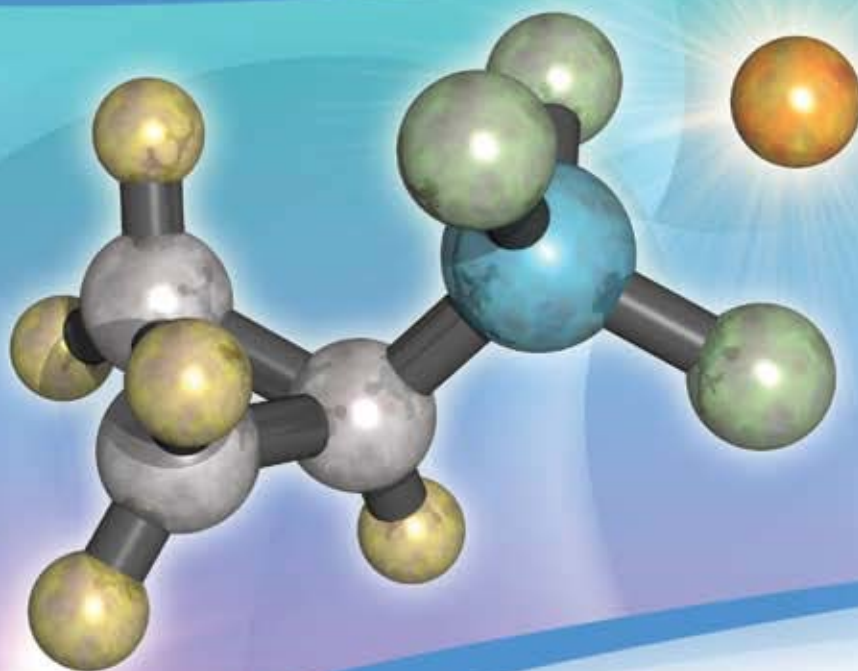


Boron Reagents for Suzuki Coupling



BORYLATION REAGENTS

BORONIC ACIDS

BORONATE ESTERS

TRIFLUOROBORATE SALTS

Potassium cyclopropyltrifluoroborate: a simple reagent for the introduction of a cyclopropyl group via Suzuki cross-coupling.

Table of Contents

Sigma-Aldrich is committed to providing the most extensive portfolio of high-quality boronic acids, boronate esters, and trifluoroborate salts for use in Suzuki coupling, and we continually expand our product listing. Within each section of this directory, products are listed by increasing molecular formula. If you are unable to find a coupling reagent for your research, "Please Bother Us" at nwallock@sial.com, or contact your local Sigma-Aldrich office (see back cover).

If viewing the electronic version, simply select from the category below to jump to that section or activate your Adobe® Bookmarks. You may also search by name, product number, molecular formula, or CAS registry number by using the "find" feature in Adobe® (Ctrl+F in Windows® or Command+F in a Macintosh® environment).

Borylation Reagents

Boronic Acids

Alkylboronic acids
Alkenylboronic acids
Arylboronic acids
N-Heteroarylboronic acids
O-Heteroarylboronic acids
S-Heteroarylboronic acids

Boronate Esters

Alkylboronate esters
Alkenylboronate esters
Allenyl- and Alkynylboronate esters
Arylboronate esters
N-Heteroarylboronate esters
O-Heteroarylboronate esters
S-Heteroarylboronate esters

Trifluoroborate Salts

Alkyltrifluoroborates
Alkenyltrifluoroborates
Aryltrifluoroborates
S-Heteroaryltrifluoroborates

Foreword



Akira Suzuki
Professor Emeritus, Hokkaido University
Sapporo, Japan

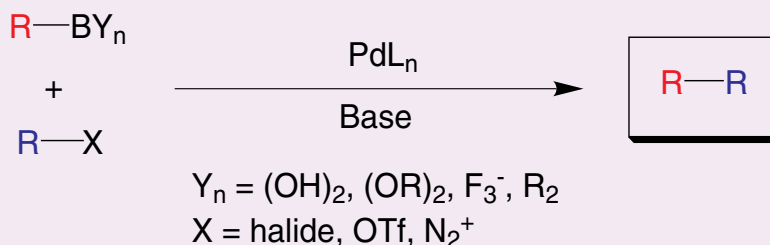
Reactions in which new carbon–carbon bonds are formed are key steps in building the complex, bioactive molecules that become developed into drugs and agrochemicals. These reactions are also vital in developing the new generation of ingeniously designed organic materials with novel electronic, optical, or mechanical properties, that are likely to play a significant role in the burgeoning area of nanotechnology.

During the past 40 years, the most important carbon–carbon bond-forming methodologies have involved the use of transition metals to mediate the reactions in a controlled and selective manner. Perhaps the most widely used of these chemistries is the cross-coupling reaction of various organoboron compounds with organic electrophiles catalyzed by a palladium complex. This reaction, Suzuki coupling, is now regarded as an integral part of innumerable synthetic routes used to build complex organic chemicals.

Suzuki coupling has many advantages. The reactants are readily available, nontoxic, and air- and water-stable. They react under mild conditions and are amenable to a variety of reaction conditions, including the use of aqueous solvents and substrate supports. The inorganic boron byproducts can be easily removed after completion of the reaction. Most important of all, the coupling proceeds with high regio- and stereoselectivity, and is little affected by steric hindrance. It does not affect other functional groups in the molecule, and can be used in one-pot strategies.

The reaction has proved to be extremely versatile. It was first carried out between alkenyl reactants, but over the years, the scope of the reaction has been extended to couple carbon atoms in aryl, alkyl, and alkynyl groups under a wide variety of conditions. Thus, the Suzuki coupling reaction is truly "carbon–carbon bond formation made easy."

The Suzuki Coupling Reaction



TO ORDER: Contact your local Sigma-Aldrich office (see back cover), or visit sigma-aldrich.com/chemicalsynthesis.

Borylation Reagents

4,4,5,5-Tetramethyl-1,3,2-dioxaborolane, sold by volume

456462

Pinacolborane
C₆H₁₃BO₂
MW: 127.98
[25015-63-8]



4,4,5,5-Tetramethyl-1,3,2-dioxaborolane, sold by weight

655856

Pinacolborane
C₆H₁₃BO₂
MW: 127.98
[25015-63-8]



4,4,5,5-Tetramethyl-1,3,2-dioxaborolane solution, 1.0 M in tetrahydrofuran

458945

Pinacolborane
C₆H₁₃BO₂
MW: 127.98
[25015-63-8]



Catecholborane solution, 1.0 M in tetrahydrofuran

225762

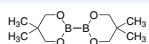
C₆H₈BO₂
MW: 119.91
[274-07-7]



Bis(neopentyl glycolato)diboron

518808

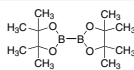
C₁₀H₂₀B₂O₄
MW: 225.89
[201733-56-4]



Bis(pinacolato)diboron

473294

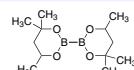
C₁₂H₂₄B₂O₄
MW: 253.94
[73183-34-3]



Bis(hexylene glycolato)diboron

525685

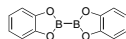
C₁₂H₂₄B₂O₄
MW: 253.94
[230299-21-5]



Bis(catecholato)diboron

473286

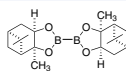
C₁₂H₈B₂O₄
MW: 237.81
[13826-27-2]



Bis[(+)-pinanediolato)diboron

527130

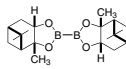
C₂₀H₃₂B₂O₄
MW: 358.09
[230299-17-9]



Bis[(-)-pinanediolato)diboron

527149

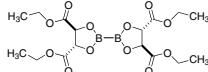
C₂₀H₃₂B₂O₄
MW: 358.09
[230299-05-5]



Bis(diethyl-D-tartrate glycolato)diboron

527157

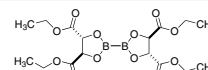
C₁₆H₂₄B₂O₁₂
MW: 429.98
[312693-46-2]



Bis(diethyl-L-tartrate glycolato)diboron

527165

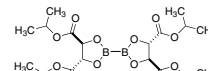
C₁₆H₂₄B₂O₁₂
MW: 429.98
[480438-20-8]



Bis(diisopropyl-D-tartrate glycolato)-diboron

527173

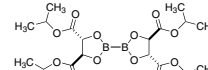
C₂₀H₃₂B₂O₁₂
MW: 486.08
[480438-21-9]



Bis(diisopropyl-L-tartrate glycolato)diboron

527181

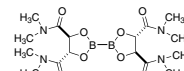
C₂₀H₃₂B₂O₁₂
MW: 486.08
[230299-10-2]



Bis(N,N,N',N'-tetramethyl-D-tartaramide glycolato)diboron

527203

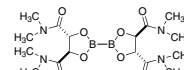
C₁₆H₂₈B₂N₄O₈
MW: 426.04
[230299-42-0]



Bis(N,N,N',N'-tetramethyl-L-tartaramide glycolato)diboron

527238

C₁₆H₂₈B₂N₄O₈
MW: 426.04
[480438-22-0]



Boric Acids

Alkylboronic acids

Methylboronic acid

165336

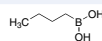
CH₃BO₂
MW: 59.86
[13061-96-6]



Butylboronic acid

163244

C₄H₁₁BO₂
MW: 101.94
[4426-47-5]



3-Methyl-1-butylboronic acid

655090

C₅H₁₃BO₂
MW: 115.97
[98139-72-1]



Cyclopropylboronic acid

597988

C₃H₇BO₂
MW: 85.90
[411235-57-9]



(2-Methylpropyl)boronic acid

346225

C₄H₁₁BO₂
MW: 101.94
[84110-40-7]



Cyclohexylboronic acid

556580

C₆H₁₃BO₂
MW: 127.98
[4441-56-9]



Isopropylboronic acid

648787

C₃H₇BO₂
MW: 87.91
[80041-89-0]



Cyclopentylboronic acid

588415

C₅H₁₁BO₂
MW: 113.95
[63076-51-7]



Phenethylboronic acid

588423

C₈H₁₁BO₂
MW: 149.98
[34420-17-2]



Cyclobutylboronic acid

646598

C₄H₉BO₂
MW: 99.92
[849052-26-2]



Neopentylboronic acid

683671

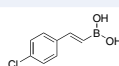
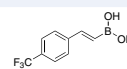
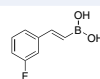
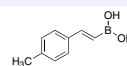
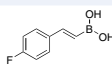
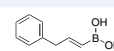
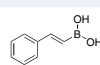
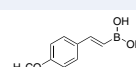
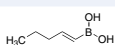
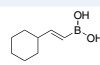
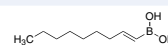
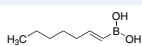
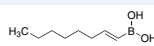
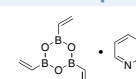
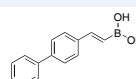
C₅H₁₃BO₂
MW: 115.97
[701261-35-0]



NEW

Ready to scale up? For competitive quotes on larger quantities or custom synthesis, contact your local Sigma-Aldrich office, or visit safglobal.com.

Alkenylboronic acids

trans-2-Chloromethylvinylboronic acid**556599**C₃H₆BClO₂
MW: 120.34
[215951-86-3]**trans-2-(4-Chlorophenyl)vinylboronic acid****523569**C₈H₈BClO₂
MW: 182.41
[154230-29-2]**trans-2-(4-(Trifluoromethyl)phenyl)vinylboronic acid****519022**C₉H₈BF₃O₂
MW: 215.96
[352525-91-8]**cis-1-Propene-1-boronic acid****572179**C₃H₇BO₂
MW: 85.90
[7547-96-8]**trans-2-(3-Fluorophenyl)vinylboronic acid****637971**C₈H₈BFO₂
MW: 165.96
[849062-22-2]**trans-2-(4-Methylphenyl)vinylboronic acid****568139**C₉H₁₁BO₂
MW: 161.99
[72316-17-7]**trans-1-Propen-1-ylboronic acid****576638**C₃H₇BO₂
MW: 85.90
[7547-97-9]**trans-2-(4-Fluorophenyl)vinylboronic acid****518972**C₈H₈BFO₂
MW: 165.96
[214907-24-1]**trans-3-Phenyl-1-propen-1-ylboronic acid****642606**C₉H₁₁BO₂
MW: 161.99
[129423-29-6]**trans-2-Buten-2-ylboronic acid****638056**C₄H₉BO₂
MW: 99.92
[125261-72-5]**trans-2-Phenylvinylboronic acid****473790**C₈H₉BO₂
MW: 147.97
[6783-05-7]**trans-2-(4-Methoxyphenyl)vinylboronic acid****518980**C₉H₁₁BO₃
MW: 177.99
[72316-18-8]**3-Methyl-2-buten-2-ylboronic acid****639079**C₅H₁₁BO₂
MW: 113.95
[870777-16-5]**1-Phenylvinylboronic acid****571350**C₈H₉BO₂
MW: 147.97
[14900-39-1]**trans-1-Penten-1-ylboronic acid****578452**C₅H₁₁BO₂
MW: 113.95
[104376-24-1]**trans-2-Cyclohexylvinylboronic acid****596256**C₈H₁₅BO₂
MW: 154.01
[37490-33-8]**trans-1-Nonenylboronic acid****579394**C₉H₁₉BO₂
MW: 170.06
[57404-77-0]**trans-1-Heptenylboronic acid****579386**C₇H₁₅BO₂
MW: 142.00**trans-1-Octen-1-ylboronic acid****521027**C₈H₁₇BO₂
MW: 156.03
[42599-16-6]**Vinylboronic anhydride pyridine complex****637998**C₁₁H₁₄B₃NO₂₃
MW: 240.67
[442850-89-7]**trans-2-(4-Biphenyl)vinylboronic acid****526045**C₁₄H₁₃BO₂
MW: 224.06
[352530-23-5]

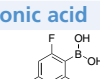
Arylboronic acids

4-Bromo-2,3,5,6-tetrafluorophenylboronic acid**593966**C₆H₂BBrF₄O₂
MW: 272.79**2,3,4,6-Tetrafluorophenylboronic acid**

NEW

680621C₆H₃BF₄O₂
MW: 193.89
[511295-00-4]**3-Bromo-2,6-difluorophenylboronic acid****557242**C₆H₄BBrF₂O₂
MW: 236.81
[352535-84-3]**Phenyl-d₅-boronic acid****517860**C₆H₂BD₅O₂
MW: 126.96
[215527-70-1]**6-Bromo-2-fluoro-3-iodophenylboronic acid**

NEW

666793C₆H₄BBrFO₂
MW: 344.71**6-Bromo-2,3-difluorophenylboronic acid****635774**C₆H₄BBrF₂O₂
MW: 236.81
[870718-10-8]**Pentafluorophenylboronic acid****465097**C₆H₂BF₅O₂
MW: 211.88
[1582-24-7]**4-Bromo-2,6-difluorophenylboronic acid****557218**C₆H₄BBrF₂O₂
MW: 236.81**2-Bromo-4,5-difluorophenylboronic acid****645281**C₆H₄BBrF₂O₂
MW: 236.81
[849062-34-6]

3,6-Dibromo-2-fluorophenylboronic acid

651087
 $C_6H_4BBr_2FO_2$
 MW: 297.71
 [870778-92-0]

**3-Bromo-5-fluorophenylboronic acid**

645346
 $C_6H_5BBrFO_2$
 MW: 218.82
 [849062-37-9]

**2-Fluoro-3-iodophenylboronic acid**

593524
 $C_6H_5BFO_2$
 MW: 265.82

**2,4-Dibromo-6-fluorophenylboronic acid**

651257
 $C_6H_4BBr_2FO_2$
 MW: 297.71
 [870778-96-4]

**3,5-Dibromophenylboronic acid**

499501
 $C_6H_5BBr_2O_2$
 MW: 279.72
 [117695-55-3]

**2-Fluoro-6-iodophenylboronic acid**

639451
 $C_6H_5BFO_2$
 MW: 265.82
 [870777-22-3]

**3,5-Dibromo-2-fluorophenylboronic acid**

661937
 $C_6H_4BBr_2FO_2$
 MW: 297.71

**3-Chloro-4-fluorophenylboronic acid**

512230
 $C_6H_5BClFO_2$
 MW: 174.37
 [144432-85-9]

**2,4-Difluorophenylboronic acid**

465070
 $C_6H_4BF_2O_2$
 MW: 157.91
 [144025-03-6]

**2,3,4-Trifluorophenylboronic acid**

524085
 $C_6H_4BF_3O_2$
 MW: 175.90
 [226396-32-3]

**3-Chloro-2-fluorophenylboronic acid**

557226
 $C_6H_5BClFO_2$
 MW: 174.37

**3,4-Difluorophenylboronic acid**

465089
 $C_6H_5BF_2O_2$
 MW: 157.91
 [168267-41-2]

**2,3,6-Trifluorophenylboronic acid**

524093
 $C_6H_4BF_3O_2$
 MW: 175.90
 [247564-71-2]

**5-Chloro-2-fluorophenylboronic acid**

557234
 $C_6H_5BClFO_2$
 MW: 174.37

**2,6-Difluorophenylboronic acid**

470791
 $C_6H_4BF_2O_2$
 MW: 157.91
 [162101-25-9]

**2,4,6-Trifluorophenylboronic acid**

524107
 $C_6H_4BF_3O_2$
 MW: 175.90
 [182482-25-3]

**2-Chloro-6-fluorophenylboronic acid**

566071
 $C_6H_5BClFO_2$
 MW: 174.37
 [313545-32-3]

**3,5-Difluorophenylboronic acid**

471925
 $C_6H_5BF_2O_2$
 MW: 157.91
 [156545-07-2]

**2,4,5-Trifluorophenylboronic acid**

524689
 $C_6H_4BF_3O_2$
 MW: 175.90
 [247564-72-3]

**3,5-Dichlorophenylboronic acid**

445207
 $C_6H_5BCl_2O_2$
 MW: 190.82
 [67492-50-6]

**2,5-Difluorophenylboronic acid**

514020
 $C_6H_5BF_2O_2$
 MW: 157.91
 [193353-34-3]

**2,3,5-Trifluorophenylboronic acid**

524697
 $C_6H_4BF_3O_2$
 MW: 175.90
 [247564-73-4]

**3,4-Dichlorophenylboronic acid**

471917
 $C_6H_5BCl_2O_2$
 MW: 190.82
 [151169-75-4]

**2,3-Difluorophenylboronic acid**

514039
 $C_6H_5BF_2O_2$
 MW: 157.91
 [121219-16-7]

**3,4,5-Trifluorophenylboronic acid**

524700
 $C_6H_4BF_3O_2$
 MW: 175.90
 [143418-49-9]

**2,5-Dichlorophenylboronic acid**

512311
 $C_6H_5BCl_2O_2$
 MW: 190.82
 [135145-90-3]

**2-Bromophenylboronic acid**

473804
 $C_6H_6BBrO_2$
 MW: 200.83
 [244205-40-1]

**3-Bromo-2-fluorophenylboronic acid**

558214
 $C_6H_5BBrFO_2$
 MW: 218.82
 [352535-97-8]

**2,3-Dichlorophenylboronic acid**

514047
 $C_6H_5BCl_2O_2$
 MW: 190.82
 [151169-74-3]

**3-Bromophenylboronic acid**

441627
 $C_6H_6BBrO_2$
 MW: 200.83
 [89598-96-9]

**5-Bromo-2-fluorophenylboronic acid**

593621
 $C_6H_5BBrFO_2$
 MW: 218.82
 [112204-57-6]

**2,4-Dichlorophenylboronic acid**

521388
 $C_6H_5BCl_2O_2$
 MW: 190.82
 [68716-47-2]

**4-Bromophenylboronic acid**

B75956
 $C_6H_6BBrO_2$
 MW: 200.83
 [5467-74-3]

**2-Bromo-6-fluorophenylboronic acid**

593737
 $C_6H_5BBrFO_2$
 MW: 218.82

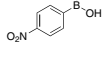
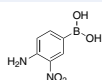
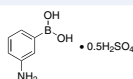
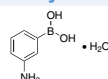
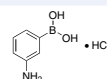
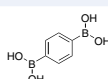
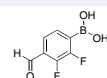
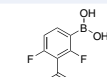
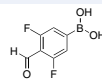
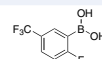
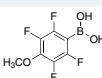
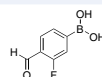
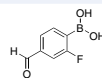
**2-Fluoro-5-iodophenylboronic acid**

593419
 $C_6H_5BFO_2$
 MW: 265.82

**2-Chlorophenylboronic acid**

445215
 $C_6H_6BClO_2$
 MW: 156.37
 [3900-89-8]



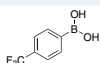
3-Chlorophenylboronic acid**417521**C₆H₆BClO₂
MW: 156.37
[63503-60-6]**4-Chlorophenylboronic acid****417548**C₆H₆BClO₂
MW: 156.37
[1679-18-1]**2-Fluorophenylboronic acid****445223**C₆H₆BF₂O₂
MW: 139.92
[1993-03-9]**3-Fluorophenylboronic acid****441643**C₆H₆BF₂O₂
MW: 139.92
[768-35-4]**4-Fluorophenylboronic acid****417556**C₆H₆BF₂O₂
MW: 139.92
[1765-93-1]**3-Iodophenylboronic acid****441678**C₆H₆BIO₂
MW: 247.83
[221037-98-5]**4-Iodophenylboronic acid****471933**C₆H₆BIO₂
MW: 247.83
[5122-99-6]**2-Nitrophenylboronic acid****673862**C₆H₆BNO₄
MW: 166.93
[5570-19-4]**3-Nitrophenylboronic acid****325104**C₆H₆BNO₄
MW: 166.93
[13331-27-6]**4-Nitrophenylboronic acid****673854**C₆H₆BNO₄
MW: 166.93
[24067-17-2]**4-Amino-3-nitrophenylboronic acid****651621**C₆H₇BN₂O₄
MW: 181.94
[89466-07-9]**Phenylboronic acid****P20009**C₆H₇BO₂
MW: 121.93
[98-80-6]**3-Mercaptophenylboronic acid****523992**C₆H₇BO₂S
MW: 153.99
[352526-01-3]**4-Mercaptophenylboronic acid****524018**C₆H₇BO₂S
MW: 153.99
[237429-33-3]**3-Hydroxyphenylboronic acid****523968**C₆H₇BO₃
MW: 137.93
[87199-18-6]**4-Hydroxyphenylboronic acid****523976**C₆H₇BO₃
MW: 137.93
[71597-85-8]**3-Aminophenylboronic acid hemisulfate****A71751**C₆H₈BNO₂ • 0.5H₂SO₄
MW: 185.98
[66472-86-4]**3-Aminophenylboronic acid monohydrate****287512**C₆H₈BNO₂ • H₂O
MW: 154.96
[206658-89-1]**3-Aminophenylboronic acid hydrochloride****410705**C₆H₈BNO₂ • HCl
MW: 173.41
[85006-23-1]**Benzene-1,4-diboronic acid****417130**C₆H₈B₂O₄
MW: 165.75
[4612-26-4]**2,3-Difluoro-4-formylphenylboronic acid****571377**C₇H₅BF₂O₃
MW: 185.92
[480424-84-8]**2,4-Difluoro-3-formylphenylboronic acid****597406**C₇H₅BF₂O₃
MW: 185.92
[870718-06-2]**3,5-Difluoro-4-formylphenylboronic acid****635782**C₇H₅BF₂O₃
MW: 185.92
[870718-11-9]**2,6-Difluoro-3-formylphenylboronic acid****652210**C₇H₅BF₂O₃
MW: 185.92
[849062-09-5]**2-Fluoro-5-(trifluoromethyl)phenylboronic acid****558184**C₇H₅BF₄O₂
MW: 207.92
[352535-96-7]**2-Fluoro-6-(trifluoromethyl)phenylboronic acid****558192**C₇H₅BF₄O₂
MW: 207.92
[313545-34-5]**2-Fluoro-3-(trifluoromethyl)phenylboronic acid****558206**C₇H₅BF₄O₂
MW: 207.92
[157834-21-4]**4-Methoxy-2,3,5,6-tetrafluorophenylboronic acid****657301**C₇H₅BF₄O₃
MW: 223.92
[871126-20-4]**3-Fluoro-4-formylphenylboronic acid****542296**C₇H₆BFO₃
MW: 167.93
[248270-25-9]**2-Fluoro-3-formylphenylboronic acid****645788**C₇H₆BFO₃
MW: 167.93
[849061-98-9]**2-Fluoro-4-formylphenylboronic acid****657344**C₇H₆BFO₃
MW: 167.93
[871126-22-6]**2-(Trifluoromethyl)phenylboronic acid****445193**C₇H₆BF₃O₂
MW: 189.93
[1423-27-4]

3-(Trifluoromethyl)phenylboronic acid**432032**

$C_7H_6BF_3O_2$
 MW: 189.93
 [1423-26-3]

**4-(Trifluoromethyl)phenylboronic acid****439320**

$C_7H_6BF_3O_2$
 MW: 189.93
 [128796-39-4]

**2-(Trifluoromethoxy)phenylboronic acid****NEW****683787**

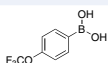
$C_7H_6BF_3O_3$
 MW: 205.93
 [175676-65-0]

**3-(Trifluoromethoxy)phenylboronic acid****510122**

$C_7H_6BF_3O_3$
 MW: 205.93
 [179113-90-7]

**4-(Trifluoromethoxy)phenylboronic acid****510130**

$C_7H_6BF_3O_3$
 MW: 205.93
 [139301-27-2]

**3-Methoxy-2,4,6-trifluorophenylboronic acid****652202**

$C_7H_6BF_3O_3$
 MW: 205.93
 [849062-08-4]

**2-Cyanophenylboronic acid****521396**

$C_7H_6BNO_2$
 MW: 146.94
 [138642-62-3]

**3-Cyanophenylboronic acid****513016**

$C_7H_6BNO_2$
 MW: 146.94
 [150255-96-2]

**4-Cyanophenylboronic acid****521418**

$C_7H_6BNO_2$
 MW: 146.94
 [126747-14-6]

**3-Bromo-5-fluoro-2-methoxyphenylboronic acid****645176**

$C_7H_7BBrFO_3$
 MW: 248.84
 [352525-85-0]

**2-Bromo-6-fluoro-3-methoxyphenylboronic acid****662011**

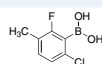
$C_7H_7BBrFO_3$
 MW: 248.84

**2-Chloro-6-fluoro-3-methylphenylboronic acid****557250**

$C_7H_7BClFO_2$
 MW: 188.39
 [352535-85-4]

**2-Chloro-6-fluoro-5-methylphenylboronic acid****557269**

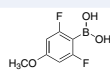
$C_7H_7BClFO_2$
 MW: 188.39
 [352535-86-5]

**5-Chloro-2-fluoro-3-methylphenylboronic acid****557277**

$C_7H_7BClFO_2$
 MW: 188.39
 [352535-87-6]

**2,6-Difluoro-4-methoxyphenylboronic acid****593060**

$C_7H_7BF_2O_3$
 MW: 187.94
 [406482-20-0]

**4,5-Difluoro-2-methoxyphenylboronic acid****645184**

$C_7H_7BF_2O_3$
 MW: 187.94
 [870777-32-5]

**2,6-Difluoro-3-methoxyphenylboronic acid****652164**

$C_7H_7BF_2O_3$
 MW: 187.94
 [870779-02-5]

**3,5-Diiodo-2-methoxyphenylboronic acid****NEW****666165**

$C_7H_7BI_2O_3$
 MW: 403.75

**2-Formylphenylboronic acid****431958**

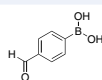
$C_7H_7BO_3$
 MW: 149.94
 [40138-16-7]

**3-Formylphenylboronic acid****441651**

$C_7H_7BO_3$
 MW: 149.94
 [87199-16-4]

**4-Formylphenylboronic acid****431966**

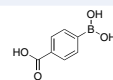
$C_7H_7BO_3$
 MW: 149.94
 [87199-17-5]

**3-Carboxyphenylboronic acid****456764**

$C_7H_7BO_4$
 MW: 165.94
 [25487-66-5]

**4-Carboxyphenylboronic acid****456772**

$C_7H_7BO_4$
 MW: 165.94
 [14047-29-1]

**3,4-(Methylenedioxy)phenylboronic acid****499994**

$C_7H_7BO_4$
 MW: 165.94
 [94839-07-3]

**2-(Bromomethyl)phenylboronic acid****NEW****679453**

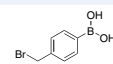
$C_7H_8BBrO_2$
 MW: 214.85
 [91983-14-1]

**3-(Bromomethyl)phenylboronic acid****NEW****679445**

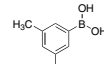
$C_7H_8BBrO_2$
 MW: 214.85
 [51323-43-4]

**4-(Bromomethyl)phenylboronic acid****NEW****679437**

$C_7H_8BBrO_2$
 MW: 214.85
 [68162-47-0]

**3-Bromo-5-methylphenylboronic acid****645311**

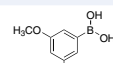
$C_7H_8BBrO_2$
 MW: 214.85
 [849062-36-8]

**5-Bromo-2-methoxyphenylboronic acid****512702**

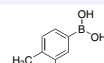
$C_7H_8BBrO_3$
 MW: 230.85
 [89694-45-1]

**3-Bromo-5-methoxyphenylboronic acid****652326**

$C_7H_8BBrO_3$
 MW: 230.85
 [849062-12-0]

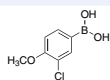
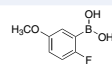
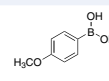
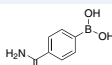
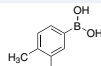
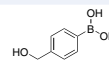
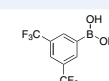
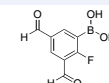
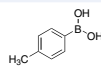
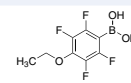
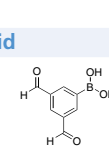
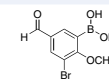
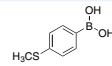
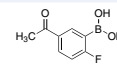
**3-Chloro-4-methylphenylboronic acid****NEW****680559**

$C_7H_8BClO_2$
 MW: 170.40
 [175883-63-3]

**2-Chloro-6-methoxyphenylboronic acid****512753**

$C_7H_8BClO_3$
 MW: 180.35
 [385370-80-9]



5-Chloro-2-methoxyphenylboronic acid**512249**C₇H₈BClO₃
MW: 186.40
[89694-48-4]**4-Fluoro-2-methoxyphenylboronic acid****564494**C₇H₈BFO₃
MW: 169.95
[179899-07-1]**2-Methoxyphenylboronic acid****445231**C₇H₉BO₃
MW: 151.96
[5720-06-9]**3-Chloro-4-methoxyphenylboronic acid****564486**C₇H₈BClO₃
MW: 186.40
[175883-60-0]**2-Fluoro-3-methoxyphenylboronic acid****594253**C₇H₈BFO₃
MW: 169.95
[352303-67-4]**3-Methoxyphenylboronic acid****441686**C₇H₉BO₃
MW: 151.96
[10365-98-7]**4-Fluoro-3-methylphenylboronic acid****483567**C₇H₈BFO₂
MW: 153.95
[139911-27-6]**2-Fluoro-5-methoxyphenylboronic acid****597112**C₇H₈BFO₃
MW: 169.95
[406482-19-7]**4-Methoxyphenylboronic acid****417599**C₇H₉BO₃
MW: 151.96
[5720-07-0]**4-Fluoro-2-methylphenylboronic acid****565652**C₇H₈BFO₂
MW: 153.95
[139911-29-8]**4-Aminocarbonylphenylboronic acid** NEW**683876**C₇H₈BNO₃
MW: 164.95
[123088-59-5]**3-(Hydroxymethyl)phenylboronic acid****512834**C₇H₉BO₃
MW: 151.96
[87199-15-3]**2-Fluoro-4-methylphenylboronic acid****567418**C₇H₈BFO₂
MW: 153.95
[170981-26-7]**4-Methyl-3-nitrophenylboronic acid****521477**C₇H₈BNO₄
MW: 180.95
[80500-27-2]**4-(Hydroxymethyl)phenylboronic acid****512338**C₇H₉BO₃
MW: 151.96
[59016-93-2]**3-Fluoro-4-methylphenylboronic acid****567426**C₇H₈BFO₂
MW: 153.95
[168267-99-0]**o-Tolylboronic acid****393606**C₇H₉BO₂
MW: 135.96
[16419-60-6]**3,5-Bis(trifluoromethyl)phenylboronic acid****471070**C₈H₅BF₆O₂
MW: 257.93
[73852-19-4]**2-Fluoro-5-methylphenylboronic acid****567434**C₇H₈BFO₂
MW: 153.95
[166328-16-1]**m-Tolylboronic acid****393614**C₇H₉BO₂
MW: 135.96
[17933-03-8]**2-Fluoro-3,5-diformylphenylboronic acid****645796**C₈H₆BFO₄
MW: 195.94
[870778-85-1]**5-Fluoro-2-methylphenylboronic acid****567442**C₇H₈BFO₂
MW: 153.95
[163517-62-2]**p-Tolylboronic acid****393622**C₇H₉BO₂
MW: 135.96
[5720-05-8]**4-Ethoxy-2,3,5,6-tetrafluorophenylboronic acid****657298**C₈H₇BF₄O₃
MW: 237.94
[871125-72-3]**5-Fluoro-2-methoxyphenylboronic acid****483540**C₇H₈BFO₃
MW: 169.95
[179897-94-0]**2-(Methylthio)phenylboronic acid****521248**C₇H₉BO₂S
MW: 168.02
[168618-42-6]**3,5-Diformylphenylboronic acid****567485**C₈H₇BO₄
MW: 177.95
[480424-62-2]**2-Fluoro-6-methoxyphenylboronic acid****512761**C₇H₈BFO₃
MW: 169.95
[78495-63-3]**3-(Methylthio)phenylboronic acid****526002**C₇H₉BO₂S
MW: 168.02
[128312-11-8]**3-Bromo-5-formyl-2-methoxyphenylboronic acid** NEW**680664**C₈H₈BBrO₄
MW: 258.86**3-Fluoro-4-methoxyphenylboronic acid****564036**C₇H₈BFO₃
MW: 169.95
[149507-26-6]**4-(Methylthio)phenylboronic acid****456802**C₇H₉BO₂S
MW: 168.02
[98546-51-1]**2-Fluoro-5-acetylphenylboronic acid****639583**C₈H₈BFO₃
MW: 181.96
[870777-29-0]

3-Acetyl-2-fluorophenylboronic acid

651249
 $C_8H_8BFO_3$
 MW: 181.96
 [870778-95-3]

**5-Fluoro-3-formyl-2-methoxyphenylboronic acid**

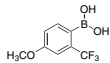
680699
 $C_8H_8BFO_4$
 MW: 197.96

**3-Ethoxy-2,4,6-trifluorophenylboronic acid**

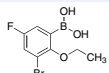
657247
 $C_8H_8BF_3O_3$
 MW: 219.95
 [871125-69-8]

**4-Methoxy-2-(trifluoromethyl)phenylboronic acid**

666912
 $C_8H_8BF_3O_3$
 MW: 219.95
 [313546-16-6]

**3-Bromo-2-ethoxy-5-fluorophenylboronic acid**

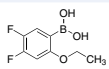
542512
 $C_8H_8BBrFO_3$
 MW: 262.87
 [352534-82-8]

**2-Bromo-3-ethoxy-6-fluorophenylboronic acid**

645710
 $C_8H_8BBrFO_3$
 MW: 262.87
 [849052-19-3]

**4,5-Difluoro-2-ethoxyphenylboronic acid**

650943
 $C_8H_8BF_2O_3$
 MW: 201.96
 [870778-87-3]

**2,6-Difluoro-3-ethoxyphenylboronic acid**

650951
 $C_8H_8BF_2O_3$
 MW: 201.96
 [849062-00-6]

**3-Vinylphenylboronic acid**

680575
 $C_8H_9BO_2$
 MW: 147.97
 [15016-43-0]

**4-Vinylphenylboronic acid**

417580
 $C_8H_9BO_2$
 MW: 147.97
 [2156-04-9]

**2-Acetylphenylboronic acid**

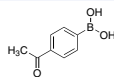
470805
 $C_8H_9BO_3$
 MW: 163.97
 [308103-40-4]

**3-Acetylphenylboronic acid**

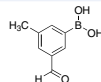
470813
 $C_8H_9BO_3$
 MW: 163.97
 [204841-19-0]

**4-Acetylphenylboronic acid**

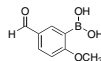
470821
 $C_8H_9BO_3$
 MW: 163.97
 [149104-90-5]

**3-Formyl-5-methylphenylboronic acid**

645338
 $C_8H_9BO_3$
 MW: 163.97
 [870777-33-6]

**5-Formyl-2-methoxyphenylboronic acid**

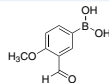
512257
 $C_8H_9BO_4$
 MW: 179.97
 [127972-02-5]

**3-Formyl-2-methoxyphenylboronic acid**

666157
 $C_8H_9BO_4$
 MW: 179.97

**3-Formyl-4-methoxyphenylboronic acid**

512869
 $C_8H_9BO_4$
 MW: 179.97
 [121124-97-8]

**2-Methoxycarbonylphenylboronic acid**

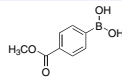
683809
 $C_8H_9BO_4$
 MW: 179.97
 [374538-03-1]

**3-Methoxycarbonylphenylboronic acid**

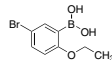
591130
 $C_8H_9BO_4$
 MW: 179.97
 [99769-19-4]

**4-Methoxycarbonylphenylboronic acid**

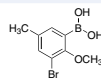
594539
 $C_8H_9BO_4$
 MW: 179.97
 [99768-12-4]

**5-Bromo-2-ethoxyphenylboronic acid**

512826
 $C_8H_9BBrO_3$
 MW: 244.88
 [352525-82-7]

**3-Bromo-2-methoxy-5-methylphenylboronic acid**

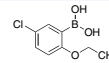
596183
 $C_8H_9BBrO_3$
 MW: 244.88
 [870717-99-0]

**5-Bromo-3-ethoxyphenylboronic acid**

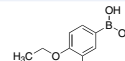
651192
 $C_8H_9BBrO_3$
 MW: 244.88
 [849062-02-8]

**5-Chloro-2-ethoxyphenylboronic acid**

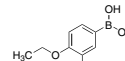
542547
 $C_8H_9BClO_3$
 MW: 200.43
 [352534-86-2]

**3-Chloro-4-ethoxyphenylboronic acid**

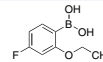
564478
 $C_8H_9BClO_3$
 MW: 200.43
 [279261-81-3]

**4-Ethoxy-3-fluorophenylboronic acid**

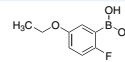
564044
 $C_8H_9BFO_3$
 MW: 183.97
 [279263-10-4]

**2-Ethoxy-4-fluorophenylboronic acid**

564508
 $C_8H_9BFO_3$
 MW: 183.97
 [480438-58-2]

**5-Ethoxy-2-fluorophenylboronic acid**

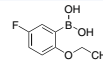
594393
 $C_8H_9BFO_3$
 MW: 183.97

**2-Ethoxy-6-fluorophenylboronic acid**

594504
 $C_8H_9BFO_3$
 MW: 183.97

**2-Ethoxy-5-fluorophenylboronic acid**

594954
 $C_8H_9BFO_3$
 MW: 183.97

**3-Ethoxy-2-fluorophenylboronic acid**

639389
 $C_8H_9BFO_3$
 MW: 183.97
 [855230-61-4]

**2-Acetamidophenylboronic acid**

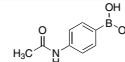
683906
 $C_8H_9BNO_3$
 MW: 178.98
 [169760-16-1]

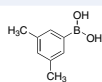
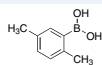
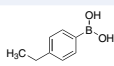
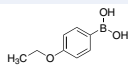
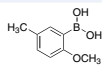
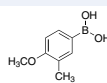
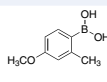
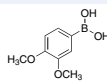
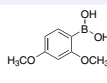
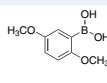
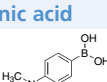
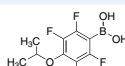
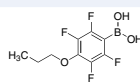
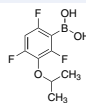
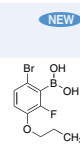
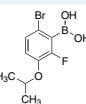
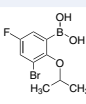
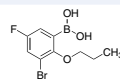
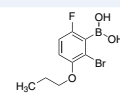
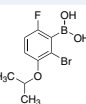
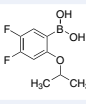
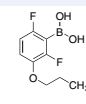
**3-Acetamidophenylboronic acid**

566012
 $C_8H_9BNO_3$
 MW: 178.98
 [78887-39-5]

**4-Acetamidophenylboronic acid**

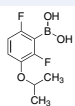
565806
 $C_8H_9BNO_3$
 MW: 178.98
 [101251-09-6]



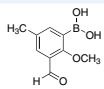
2,6-Dimethylphenylboronic acid**480061**C₈H₁₁BO₂
MW: 149.98
[100379-00-8]**3,5-Dimethylphenylboronic acid****480088**C₈H₁₁BO₂
MW: 149.98
[172975-69-8]**2,3-Dimethylphenylboronic acid****483508**C₈H₁₁BO₂
MW: 149.98
[183158-34-1]**2,5-Dimethylphenylboronic acid****483516**C₈H₁₁BO₂
MW: 149.98
[85199-06-0]**2-Ethylphenylboronic acid****521523**C₈H₁₁BO₂
MW: 149.98
[90002-36-1]**4-Ethylphenylboronic acid****499536**C₈H₁₁BO₂
MW: 149.98
[63139-21-9]**3-(Ethylthio)phenylboronic acid****597295**C₈H₁₁BO₂S
MW: 182.05
[870718-05-1]**2-Ethoxyphenylboronic acid****455520**C₈H₁₁BO₃
MW: 165.98
[213211-69-9]**3-Ethoxyphenylboronic acid****441635**C₈H₁₁BO₃
MW: 165.98
[90555-66-1]**4-Ethoxyphenylboronic acid****455539**C₈H₁₁BO₃
MW: 165.98
[22237-13-4]**2-Methoxy-5-methylphenylboronic acid****567507**C₈H₁₁BO₃
MW: 165.98
[127972-00-3]**4-Methoxy-3-methylphenylboronic acid****595500**C₈H₁₁BO₃
MW: 165.98
[175883-62-2]**4-Methoxy-2-methylphenylboronic acid****639370**C₈H₁₁BO₃
MW: 165.98
[208399-66-0]**2,6-Dimethoxyphenylboronic acid****480096**C₈H₁₁BO₄
MW: 181.98
[23112-96-1]**3,4-Dimethoxyphenylboronic acid****480118**C₈H₁₁BO₄
MW: 181.98
[122775-35-3]**2,4-Dimethoxyphenylboronic acid****483486**C₈H₁₁BO₄
MW: 181.98
[133730-34-4]**2,5-Dimethoxyphenylboronic acid****483494**C₈H₁₁BO₄
MW: 181.98
[107099-99-0]**2,3-Dimethoxyphenylboronic acid****557730**C₈H₁₁BO₄
MW: 181.98
[40972-86-9]**3-(Dimethylamino)phenylboronic acid****680532**C₈H₁₂BNO₂
MW: 165.00
[178752-79-9]**4-(Dimethylamino)phenylboronic acid****483532**C₈H₁₂BNO₂
MW: 165.00
[28611-39-4]**4-Isopropoxy-2,3,5,6-tetrafluorophenylboronic acid****657263**C₉H₉BF₄O₃
MW: 251.97
[871126-28-2]**4-Propoxy-2,3,5,6-tetrafluorophenylboronic acid****657271**C₉H₉BF₄O₃
MW: 251.97
[871125-71-2]**3-Isopropoxy-2,4,6-trifluorophenylboronic acid****657360**C₉H₁₀BF₃O₃
MW: 233.98
[871125-73-4]**3-Propoxy-2,4,6-trifluorophenylboronic acid****657379**C₉H₁₀BF₃O₃
MW: 233.98
[871125-70-1]**6-Bromo-2-fluoro-3-propoxyphenylboronic acid****666882**C₉H₁₁BrFO₃
MW: 276.90**6-Bromo-2-fluoro-3-isopropoxyphenylboronic acid****666890**C₉H₁₁BrFO₃
MW: 276.90**3-Bromo-5-fluoro-2-isopropoxyphenylboronic acid****542520**C₉H₁₁BrFO₃
MW: 276.90
[352534-84-0]**3-Bromo-5-fluoro-2-propoxyphenylboronic acid****565784**C₉H₁₁BrFO₃
MW: 276.90
[868272-84-8]**2-Bromo-6-fluoro-3-propoxyphenylboronic acid****645729**C₉H₁₁BrFO₃
MW: 276.90
[849052-20-6]**2-Bromo-6-fluoro-3-isopropoxyphenylboronic acid****661996**C₉H₁₁BrFO₃
MW: 276.90**4,5-Difluoro-2-isopropoxyphenylboronic acid****680702**C₉H₁₁BF₂O₃
MW: 215.99**2,6-Difluoro-3-propoxyphenylboronic acid****652172**C₉H₁₁BF₂O₃
MW: 215.99
[849062-14-2]

2,6-Difluoro-3-isopropoxyphenylboronic acid

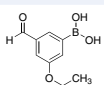
652180
 $C_9H_{11}BF_2O_3$
 MW: 215.99
 [849062-04-0]

**3-Formyl-2-methoxy-5-methylphenylboronic acid**

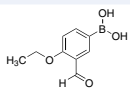
567337
 $C_9H_{11}BO_4$
 MW: 193.99
 [480424-55-3]

**3-Ethoxy-5-formylphenylboronic acid**

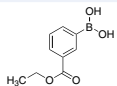
652342
 $C_9H_{11}BO_4$
 MW: 193.99

**4-Ethoxy-3-formylphenylboronic acid**

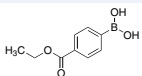
677817
 $C_9H_{11}BO_4$
 MW: 193.99
 [480424-63-3]

**3-Ethoxycarbonylphenylboronic acid**

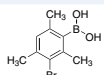
574651
 $C_9H_{11}BO_4$
 MW: 193.99
 [4334-87-6]

**4-Ethoxycarbonylphenylboronic acid**

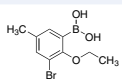
574643
 $C_9H_{11}BO_4$
 MW: 193.99
 [4334-88-7]

**3-Bromo-2,4,6-trimethylphenylboronic acid**

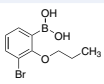
593850
 $C_9H_{12}BBrO_2$
 MW: 242.91

**3-Bromo-2-ethoxy-5-methylphenylboronic acid**

596299
 $C_9H_{12}BBrO_3$
 MW: 258.90
 [870718-00-6]

**3-Bromo-2-propoxyphenylboronic acid**

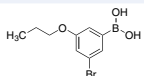
597074
 $C_9H_{12}BBrO_3$
 MW: 258.90
 [848779-86-2]

**3-Bromo-2-isopropoxyphenylboronic acid**

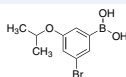
597171
 $C_9H_{12}BBrO_3$
 MW: 258.90
 [870718-04-0]

**3-Bromo-5-propoxyphenylboronic acid**

657514
 $C_9H_{12}BBrO_3$
 MW: 258.90
 [871126-27-1]

**3-Bromo-5-isopropoxyphenylboronic acid**

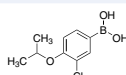
657735
 $C_9H_{12}BBrO_3$
 MW: 258.90
 [871125-81-4]

**5-Chloro-2-isopropoxyphenylboronic acid**

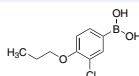
542555
 $C_9H_{12}BClO_3$
 MW: 214.45
 [352534-87-3]

**3-Chloro-4-isopropoxyphenylboronic acid**

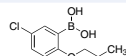
564443
 $C_9H_{12}BClO_3$
 MW: 214.45
 [480438-56-0]

**3-Chloro-4-propoxyphenylboronic acid**

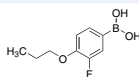
564451
 $C_9H_{12}BClO_3$
 MW: 214.45
 [480438-57-1]

**5-Chloro-2-propoxyphenylboronic acid**

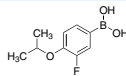
645192
 $C_9H_{12}BClO_3$
 MW: 214.45
 [849062-29-9]

**3-Fluoro-4-propoxyphenylboronic acid**

564052
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [192376-68-4]

**3-Fluoro-4-isopropoxyphenylboronic acid**

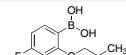
564060
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [480438-54-8]

**4-Fluoro-2-isopropoxyphenylboronic acid**

564516
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [480438-59-3]

**4-Fluoro-2-propoxyphenylboronic acid**

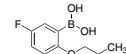
564524
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [480438-60-6]

**5-Fluoro-2-isopropoxyphenylboronic acid**

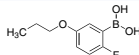
564559
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [480438-63-9]

**5-Fluoro-2-propoxyphenylboronic acid**

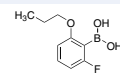
565776
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [480438-73-1]

**2-Fluoro-5-propoxyphenylboronic acid**

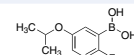
594288
 $C_9H_{12}BF_2O_3$
 MW: 198.00

**2-Fluoro-6-propoxyphenylboronic acid**

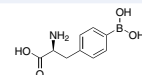
639419
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [870777-18-7]

**2-Fluoro-5-isopropoxyphenylboronic acid**

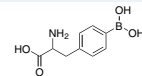
645206
 $C_9H_{12}BF_2O_3$
 MW: 198.00
 [849062-30-2]

**4-Borono-L-phenylalanine**

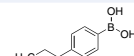
17755
 $C_9H_{12}BNO_4$
 MW: 209.01
 [76410-58-7]

**4-Borono-DL-phenylalanine**

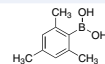
512680
 $C_9H_{12}BNO_4$
 MW: 209.01
 [90580-64-6]

**4-Propylphenylboronic acid**

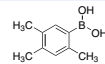
521507
 $C_9H_{13}BO_2$
 MW: 164.01
 [134150-01-9]

**2,4,6-Trimethylphenylboronic acid**

542318
 $C_9H_{13}BO_2$
 MW: 164.01
 [5980-97-2]

**2,4,5-Trimethylphenylboronic acid**

542326
 $C_9H_{13}BO_2$
 MW: 164.01
 [352534-80-6]

**2-Isopropoxyphenylboronic acid**

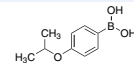
542466
 $C_9H_{13}BO_3$
 MW: 180.01
 [138008-97-6]

**3-Isopropoxyphenylboronic acid**

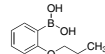
542458
 $C_9H_{13}BO_3$
 MW: 180.01
 [216485-86-8]

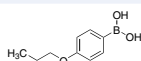
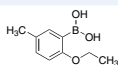
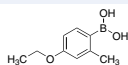
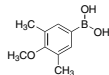
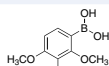
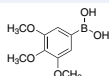
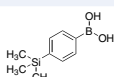
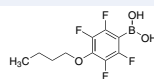
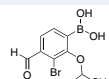
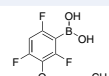
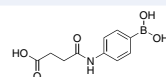
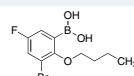
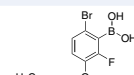
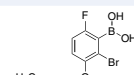
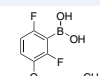
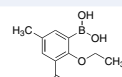
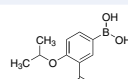
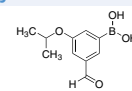
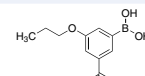
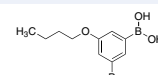
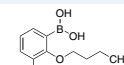
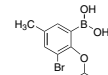
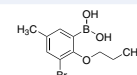
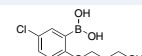
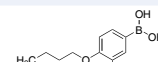
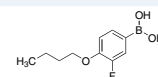
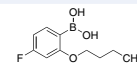
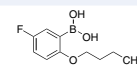
**4-Isopropoxyphenylboronic acid**

542474
 $C_9H_{13}BO_3$
 MW: 180.01
 [153624-46-5]

**2-Propoxyphenylboronic acid**

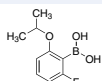
557722
 $C_9H_{13}BO_3$
 MW: 180.01
 [134896-34-7]



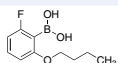
3-Propoxyphenylboronic acid**557714**C₉H₁₃BO₃
MW: 180.01
[149557-18-6]**4-Propoxyphenylboronic acid****557706**C₉H₁₃BO₃
MW: 180.01
[186497-67-6]**2-Ethoxy-5-methylphenylboronic acid****567493**C₉H₁₃BO₃
MW: 180.01
[123291-97-4]**4-Ethoxy-2-methylphenylboronic acid****594830**C₉H₁₃BO₃
MW: 180.01
[313545-31-2]**3,5-Dimethyl-4-methoxyphenylboronic acid****598062**C₉H₁₃BO₃
MW: 180.01
[301699-39-8]**2,3,4-Trimethoxyphenylboronic acid****512281**C₉H₁₃BO₅
MW: 212.01
[118062-05-8]**3,4,5-Trimethoxyphenylboronic acid****651648**C₉H₁₃BO₅
MW: 212.01
[182163-96-8]**3-(Trimethylsilyl)phenylboronic acid****523666**C₉H₁₅BO₂Si
MW: 194.11
[177171-16-3]**4-(Trimethylsilyl)phenylboronic acid****523674**C₉H₁₅BO₂Si
MW: 194.11
[17865-11-1]**1-Naphthylboronic acid****N257**C₁₀H₉BO₂
MW: 171.99
[13922-41-3]**2-Naphthylboronic acid****480134**C₁₀H₉BO₂
MW: 171.99
[32316-92-0]**4-Butoxy-2,3,5,6-tetrafluorophenylboronic acid****657255**C₁₀H₁₁BF₄O₃
MW: 266.00
[871126-19-1]**Ferroceneboronic acid****455547**C₁₀H₁₁BFeO₂
MW: 229.85
[12152-94-2]**3-Bromo-2-isopropoxy-5-formylphenylboronic acid****684619**C₁₀H₁₂BBro₃
MW: 286.91**3-Butoxy-2,4,6-trifluorophenylboronic acid****657352**C₁₀H₁₂BF₃O₃
MW: 248.01
[871126-23-7]**N-(4-Phenylboronic)succinamic acid****578754**C₁₀H₁₂BNO₅
MW: 237.02
[480424-95-1]**3-Bromo-2-butoxy-5-fluorophenylboronic acid****542539**C₁₀H₁₃BBroF₃
MW: 290.92
[352534-85-1]**6-Bromo-3-butoxy-2-fluorophenylboronic acid****661953**C₁₀H₁₃BBroF₃
MW: 290.92**2-Bromo-3-butoxy-6-fluorophenylboronic acid****661945**C₁₀H₁₃BBroF₃
MW: 290.92**3-Butoxy-2,6-difluorophenylboronic acid****652199**C₁₀H₁₃BF₂O₃
MW: 230.02
[849062-15-3]**2-Ethoxy-3-formyl-5-methylphenylboronic acid****567329**C₁₀H₁₃BO₄
MW: 208.02
[480424-54-2]**3-Formyl-4-isopropoxyphenylboronic acid****666742**C₁₀H₁₃BO₄
MW: 208.02**3-Formyl-5-isopropoxyphenylboronic acid****657557**C₁₀H₁₃BO₄
MW: 208.02
[871125-79-0]**3-Formyl-5-propoxyphenylboronic acid****657638**C₁₀H₁₃BO₄
MW: 208.02
[871125-80-3]**3-Bromo-5-butoxyphenylboronic acid****661961**C₁₀H₁₄BBro₃
MW: 272.93**3-Bromo-2-butoxyphenylboronic acid****592129**C₁₀H₁₄BBro₃
MW: 272.93
[480425-34-1]**3-Bromo-2-isopropoxy-5-methylphenylboronic acid****596620**C₁₀H₁₄BBro₃
MW: 272.93
[870718-01-7]**3-Bromo-5-methyl-2-propoxyphenylboronic acid****596744**C₁₀H₁₄BBro₃
MW: 272.93
[870718-02-8]**2-Butoxy-5-chlorophenylboronic acid****542563**C₁₀H₁₄BClO₃
MW: 228.48
[352534-88-4]**4-Butoxy-3-chlorophenylboronic acid****564435**C₁₀H₁₄BClO₃
MW: 228.48
[480438-55-9]**4-Butoxy-3-fluorophenylboronic acid****564079**C₁₀H₁₄BF₃O₃
MW: 212.03
[156487-13-7]**2-Butoxy-4-fluorophenylboronic acid****564532**C₁₀H₁₄BFO₃
MW: 212.03
[480438-61-7]**2-Butoxy-5-fluorophenylboronic acid****564540**C₁₀H₁₄BFO₃
MW: 212.03
[480438-62-8]

2-Fluoro-6-isopropoxyphenylboronic acid**639397**

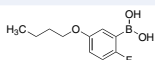
$C_{10}H_{14}BFO_3$
MW: 212.03
[870777-17-6]

**2-Butoxy-6-fluorophenylboronic acid****639427**

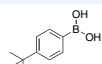
$C_{10}H_{14}BFO_3$
MW: 212.03
[870777-19-8]

**5-Butoxy-2-fluorophenylboronic acid****645230**

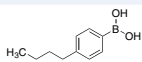
$C_{10}H_{14}BFO_3$
MW: 212.03
[849062-31-3]

**4-tert-Butylphenylboronic acid****480053**

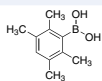
$C_{10}H_{15}BO_2$
MW: 178.04
[123324-71-0]

**4-Butylphenylboronic acid****521493**

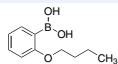
$C_{10}H_{15}BO_2$
MW: 178.04
[145240-28-4]

**2,3,5,6-Tetramethylphenylboronic acid****521515**

$C_{10}H_{15}BO_2$
MW: 178.04
[197223-36-2]

**2-Butoxyphenylboronic acid****542482**

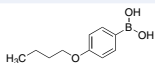
$C_{10}H_{15}BO_3$
MW: 194.04
[91129-69-0]

**3-Butoxyphenylboronic acid****542490**

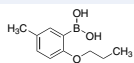
$C_{10}H_{15}BO_3$
MW: 194.04
[352534-81-7]

**4-Butoxyphenylboronic acid****542504**

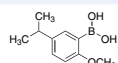
$C_{10}H_{15}BO_3$
MW: 194.04
[105365-51-3]

**5-Methyl-2-propoxyphenylboronic acid****565644**

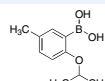
$C_{10}H_{15}BO_3$
MW: 194.04
[480438-70-8]

**5-Isopropyl-2-methoxyphenylboronic acid****680583**

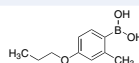
$C_{10}H_{15}BO_3$
MW: 194.04
[216393-63-4]

**2-Isopropoxy-5-methylphenylboronic acid****565660**

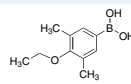
$C_{10}H_{15}BO_3$
MW: 194.04
[480438-71-9]

**4-Propoxy-2-methylphenylboronic acid****594725**

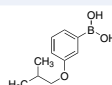
$C_{10}H_{15}BO_3$
MW: 194.04

**3,5-Dimethyl-4-ethoxyphenylboronic acid****597945**

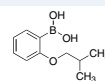
$C_{10}H_{15}BO_3$
MW: 194.04
[850568-59-1]

**3-Isobutoxyphenylboronic acid****645737**

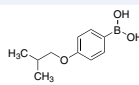
$C_{10}H_{15}BO_3$
MW: 194.04
[849052-21-7]

**2-Isobutoxyphenylboronic acid****645745**

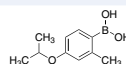
$C_{10}H_{15}BO_3$
MW: 194.04
[833486-92-3]

**4-Isobutoxyphenylboronic acid****639605**

$C_{10}H_{15}BO_3$
MW: 194.04
[153624-44-3]

**4-Isopropoxy-2-methylphenylboronic acid****657328**

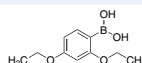
$C_{10}H_{15}BO_3$
MW: 194.04
[871126-21-5]

**2-Isopropoxy-6-methoxyphenylboronic acid****650978**

$C_{10}H_{15}BO_4$
MW: 210.03
[870778-88-4]

**2,4-Diethoxyphenylboronic acid****680680**

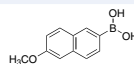
$C_{10}H_{15}BO_4$
MW: 210.03

**4-Methyl-1-naphthaleneboronic acid****521450**

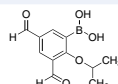
$C_{11}H_{11}BO_2$
MW: 186.01
[103986-53-4]

**6-Methoxy-2-naphthaleneboronic acid****521892**

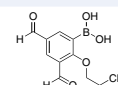
$C_{11}H_{11}BO_3$
MW: 202.01
[156641-98-4]

**3,5-Diformyl-2-isopropoxyphenylboronic acid****680648**

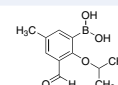
$C_{11}H_{13}BO_5$
MW: 236.03

**3,5-Diformyl-2-propoxyphenylboronic acid****684546**

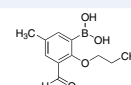
$C_{11}H_{13}BO_5$
MW: 236.03

**3-Formyl-2-isopropoxy-5-methylphenylboronic acid****567302**

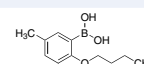
$C_{11}H_{15}BO_4$
MW: 222.05
[480424-52-0]

**3-Formyl-5-methyl-2-propoxyphenylboronic acid****661988**

$C_{11}H_{15}BO_4$
MW: 222.05
[480424-53-1]

**3-Bromo-2-butoxy-5-methylphenylboronic acid****596841**

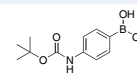
$C_{11}H_{16}BrO_3$
MW: 286.96
[870718-03-9]

**3-(N-Boc-amino)phenylboronic acid****578851**

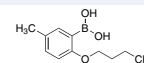
$C_{11}H_{16}BNO_4$
MW: 237.06
[380430-68-2]

**4-(N-Boc-amino)phenylboronic acid****565814**

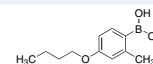
$C_{11}H_{16}BNO_4$
MW: 237.06
[380430-49-9]

**2-Butoxy-5-methylphenylboronic acid****565679**

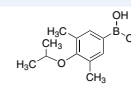
$C_{11}H_{17}BO_3$
MW: 208.06
[480438-72-0]

**4-Butoxy-2-methylphenylboronic acid****594628**

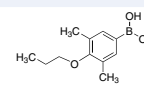
$C_{11}H_{17}BO_3$
MW: 208.06

**3,5-Dimethyl-4-isopropoxyphenylboronic acid****597732**

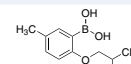
$C_{11}H_{17}BO_3$
MW: 208.06
[849062-16-4]

**3,5-Dimethyl-4-propoxyphenylboronic acid****597848**

$C_{11}H_{17}BO_3$
MW: 208.06
[357611-51-9]

**2-Isobutoxy-5-methylphenylboronic acid****651230**

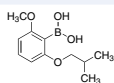
$C_{11}H_{17}BO_3$
MW: 208.06
[870778-94-2]



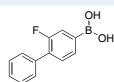
2-Isobutoxy-6-methoxyphenylboronic acid

NEW

662003
 $C_{11}H_{17}BO_4$
 MW: 224.06

**2-Fluoro-4-biphenylboronic acid**

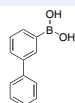
512109
 $C_{12}H_{10}BFO_2$
 MW: 216.02
 [178305-99-2]

**2-Biphenylboronic acid**

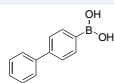
542202
 $C_{12}H_{11}BO_2$
 MW: 198.03
 [4688-76-0]

**3-Biphenylboronic acid**

542199
 $C_{12}H_{11}BO_2$
 MW: 198.03
 [5122-95-2]

**4-Biphenylboronic acid**

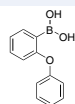
483451
 $C_{12}H_{11}BO_2$
 MW: 198.03
 [5122-94-1]

**Acenaphthene-5-boronic acid**

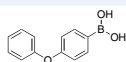
567477
 $C_{12}H_{11}BO_2$
 MW: 198.03
 [183158-33-0]

**2-Phenoxyphenylboronic acid**

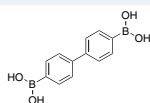
521485
 $C_{12}H_{11}BO_3$
 MW: 214.02
 [108238-09-1]

**4-Phenoxyphenylboronic acid**

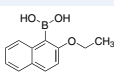
480142
 $C_{12}H_{11}BO_3$
 MW: 214.02
 [51067-38-0]

**4,4'-Biphenyldiboronic acid**

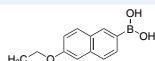
456799
 $C_{12}H_{12}B_2O_4$
 MW: 241.84
 [4151-80-8]

**2-Ethoxy-1-naphthaleneboronic acid**

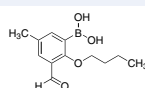
521884
 $C_{12}H_{13}BO_3$
 MW: 216.04
 [148345-64-6]

**6-Ethoxy-2-naphthaleneboronic acid**

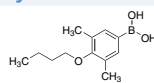
521906
 $C_{12}H_{13}BO_3$
 MW: 216.04
 [352525-98-5]

**2-Butoxy-3-formyl-5-methylphenylboronic acid**

567299
 $C_{12}H_{17}BO_4$
 MW: 236.07
 [480424-51-9]

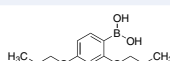
**4-Butoxy-3,5-dimethylphenylboronic acid**

597511
 $C_{12}H_{19}BO_3$
 MW: 222.09
 [845551-41-9]

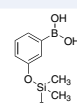
**2,4-Dipropoxyphenylboronic acid**

NEW

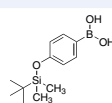
680672
 $C_{12}H_{19}BO_4$
 MW: 238.09
 [150145-25-8]

**3-(tert-Butyldimethylsilyloxy)phenylboronic acid**

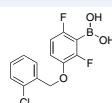
524034
 $C_{12}H_{21}BO_3Si$
 MW: 252.19
 [261621-12-9]

**4-(tert-Butyldimethylsilyloxy)phenylboronic acid**

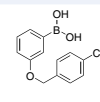
524042
 $C_{12}H_{21}BO_3Si$
 MW: 252.19
 [159191-56-7]

**2,6-Difluoro-3-(2'-chlorobenzoyloxy)phenylboronic acid**

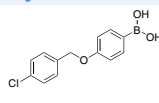
651362
 $C_{13}H_{10}BClF_2O_3$
 MW: 298.48
 [870778-99-7]

**3-(4'-Chlorobenzoyloxy)phenylboronic acid**

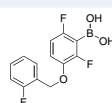
651036
 $C_{13}H_{10}BClO_3$
 MW: 260.48
 [870778-90-8]

**4-(4'-Chlorobenzoyloxy)phenylboronic acid**

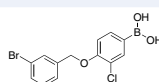
651044
 $C_{13}H_{10}BClO_3$
 MW: 260.48
 [870778-91-9]

**2,6-Difluoro-3-(2'-fluorobenzoyloxy)phenylboronic acid**

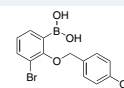
651346
 $C_{13}H_{10}BF_3O_3$
 MW: 282.02
 [836615-83-9]

**4-(3-Bromobenzoyloxy)-3-chlorophenylboronic acid**

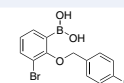
639524
 $C_{13}H_{11}BBrClO_3$
 MW: 341.39
 [849062-25-5]

**3-Bromo-2-(4'-chlorobenzoyloxy)phenylboronic acid**

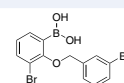
645842
 $C_{13}H_{11}BBrClO_3$
 MW: 341.39
 [849052-23-9]

**3-Bromo-2-(4'-fluorobenzoyloxy)phenylboronic acid**

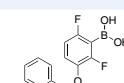
645761
 $C_{13}H_{11}BBrFO_3$
 MW: 324.94
 [849052-22-8]

**3-Bromo-2-(3'-bromobenzoyloxy)phenylboronic acid**

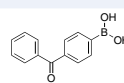
639567
 $C_{13}H_{11}BBr_2O_3$
 MW: 385.84
 [849062-27-7]

**3-Benzyloxy-2,6-difluorophenylboronic acid**

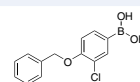
635707
 $C_{13}H_{11}BF_2O_3$
 MW: 264.03
 [870718-07-3]

**4-Benzoylphenylboronic acid**

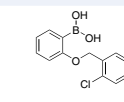
526010
 $C_{13}H_{11}BO_3$
 MW: 226.04
 [268218-94-6]

**4-Benzyloxy-3-chlorophenylboronic acid**

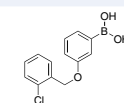
595616
 $C_{13}H_{12}BClO_3$
 MW: 262.50
 [845551-44-2]

**2-(2'-Chlorobenzoyloxy)phenylboronic acid**

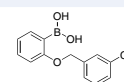
639443
 $C_{13}H_{12}BClO_3$
 MW: 262.50
 [870777-21-2]

**3-(2-Chlorobenzoyloxy)phenylboronic acid**

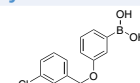
635731
 $C_{13}H_{12}BClO_3$
 MW: 262.50
 [845551-45-3]

**2-(3'-Chlorobenzoyloxy)phenylboronic acid**

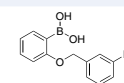
645249
 $C_{13}H_{12}BClO_3$
 MW: 262.50
 [849062-32-4]

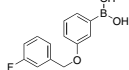
**3-(3'-Chlorobenzoyloxy)phenylboronic acid**

645265
 $C_{13}H_{12}BClO_3$
 MW: 262.50
 [849062-33-5]

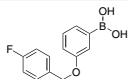
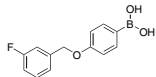
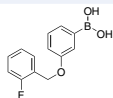
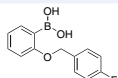
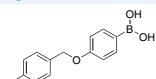
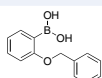
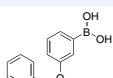
**2-(3'-Fluorobenzoyloxy)phenylboronic acid**

657492
 $C_{13}H_{12}BFO_3$
 MW: 246.04
 [871126-24-8]

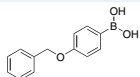
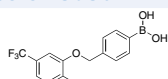
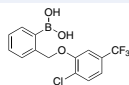
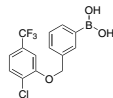


3-(3-Fluorobenzoyloxy)phenylboronic acid**661872**C₁₃H₁₂BO₃
MW: 246.04**3-(4'-Fluorobenzoyloxy)phenylboronic acid**

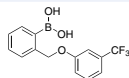
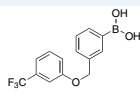
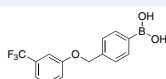
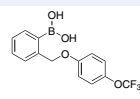
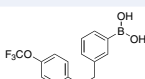
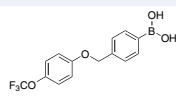
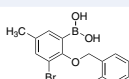
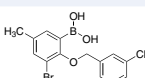
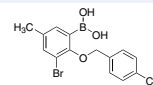
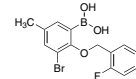
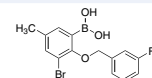
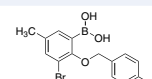
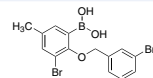
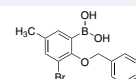
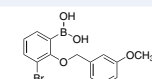
NEW

666777C₁₃H₁₂BO₃
MW: 246.04**4-(3-Fluorobenzoyloxy)phenylboronic acid****661864**C₁₃H₁₂BO₃
MW: 246.04**3-(2'-Fluorobenzoyloxy)phenylboronic acid****652113**C₁₃H₁₂BO₃
MW: 246.04
[849062-13-1]**2-(4'-Fluorobenzoyloxy)phenylboronic acid****652105**C₁₃H₁₂BO₃
MW: 246.04
[870779-01-4]**4-(4'-Fluorobenzoyloxy)phenylboronic acid****658073**C₁₃H₁₂BO₃
MW: 246.04
[871125-82-5]**2-Benzyloxyphenylboronic acid****521337**C₁₃H₁₃BO₃
MW: 228.05
[190661-29-1]**3-Benzyloxyphenylboronic acid****526339**C₁₃H₁₃BO₃
MW: 228.05
[156682-54-1]**4-Benzyloxyphenylboronic acid**

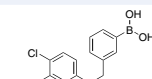
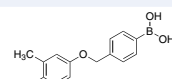
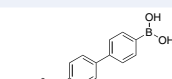
NEW

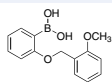
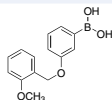
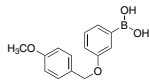
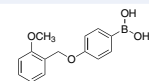
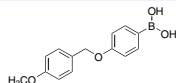
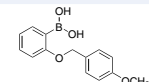
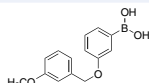
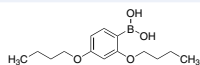
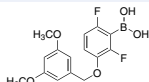
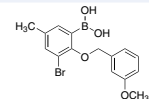
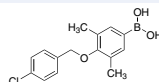
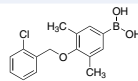
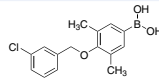
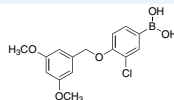
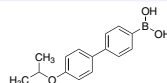
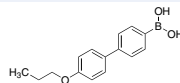
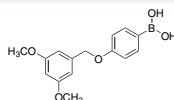
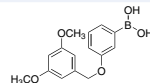
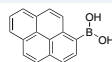
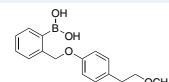
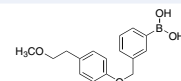
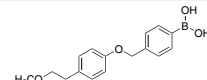
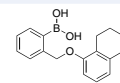
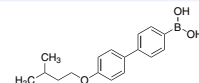
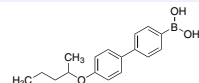
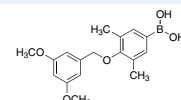
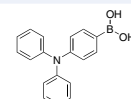
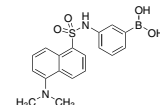
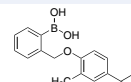
666920C₁₃H₁₃BO₃
MW: 228.05
[146631-00-7]**10-Bromoanthracene-9-boronic acid****595756**C₁₄H₁₀BBrO₂
MW: 300.94
[641144-16-3]**4-[(2'-Chloro-5'-(trifluoromethyl)-phenoxy)methyl]phenylboronic acid****652237**C₁₄H₁₁BClF₃O₃
MW: 330.49
[849062-05-1]**2-[(2'-Chloro-5'-(trifluoromethyl)-phenoxy)methyl]phenylboronic acid****652245**C₁₄H₁₁BClF₃O₃
MW: 330.49
[849062-11-9]**3-[(2'-Chloro-5'-(trifluoromethyl)-phenoxy)methyl]phenylboronic acid****657530**C₁₄H₁₁BClF₃O₃
MW: 330.49
[871126-25-9]**2-[(3'-(Trifluoromethyl)-phenoxy)methyl]phenylboronic acid**

NEW

666866C₁₄H₁₂BF₃O₃
MW: 296.05**3-(3'-(Trifluoromethyl)phenoxy)methyl-phenylboronic acid****651311**C₁₄H₁₂BF₃O₃
MW: 296.05
[870778-98-6]**4-(3'-(Trifluoromethyl)phenoxy)methyl-phenylboronic acid****651222**C₁₄H₁₂BF₃O₃
MW: 296.05
[849062-03-9]**2-[(4'-(Trifluoromethoxy)phenoxy)methyl]phenylboronic acid****652156**C₁₄H₁₂BF₃O₄
MW: 312.05
[849062-07-3]**3-[(4'-(Trifluoromethoxy)phenoxy)methyl]phenylboronic acid****652075**C₁₄H₁₂BF₃O₄
MW: 312.05
[849062-06-2]**4-[4'-(Trifluoromethoxy)phenoxy)methyl]phenylboronic acid****651303**C₁₄H₁₂BF₃O₄
MW: 312.05
[870778-97-5]**3-Bromo-2-(2'-chlorobenzoyloxy)-5-methylphenylboronic acid****645486**C₁₄H₁₃BBrClO₃
MW: 355.42
[849052-17-1]**3-Bromo-2-(3'-chlorobenzoyloxy)-5-methylphenylboronic acid****645494**C₁₄H₁₃BBrClO₃
MW: 355.42
[870778-83-9]**3-Bromo-2-(4'-chlorobenzoyloxy)-5-methylphenylboronic acid****645508**C₁₄H₁₃BBrClO₃
MW: 355.42
[849052-18-2]**3-Bromo-2-(2'-fluorobenzoyloxy)-5-methylphenylboronic acid****645400**C₁₄H₁₃BBrFO₃
MW: 338.96
[849062-18-6]**3-Bromo-2-(3'-fluorobenzoyloxy)-5-methylphenylboronic acid****645419**C₁₄H₁₃BBrFO₃
MW: 338.96
[849062-40-4]**3-Bromo-2-(4'-fluorobenzoyloxy)-5-methylphenylboronic acid****645427**C₁₄H₁₃BBrFO₃
MW: 338.96
[849062-41-5]**3-Bromo-2-(3'-bromobenzoyloxy)-5-methylphenylboronic acid****645478**C₁₄H₁₃BBr₂O₃
MW: 399.87
[849052-16-0]**2-Benzyloxy-3-bromo-5-methylphenylboronic acid****639435**C₁₄H₁₄BBrO₃
MW: 320.97
[870777-20-1]**3-Bromo-2-(3'-methoxybenzoyloxy)phenylboronic acid****645850**C₁₄H₁₄BBrO₄
MW: 336.97
[849052-24-0]**3-[(4-Chloro-3-methylphenoxy)methyl]phenylboronic acid**

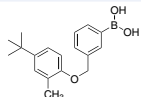
NEW

680605C₁₄H₁₄BClO₃
MW: 276.52**4-[(4-Chloro-3-methylphenoxy)methyl]phenylboronic acid****645931**C₁₄H₁₄BClO₃
MW: 276.52
[849052-25-1]**4-Ethoxybiphenyl-4'-boronic acid****567531**C₁₄H₁₅BO₃
MW: 242.08
[182344-29-2]

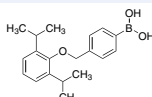
2-(2'-Methoxybenzyloxy)phenylboronic acid**657409**C₁₄H₁₅BO₄
MW: 258.08
[871125-76-7]**3-(2-Methoxybenzyloxy)phenylboronic acid****662089**C₁₄H₁₅BO₄
MW: 258.08**3-(4'-Methoxybenzyloxy)phenylboronic acid****666750**C₁₄H₁₅BO₄
MW: 258.08**4-(2'-Methoxybenzyloxy)phenylboronic acid****657387**C₁₄H₁₅BO₄
MW: 258.08
[871125-74-5]**4-(4'-Methoxybenzyloxy)phenylboronic acid****666769**C₁₄H₁₅BO₄
MW: 258.08**2-(4'-Methoxybenzyloxy)phenylboronic acid****657417**C₁₄H₁₅BO₄
MW: 258.08
[871125-77-8]**3-(3'-Methoxybenzyloxy)phenylboronic acid****657395**C₁₄H₁₅BO₄
MW: 258.08
[871125-75-6]**2,4-Dibutoxyphenylboronic acid****650994**C₁₄H₂₃BO₄
MW: 266.14
[870778-89-5]**2,6-Difluoro-3-(3',5'-dimethoxybenzyloxy)phenylboronic acid****651109**C₁₅H₁₅BF₂O₅
MW: 324.08
[849062-01-7]**3-Bromo-5-methyl-2-(3'-methoxybenzyloxy)phenylboronic acid****639478**C₁₅H₁₆BBrO₄
MW: 351.00
[849062-23-3]**4-(4'-Chlorobenzyloxy)-3,5-dimethylphenylboronic acid****645354**C₁₅H₁₆BClO₃
MW: 290.55
[849062-38-0]**4-(2'-Chlorobenzyloxy)-3,5-dimethylphenylboronic acid****645435**C₁₅H₁₆BClO₃
MW: 290.55
[849052-15-9]**4-(3'-Chlorobenzyloxy)-3,5-dimethylphenylboronic acid****645443**C₁₅H₁₆BClO₃
MW: 290.55
[849062-21-1]**3-Chloro-4-(3',5'-dimethoxybenzyloxy)phenylboronic acid****639516**C₁₅H₁₆BClO₅
MW: 322.55
[849062-24-4]**4-(4'-Isopropoxyphenyl)phenylboronic acid****595179**C₁₅H₁₇BO₃
MW: 256.10
[870717-98-9]**4-(4'-Propoxyphenyl)phenylboronic acid****595284**C₁₅H₁₇BO₃
MW: 256.10
[849062-20-0]**4-(3,5-Dimethoxybenzyloxy)phenylboronic acid****635758**C₁₅H₁₇BO₅
MW: 288.10
[870718-08-4]**3-(3,5-Dimethoxybenzyloxy)phenylboronic acid****635766**C₁₅H₁₇BO₅
MW: 288.10
[870718-09-5]**Pyrene-1-boronic acid****542873**C₁₆H₁₁BO₂
MW: 246.07
[164461-18-1]**2-[(4'-(2-Methoxyethyl)phenoxy)methyl]phenylboronic acid****658065**C₁₆H₁₉BO₄
MW: 286.13
[871126-29-3]**3-[(4'-(2-Methoxyethyl)phenoxy)methyl]phenylboronic acid****657506**C₁₆H₁₉BO₄
MW: 286.13
[871126-26-0]**4-[(4'-(2-Methoxyethyl)phenoxy)methyl]phenylboronic acid****652083**C₁₆H₁₉BO₄
MW: 286.13
[870779-00-3]**2-[(1',2',3',4'-Tetrahydro-5-naphthoxy)methyl]phenylboronic acid****652229**C₁₇H₁₉BO₃
MW: 282.14
[849062-10-8]**4-(4-Isopentyloxyphenyl)phenylboronic acid****662070**C₁₇H₂₁BO₃
MW: 284.16**4-(4'-(2-Pentyloxy)phenyl)phenylboronic acid****666815**C₁₇H₂₁BO₃
MW: 284.16**3,5-Dimethyl-4-(3',5'-dimethoxybenzyloxy)phenylboronic acid****652121**C₁₇H₂₁BO₅
MW: 316.16**4-(Diphenylamino)phenylboronic acid****647292**C₁₈H₁₆BN₂O₂
MW: 289.14
[201802-67-7]**3-(Dansylamino)phenylboronic acid****D0520**C₁₈H₁₉BN₂O₄S
MW: 370.23
[75806-94-9]**2-[(4-tert-Butyl-2-methylphenoxy)methyl]phenylboronic acid****662046**C₁₈H₂₃BO₃
MW: 298.18

3-((4'-*tert*-Butyl-2'-methylphenoxy)-methyl)phenylboronic acid NEW

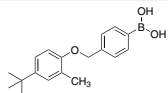
666858
 $C_{18}H_{23}BO_3$
 MW: 298.18

**4-[(2',6'-Diisopropylphenoxy)methyl]phenylboronic acid**

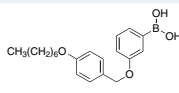
662038
 $C_{19}H_{25}BO_3$
 MW: 312.21

**4-((4'-*tert*-Butyl-2-methylphenoxy)-methyl)phenylboronic acid** NEW

680591
 $C_{18}H_{23}BO_3$
 MW: 298.18

**3-(4'-Heptyloxyphenoxy)methyl)phenylboronic acid**

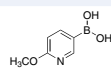
651214
 $C_{20}H_{27}BO_4$
 MW: 342.24
 [870778-93-1]

**N-Heteroarylboronic acids****2-Bromopyridine-5-boronic acid** NEW

666556
 $C_5H_5BBrNO_2$
 MW: 201.81
 [223463-14-7]

**2-Methoxypyridine-5-boronic acid**

637610
 $C_6H_8BNO_3$
 MW: 152.94
 [163105-89-3]

**N-Boc-pyrrole-2-boronic acid**

15047
 $C_9H_{14}BNO_4$
 MW: 211.02
 [135884-31-0]

**2-Chloropyridine-5-boronic acid**

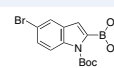
637386
 $C_5H_5BClNO_2$
 MW: 157.36
 [444120-91-6]

**2-Methoxypyridine-3-boronic acid** NEW

684589
 $C_6H_8BNO_3$
 MW: 152.94
 [163105-90-6]

**1-Boc-5-bromindole-2-boronic acid**

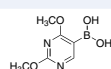
637394
 $C_{13}H_{15}BBrNO_4$
 MW: 339.98
 [475102-13-7]

**2-Chloropyridine-4-boronic acid** NEW

666513
 $C_5H_5BClNO_2$
 MW: 157.36
 [458532-96-2]

**2,4-Dimethoxypyrimidine-5-boronic acid** NEW

666491
 $C_6H_9BN_2O_4$
 MW: 183.96
 [89641-18-9]

**1-(Phenylsulfonyl)indole-2-boronic acid**

563862
 $C_{14}H_{12}BNO_4S$
 MW: 301.13
 [342404-46-0]

**2-Fluoropyridine-5-boronic acid**

639184
 $C_5H_5BFNO_2$
 MW: 140.91
 [351019-18-6]

**5-Indoleboronic acid**

666467
 $C_8H_8BNO_2$
 MW: 160.97
 [144104-59-6]

**1-(Phenylsulfonyl)indole-3-boronic acid**

563870
 $C_{14}H_{12}BNO_4S$
 MW: 301.13
 [129271-98-3]

**3-Pyridineboronic acid**

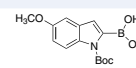
512125
 $C_5H_6BNO_2$
 MW: 122.92
 [1692-25-7]

**6-Indoleboronic acid** NEW

666459
 $C_8H_8BNO_2$
 MW: 160.97
 [147621-18-9]

**1-Boc-5-methoxyindole-2-boronic acid** NEW

680516
 $C_{14}H_{18}BNO_5$
 MW: 291.11
 [290331-71-4]

**4-Pyridineboronic acid**

634492
 $C_5H_6BNO_2$
 MW: 122.92
 [1692-15-5]

**8-Quinolineboronic acid**

542865
 $C_9H_8BNO_2$
 MW: 172.98
 [86-58-8]

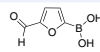


O-Heteroarylboronic acids

2-Furanboronic acid

464910C₄H₅BO₃
MW: 111.89
[13331-23-2]

5-Formylfuran-2-boronic acid

512346C₅H₅BO₄
MW: 139.90
[27329-70-0]

1,4-Benzodioxane-6-boronic acid

635995C₈H₉BO₄
MW: 179.97
[164014-95-3]

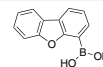
3-Furanboronic acid

512168C₄H₅BO₃
MW: 111.89
[55552-70-0]

2-Benzofuranboronic acid

499943C₈H₇BO₃
MW: 161.95
[98437-24-2]

4-Dibenzofuranboronic acid

499951C₁₂H₉BO₃
MW: 212.01
[100124-06-9]

S-Heteroarylboronic acids

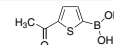
5-Bromothiophene-2-boronic acid

557684C₄H₄BBrO₂S
MW: 206.85
[162607-17-2]

2-Formylthiophene-3-boronic acid

499900C₅H₅BO₃S
MW: 155.97
[4347-31-3]

5-Acetylthiophene-2-boronic acid

499927C₆H₇BO₃S
MW: 169.99
[206551-43-1]

5-Chlorothiophene-2-boronic acid

499935C₄H₄BClO₂S
MW: 162.40
[162607-18-3]

3-Formylthiophene-2-boronic acid

499919C₅H₅BO₃S
MW: 155.97
[17303-83-2]

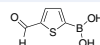
2-Benzo[b]thiopheneboronic acid

499978C₈H₇BO₂S
MW: 178.02
[98437-23-1]

2-Thiopheneboronic acid

436836C₄H₅BO₂S
MW: 127.96
[6165-68-0]

5-Formylthiophene-2-boronic acid

514055C₅H₅BO₃S
MW: 155.97
[4347-33-5]

3-Benzo[b]thiopheneboronic acid

512117C₈H₇BO₂S
MW: 178.02
[113893-08-6]

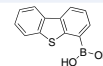
3-Thiopheneboronic acid

436844C₄H₅BO₂S
MW: 127.96
[6165-69-1]

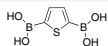
4-Methylthiophene-3-boronic acid

542393C₅H₇BO₂S
MW: 141.98
[177735-11-4]

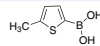
4-Dibenzothiopheneboronic acid

499986C₁₂H₉BO₂S
MW: 228.07
[108847-20-7]

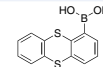
2,5-Thiophenediboronic acid

470317C₄H₆B₂O₄S
MW: 171.78
[26076-46-0]

5-Methylthiophene-2-boronic acid

512192C₅H₇BO₂S
MW: 141.98
[162607-20-7]

1-Thianthreneboronic acid

512214C₁₂H₉BO₂S₂
MW: 260.14
[108847-76-3]

Boronate Esters

Alkylboronate esters

Dichloromethyldiisopropoxyborane

324485

$C_7H_{15}BCl_2O_2$
MW: 212.91
[62260-99-5]



Diisopropoxymethylborane

283681

$C_7H_{17}BO_2$
MW: 144.02
[86595-27-9]



(3R)-(+)-Tetrahydrofuranylboronic acid diethanolamine ester

376584

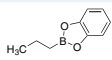
$C_8H_{16}BNO_3$
MW: 185.03
[100858-40-0]



1-Propylboronic acid catechol ester

464511

$C_9H_{11}BO_2$
MW: 161.99
[40218-49-3]



Cyclopropylboronic acid pinacol ester

659851

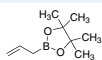
$C_9H_{17}BO_2$
MW: 168.04
[126689-01-8]



Allylboronic acid pinacol ester

324647

$C_9H_{17}BO_2$
MW: 168.04
[72824-04-5]



Alkenylboronate esters

Vinylboronic acid pinacol ester

633348

$C_8H_{15}BO_2$
MW: 154.01
[75927-49-0]



Vinylboronic acid 2-methyl-2,4-pentanediol ester

673641

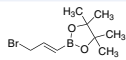
$C_8H_{15}BO_2$
MW: 154.01
[4627-10-5]



trans-3-Bromo-1-propen-1-ylboronic acid pinacol ester

643785

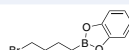
$C_9H_{16}BBrO_2$
MW: 246.94
[870777-31-4]



4-Bromo-1-butylboronic acid catechol ester

513687

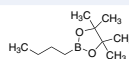
$C_{10}H_{12}BBrO_2$
MW: 254.92
[142172-51-8]



n-Butylboronic acid pinacol ester

680923

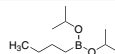
$C_{10}H_{21}BO_2$
MW: 184.08
[69190-62-1]



Butyldiisopropoxyborane

277975

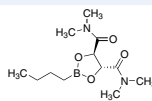
$C_{10}H_{23}BO_2$
MW: 186.10
[86595-32-6]



Butylboronic acid N,N,N'-tetramethyl-L-tartaric acid diamide ester

518379

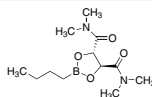
$C_{12}H_{23}BN_2O_4$
MW: 270.13
[161344-85-0]



Butylboronic acid N,N,N'-tetramethyl-D-tartaric acid diamide ester

594806

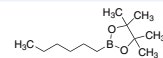
$C_{12}H_{23}BN_2O_4$
MW: 270.13
[161344-84-9]



Hexylboronic acid pinacol ester

673668

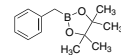
$C_{12}H_{25}BO_2$
MW: 212.14
[86308-26-1]



Benzylboronic acid pinacol ester

659207

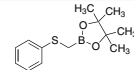
$C_{13}H_{19}BO_2$
MW: 218.10
[87100-28-5]



Phenylsulfanylmethylboronic acid pinacol ester

568155

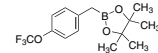
$C_{13}H_{19}BO_2S$
MW: 250.16
[66080-23-7]



4-(Trifluoromethoxy)benzylboronic acid pinacol ester

662879

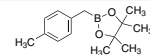
$C_{14}H_{18}BF_3O_3$
MW: 302.10
[475250-46-5]



4-Methylbenzylboronic acid pinacol ester

663298

$C_{14}H_{21}BO_2$
MW: 232.13
[356570-52-0]



Vinylboronic acid pinacol ester

633348

$C_8H_{15}BO_2$
MW: 154.01
[75927-49-0]



Vinylboronic acid 2-methyl-2,4-pentanediol ester

673641

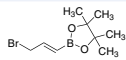
$C_8H_{15}BO_2$
MW: 154.01
[4627-10-5]



trans-3-Bromo-1-propen-1-ylboronic acid pinacol ester

643785

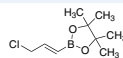
$C_9H_{16}BBrO_2$
MW: 246.94
[870777-31-4]



trans-3-Chloro-1-propen-1-ylboronic acid pinacol ester

680249

$C_9H_{16}BClO_2$
MW: 202.49



Isopropenylboronic acid pinacol ester

663212

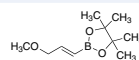
$C_9H_{17}BO_2$
MW: 168.04
[126726-62-3]



trans-3-Methoxy-1-propenylboronic acid pinacol ester

592935

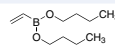
$C_{10}H_{19}BO_3$
MW: 198.07
[165059-42-7]



Vinylboronic acid dibutyl ester

568147

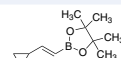
$C_{10}H_{21}BO_2$
MW: 184.08
[6336-45-4]



trans-2-Cyclopropylvinylboronic acid pinacol ester

653942

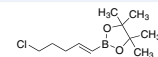
$C_{11}H_{19}BO_2$
MW: 194.08
[849061-99-0]



trans-5-Chloro-1-penten-1-ylboronic acid pinacol ester

652067

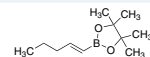
$C_{11}H_{20}BClO_2$
MW: 230.54
[154820-95-8]

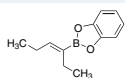
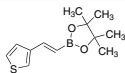
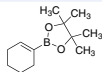
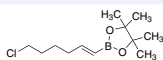
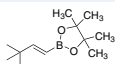
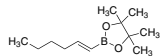
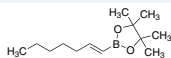
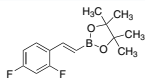
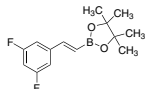
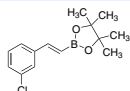
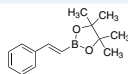
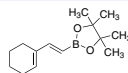
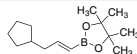
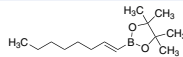
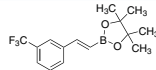
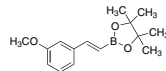
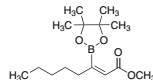
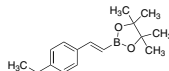
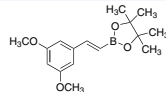
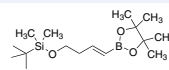
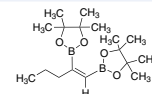
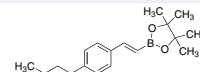
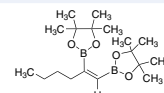
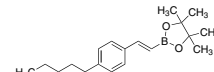
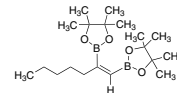
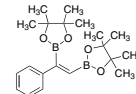
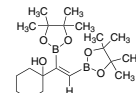
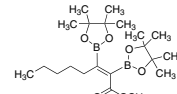
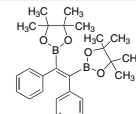


trans-1-Penten-1-ylboronic acid pinacol ester

665169

$C_{11}H_{21}BO_2$
MW: 196.09
[161395-96-6]



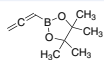
(Z)-3-Hexenyl-3-boronic acid catechol ester**575526**C₁₂H₁₅BO₂
MW: 202.06
[37490-28-1]**trans-2-(Thiophen-3-yl)vinylboronic acid pinacol ester** NEW**680257**C₁₂H₁₇BO₂S
MW: 236.14**1-Cyclohexen-1-ylboronic acid pinacol ester****650277**C₁₂H₂₁BO₂
MW: 208.10
[141091-37-4]**trans-6-Chloro-1-hexen-1-ylboronic acid pinacol ester****653950**C₁₂H₂₂BClO₂
MW: 244.57
[197313-32-9]**trans-(3,3-Dimethylbuten-1-yl)boronic acid pinacol ester****667277**C₁₂H₂₃BO₂
MW: 210.12**trans-1-Hexen-1-ylboronic acid pinacol ester****663743**C₁₂H₂₃BO₂
MW: 210.12
[126688-97-9]**trans-1-Hepten-1-ylboronic acid pinacol ester****662992**C₁₃H₂₅BO₂
MW: 224.15
[169339-75-7]**trans-2-(2,4-Difluorophenyl)vinylboronic acid pinacol ester****664871**C₁₄H₁₇BF₂O₂
MW: 266.09
[736987-78-3]**trans-2-(3,5-Difluorophenyl)vinylboronic acid pinacol ester** NEW**669199**C₁₄H₁₇BF₂O₂
MW: 266.09**trans-2-(3-Chlorophenyl)vinylboronic acid pinacol ester****665533**C₁₄H₁₈BClO₂
MW: 264.56
[871125-84-7]**1-Phenylvinylboronic acid pinacol ester****659193**C₁₄H₁₉BO₂
MW: 230.11
[78782-27-1]**trans-2-Phenylvinylboronic acid pinacol ester****570214**C₁₄H₁₉BO₂
MW: 230.11
[83947-56-2]**trans-2-(1-Cyclohexenyl)vinylboronic acid pinacol ester** NEW**676373**C₁₄H₂₃BO₂
MW: 234.14**trans-3-(Cyclopentyl)propenylboronic acid pinacol ester****667013**C₁₄H₂₅BO₂
MW: 236.16**trans-1-Octen-1-ylboronic acid pinacol ester****663050**C₁₄H₂₇BO₂
MW: 238.17
[83947-55-1]**trans-2-(3-(Trifluoromethyl)phenyl)vinylboronic acid pinacol ester** NEW**677639**C₁₅H₁₈BF₃O₂
MW: 298.11**trans-2-(3-Methoxyphenyl)vinylboronic acid pinacol ester****656909**C₁₅H₂₁BO₃
MW: 260.14
[871125-68-7]**Methyl (Z)-oct-2-enoate-3-boronic acid pinacol ester****540625**C₁₅H₂₇BO₄
MW: 282.18
[352534-74-8]**trans-2-(4-Ethylphenyl)vinylboronic acid pinacol ester****662798**C₁₆H₂₃BO₂
MW: 258.16
[870717-91-2]**trans-2-(3,5-Dimethoxyphenyl)vinylboronic acid pinacol ester** NEW**676160**C₁₆H₂₃BO₄
MW: 290.16**trans-4-(tert-Butyldimethylsiloxy)-1-buten-1-ylboronic acid****588849**C₁₆H₃₃BO₂Si
MW: 312.33
[480425-30-7]**(E)-1-Pentene-1,2-diboronic acid bis(pinacol) ester****529907**C₁₇H₃₂B₂O₄
MW: 322.06
[307531-75-5]**2-(4-Butylphenyl)vinylboronic acid pinacol ester****592870**C₁₈H₂₇BO₂
MW: 286.22**(E)-1-Hexene-1,2-diboronic acid bis(pinacol) ester****527602**C₁₈H₃₄B₂O₄
MW: 336.08
[312693-52-0]**trans-2-(4-Pentylphenyl)vinylboronic acid pinacol ester****592986**C₁₉H₂₉BO₂
MW: 300.24**(E)-1-Heptene-1,2-diboronic acid bis(pinacol) ester****529893**C₁₉H₃₆B₂O₄
MW: 350.11
[307531-74-4]**(E)-1-Phenyl-1,2-ethylenediboronic acid bis(pinacol) ester****527572**C₂₀H₃₀B₂O₄
MW: 356.07
[173603-23-1]**(E)-1-(1-Hydroxycyclohexyl)ethene-1,2-diboronic acid bis(pinacol) ester****527610**C₂₀H₃₆B₂O₅
MW: 378.12
[264144-69-6]**Methyl (Z)-oct-2-enoate-2,3-diboronic acid bis(pinacol) ester****527599**C₂₁H₃₈B₂O₆
MW: 408.14
[312693-51-9]**(Z)-1,2-Diphenyl-1,2-ethylenediboronic acid bis(pinacol) ester****527580**C₂₆H₃₄B₂O₄
MW: 432.17
[221006-76-4]

Allenyl- and Alkynylboronate esters

Allenylboronic acid pinacol ester

NEW

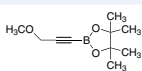
678554
 $C_9H_{15}BO_2$
 MW: 166.03



3-Methoxy-1-propyn-1-ylboronic acid pinacol ester

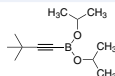
NEW

674710
 $C_{10}H_{17}BO_3$
 MW: 196.05



(2-tert-Butyl-1-ethynyl)diisopropoxyborane

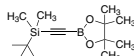
639192
 $C_{12}H_{23}BO_2$
 MW: 210.12
 [121021-24-7]



2-((tert-Butyldimethylsilyl)-ethynyl)boronic acid pinacol ester

NEW

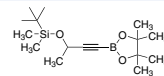
676659
 $C_{14}H_{27}BO_2Si$
 MW: 266.26



3-(tert-Butyldimethylsilyloxy)-1-butyne-1-ylboronic acid pinacol ester

NEW

674729
 $C_{16}H_{31}BO_3Si$
 MW: 310.31



Arylboronate esters

Phenylboronic acid 1,3-propanediol ester

341339
 $C_9H_{11}BO_2$
 MW: 161.99
 [4406-77-3]



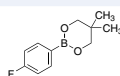
2-Cyanophenylboronic acid 1,3-propanediol ester

653934
 $C_{10}H_{10}BNO_2$
 MW: 187.00
 [172732-52-4]



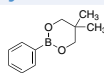
4-Fluorophenylboronic acid neopentylglycol ester

632686
 $C_{11}H_{14}BF_2O_2$
 MW: 208.04
 [225916-39-2]



Phenylboronic acid neopentylglycol ester

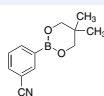
632678
 $C_{11}H_{15}BO_2$
 MW: 190.05
 [5123-13-7]



(3-Cyanophenyl)boronic acid neopentylglycol ester

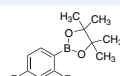
NEW

651168
 $C_{12}H_{14}BNO_2$
 MW: 215.06
 [214360-45-9]



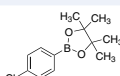
2,4-Difluorophenylboronic acid pinacol ester

632694
 $C_{12}H_{15}BF_2O_2$
 MW: 240.05
 [288101-48-4]



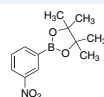
4-Chlorophenylboronic acid pinacol ester

632724
 $C_{12}H_{16}BClO_2$
 MW: 238.52
 [195062-61-4]



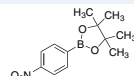
3-Nitrophenylboronic acid pinacol ester

632708
 $C_{12}H_{16}BNO_4$
 MW: 249.07
 [68716-48-3]



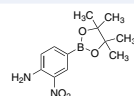
4-Nitrophenylboronic acid pinacol ester

643890
 $C_{12}H_{16}BNO_4$
 MW: 249.07
 [171364-83-3]



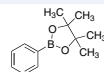
4-Amino-3-nitrophenylboronic acid pinacol ester

651613
 $C_{12}H_{17}BN_2O_4$
 MW: 264.09
 [833486-94-5]



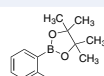
Phenylboronic acid pinacol ester

647098
 $C_{12}H_{17}BO_2$
 MW: 204.07
 [24388-23-6]



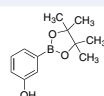
2-Hydroxyphenylboronic acid pinacol ester

522554
 $C_{12}H_{17}BO_3$
 MW: 220.07
 [269409-97-4]



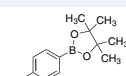
3-Hydroxyphenylboronic acid pinacol ester

522562
 $C_{12}H_{17}BO_3$
 MW: 220.07
 [214360-76-6]



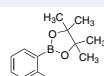
4-Hydroxyphenylboronic acid pinacol ester

522570
 $C_{12}H_{17}BO_3$
 MW: 220.07
 [269409-70-3]



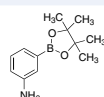
2-Aminophenylboronic acid pinacol ester

576557
 $C_{12}H_{18}BNO_2$
 MW: 219.09
 [191171-55-8]



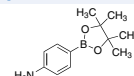
3-Aminophenylboronic acid pinacol ester

574686
 $C_{12}H_{18}BNO_2$
 MW: 219.09
 [210907-84-9]



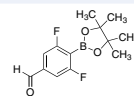
4-Aminophenylboronic acid pinacol ester

518751
 $C_{12}H_{18}BNO_2$
 MW: 219.09
 [214360-73-3]



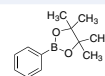
2,6-Difluoro-4-formylphenylboronic acid pinacol ester

663514
 $C_{13}H_{15}BF_2O_3$
 MW: 268.06
 [870717-92-3]



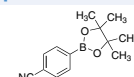
3-Cyanophenylboronic acid pinacol ester

578401
 $C_{13}H_{16}BNO_2$
 MW: 229.08
 [214360-46-0]



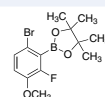
4-Cyanophenylboronic acid pinacol ester

527556
 $C_{13}H_{16}BNO_2$
 MW: 229.08
 [171364-82-2]



6-Bromo-2-fluoro-3-methoxyphenylboronic acid pinacol ester

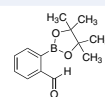
667560
 $C_{13}H_{17}BBrFO_3$
 MW: 330.99



2-Formylphenylboronic acid pinacol ester

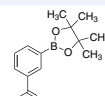
NEW

683817
 $C_{13}H_{17}BO_3$
 MW: 232.08
 [380151-85-9]



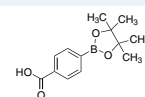
3-Carboxyphenylboronic acid pinacol ester

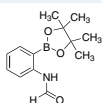
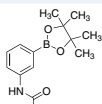
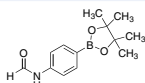
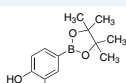
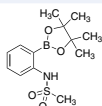
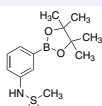
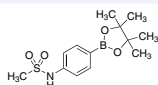
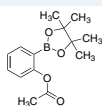
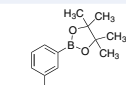
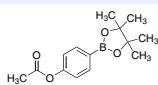
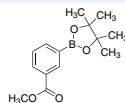
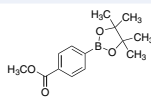
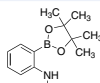
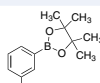
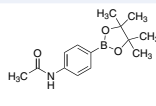
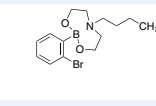
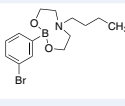
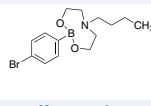
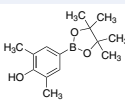
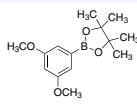
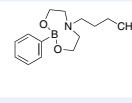
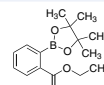
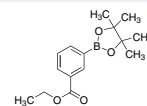
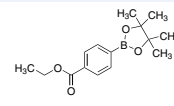
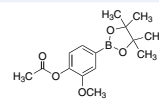
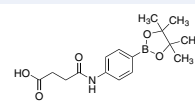
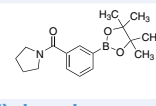
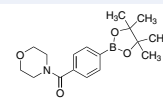
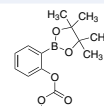
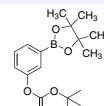
574694
 $C_{13}H_{17}BO_4$
 MW: 248.08
 [269409-73-6]



4-Carboxyphenylboronic acid pinacol ester

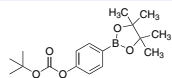
513490
 $C_{13}H_{17}BO_4$
 MW: 248.08
 [180516-87-4]



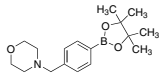
2-(Formylamino)phenylboronic acid pinacol ester**595098**C₁₃H₁₈BNO₃
MW: 247.10
[480425-36-3]**3-(Formylamino)phenylboronic acid pinacol ester****595209**C₁₃H₁₈BNO₃
MW: 247.10
[480425-37-4]**4-(Formylamino)phenylboronic acid pinacol ester****578746**C₁₃H₁₈BNO₃
MW: 247.10
[480424-94-0]**4-Hydroxy-3-methoxyphenylboronic acid pinacol ester****518786**C₁₃H₁₉BO₄
MW: 250.10
[269410-22-2]**2-(Methanesulfonylamino)phenylboronic acid pinacol ester****636045**C₁₃H₂₀BNO₄S
MW: 297.18
[380430-60-4]**3-(Methanesulfonylamino)phenylboronic acid pinacol ester****636037**C₁₃H₂₀BNO₄S
MW: 297.18
[305448-92-4]**4-(Methanesulfonylamino)phenylboronic acid pinacol ester****632716**C₁₃H₂₀BNO₄S
MW: 297.18
[616880-14-9]**2-Acetoxyphenylboronic acid pinacol ester****570184**C₁₄H₁₉BO₄
MW: 262.11
[480424-68-8]**3-Acetoxyphenylboronic acid pinacol ester****570192**C₁₄H₁₉BO₄
MW: 262.11
[480424-69-9]**4-Acetoxyphenylboronic acid pinacol ester****570206**C₁₄H₁₉BO₄
MW: 262.11
[480424-70-2]**3-Methoxycarbonylphenylboronic acid pinacol ester****594652**C₁₄H₁₉BO₄
MW: 262.11
[480425-35-2]**4-Methoxycarbonylphenylboronic acid pinacol ester****594768**C₁₄H₁₉BO₄
MW: 262.11
[171364-80-0]**2-Acetamidophenylboronic acid pinacol ester****594989**C₁₄H₂₀BNO₃
MW: 261.12
[380430-61-5]**3-Acetamidophenylboronic acid pinacol ester****578088**C₁₄H₂₀BNO₃
MW: 261.12
[480424-93-9]**4-Acetamidophenylboronic acid pinacol ester****518778**C₁₄H₂₀BNO₃
MW: 261.12
[214360-60-8]**2-Bromophenylboronic acid N-butyl-diethanolamine ester****684570**C₁₄H₂₁BBrNO₂
MW: 326.04**3-Bromophenylboronic acid N-butyl-diethanolamine ester****680486**C₁₄H₂₁BBrNO₂
MW: 326.04**4-Bromophenylboronic acid N-butyl-diethanolamine ester****680494**C₁₄H₂₁BBrNO₂
MW: 326.04**4-Hydroxy-3,5-dimethylphenylboronic acid pinacol ester****518794**C₁₄H₂₁BO₃
MW: 248.13
[269410-25-5]**3,5-Dimethoxyphenylboronic acid pinacol ester****633909**C₁₄H₂₁BO₄
MW: 264.13
[365564-07-4]**Phenylboronic acid N-butyl-diethanolamine ester****680478**C₁₄H₂₂BNO₂
MW: 247.14**2-Ethoxycarbonylphenylboronic acid pinacol ester****570168**C₁₅H₂₁BO₄
MW: 276.14
[269409-99-6]**3-Ethoxycarbonylphenylboronic acid pinacol ester****527548**C₁₅H₂₁BO₄
MW: 276.14
[269410-00-6]**4-Ethoxycarbonylphenylboronic acid pinacol ester****527564**C₁₅H₂₁BO₄
MW: 276.14
[195062-62-5]**4-Acetoxy-3-methoxyphenylboronic acid pinacol ester****594474**C₁₅H₂₁BO₅
MW: 292.14**4-(Succinylamino)phenylboronic acid pinacol ester****578762**C₁₆H₂₂BNO₅
MW: 319.16
[480424-98-4]**3-(Pyrrolidine-1-carbonyl)phenylboronic acid pinacol ester****680222**C₁₇H₂₄BNO₃
MW: 301.19**4-(Morpholine-4-carbonyl)phenylboronic acid pinacol ester****677809**C₁₇H₂₄BNO₄
MW: 317.19
[656239-38-2]**2-(tert-Butoxycarbonyloxy)phenylboronic acid pinacol ester****570222**C₁₇H₂₅BO₅
MW: 320.19
[480424-71-3]**3-(tert-Butoxycarbonyloxy)phenylboronic acid pinacol ester****565822**C₁₇H₂₅BO₅
MW: 320.19
[480438-74-2]

4-(tert-Butoxycarbonyloxy)phenylboronic acid pinacol ester

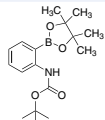
565946
 $C_{17}H_{25}BO_5$
 MW: 320.19
 [480438-75-3]

**4-(4-Morpholinomethyl)phenylboronic acid pinacol ester**

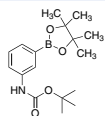
680230
 $C_{17}H_{26}BNO_3$
 MW: 303.20
 [364794-79-6]

**2-(N-Boc-amino)phenylboronic acid pinacol ester**

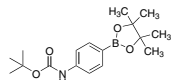
594865
 $C_{17}H_{26}BNO_4$
 MW: 319.20
 [159624-15-4]

**3-(N-Boc-amino)phenylboronic acid pinacol ester**

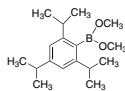
578843
 $C_{17}H_{26}BNO_4$
 MW: 319.20
 [330793-09-4]

**4-(N-Boc-amino)phenylboronic acid pinacol ester**

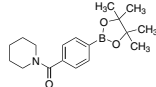
562351
 $C_{17}H_{26}BNO_4$
 MW: 319.20
 [330793-01-6]

**2,4,6-Triisopropylphenylboronic acid dimethyl ester**

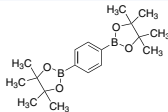
682683
 $C_{17}H_{29}BO_2$
 MW: 276.22
 [145434-22-6]

**4-(Piperidine-1-carbonyl)phenylboronic acid pinacol ester**

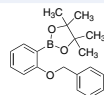
677787
 $C_{18}H_{26}BNO_3$
 MW: 315.21

**1,4-Benzenediboronic acid bis(pinacol) ester**

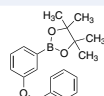
663816
 $C_{18}H_{28}B_2O_4$
 MW: 330.03
 [99770-93-1]

**2-(Benzyloxy)phenylboronic acid pinacol ester**

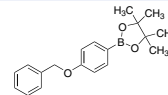
594245
 $C_{19}H_{23}BO_3$
 MW: 310.20

**3-(Benzyloxy)phenylboronic acid pinacol ester**

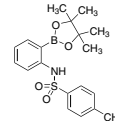
594369
 $C_{19}H_{23}BO_3$
 MW: 310.20

**4-(Benzyloxy)phenylboronic acid pinacol ester**

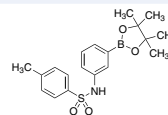
594148
 $C_{19}H_{23}BO_3$
 MW: 310.20

**2-(p-Toluenesulfonylamino)phenylboronic acid pinacol ester**

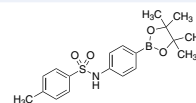
636290
 $C_{19}H_{24}BNO_4S$
 MW: 373.27
 [796061-07-9]

**3-(p-Toluenesulfonylamino)phenylboronic acid pinacol ester**

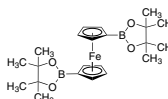
636312
 $C_{19}H_{24}BNO_4S$
 MW: 373.27
 [796061-08-0]

**4-(p-Toluenesulfonylamino)phenylboronic acid pinacol ester**

636304
 $C_{19}H_{24}BNO_4S$
 MW: 373.27
 [674776-54-6]

**1,1'-Ferrocenediboronic acid bis(pinacol) ester**

643750
 $C_{22}H_{32}B_2O_4Fe$
 MW: 437.95
 [737776-93-1]

**N-Heteroarylboronate esters****3-Pyridineboronic acid 1,3-propanediol ester**

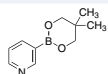
631566
 $C_8H_{10}BNO_2$
 MW: 162.98
 [131534-65-1]

**4-Pyrazoleboronic acid pinacol ester**

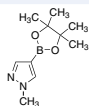
525057
 $C_9H_{15}BN_2O_2$
 MW: 194.04
 [269410-08-4]

**3-Pyridineboronic acid neopentylglycol ester**

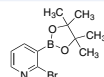
642622
 $C_{10}H_{14}BNO_2$
 MW: 191.03
 [845885-86-1]

**1-Methylpyrazole-4-boronic acid pinacol ester**

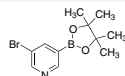
595314
 $C_{10}H_{17}BN_2O_2$
 MW: 208.07
 [761446-44-0]

**2-Bromopyridine-3-boronic acid pinacol ester**

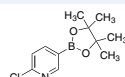
654264
 $C_{11}H_{15}BBrNO_2$
 MW: 283.96
 [452972-12-2]

**5-Bromopyridine-3-boronic acid pinacol ester**

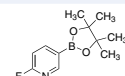
674567
 $C_{11}H_{15}BBrNO_2$
 MW: 283.96
 [452972-13-3]

**6-Chloropyridine-3-boronic acid pinacol ester**

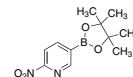
659843
 $C_{11}H_{15}BClNO_2$
 MW: 239.51
 [444120-94-9]

**2-Fluoropyridine-5-boronic acid pinacol ester**

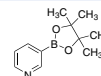
592544
 $C_{11}H_{15}BFNO_2$
 MW: 223.05
 [444120-95-0]

**2-Nitropyridine-5-boronic acid pinacol ester**

674958
 $C_{11}H_{15}BN_2O_4$
 MW: 250.06

**3-Pyridineboronic acid pinacol ester**

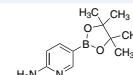
576565
 $C_{11}H_{16}BNO_2$
 MW: 205.06
 [329214-79-1]

**4-Pyridineboronic acid pinacol ester**

578770
 $C_{11}H_{16}BNO_2$
 MW: 205.06
 [181219-01-2]

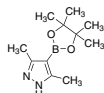
**2-Aminopyridine-5-boronic acid pinacol ester**

640379
 $C_{11}H_{17}BN_2O_2$
 MW: 220.08
 [827614-64-2]

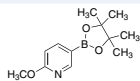


3,5-Dimethylpyrazole-4-boronic acid pinacol ester

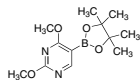
636010
 $C_{11}H_{19}BN_2O_2$
 MW: 222.09
 [857530-80-4]

**2-Methoxypyridine-5-boronic acid pinacol ester**

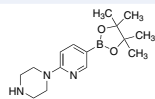
636029
 $C_{12}H_{18}BNO_3$
 MW: 235.09
 [445264-61-9]

**2,4-Dimethoxypyrimidine-5-boronic acid pinacol ester** NEW

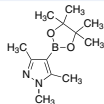
683841
 $C_{12}H_{19}BN_2O_4$
 MW: 266.10

**6-(Piperazin-1-yl)pyridine-3-boronic acid pinacol ester**

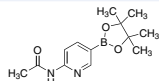
654329
 $C_{12}H_{21}BN_2O_2$
 MW: 236.12
 [871125-86-9]

**1,3,5-Trimethyl-1H-pyrazole-4-boronic acid pinacol ester** NEW

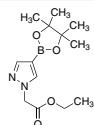
654396
 $C_{12}H_{21}BN_2O_2$
 MW: 236.12

**2-Acetamidopyridine-5-boronic acid pinacol ester** NEW

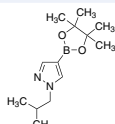
683892
 $C_{13}H_{19}BN_2O_3$
 MW: 262.11
 [904326-87-0]

**1-(Ethoxycarbonylmethyl)-1H-pyrazole-4-boronic acid pinacol ester** NEW

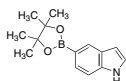
683566
 $C_{13}H_{21}BN_2O_4$
 MW: 280.13
 [864754-16-5]

**1-Isobutylpyrazole-4-boronic acid pinacol ester**

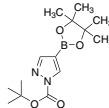
642614
 $C_{13}H_{23}BN_2O_2$
 MW: 250.14
 [827614-66-4]

**5-Indoleboronic acid pinacol ester**

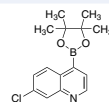
578835
 $C_{14}H_{18}BNO_2$
 MW: 243.11
 [269410-24-4]

**1-Boc-pyrazole-4-boronic acid pinacol ester**

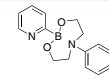
632732
 $C_{14}H_{23}BN_2O_4$
 MW: 294.15
 [552846-17-0]

**7-Chloroquinoline-4-boronic acid pinacol ester**

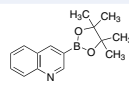
658596
 $C_{15}H_{17}BClNO_2$
 MW: 289.56
 [871125-83-6]

**2-Pyridineboronic acid N-phenyl-diethanolamine ester**

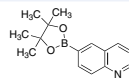
647284
 $C_{15}H_{17}BN_2O_2$
 MW: 268.12
 [662138-96-7]

**3-Quinolineboronic acid pinacol ester**

647950
 $C_{15}H_{18}BNO_2$
 MW: 255.12
 [171364-85-5]

**6-Quinolineboronic acid pinacol ester**

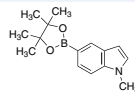
641618
 $C_{15}H_{18}BNO_2$
 MW: 255.12
 [406463-06-7]

**4-Isoquinolineboronic acid pinacol ester**

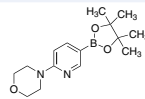
648795
 $C_{15}H_{18}BNO_2$
 MW: 255.12
 [685103-98-4]

**1-Methylindole-5-boronic acid pinacol ester**

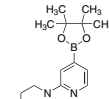
640395
 $C_{15}H_{20}BNO_2$
 MW: 257.14
 [837392-62-8]

**6-(Morpholin-4-yl)pyridine-3-boronic acid pinacol ester**

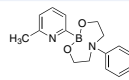
654310
 $C_{15}H_{23}BN_2O_3$
 MW: 290.17
 [485799-04-0]

**2-(1-Piperazinyl)pyridine-4-boronic acid pinacol ester**

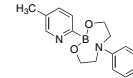
654361
 $C_{15}H_{24}BN_3O_2$
 MW: 289.18
 [871125-87-0]

**6-Methylpyridine-2-boronic acid N-phenyl-diethanolamine ester**

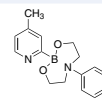
648817
 $C_{16}H_{19}BN_2O_2$
 MW: 282.15
 [872054-59-6]

**5-Methylpyridine-2-boronic acid N-phenyl-diethanolamine ester**

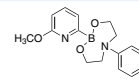
648825
 $C_{16}H_{19}BN_2O_2$
 MW: 282.15
 [872054-54-1]

**4-Methylpyridine-2-boronic acid N-phenyl-diethanolamine ester**

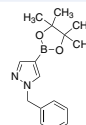
648841
 $C_{16}H_{19}BN_2O_2$
 MW: 282.15
 [849100-03-4]

**6-Methoxypyridine-2-boronic acid N-phenyl-diethanolamine ester**

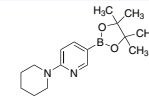
649155
 $C_{16}H_{19}BN_2O_3$
 MW: 298.14
 [872054-59-6]

**1-Benzylpyrazole-4-boronic acid pinacol ester**

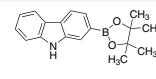
636002
 $C_{16}H_{21}BN_2O_2$
 MW: 284.16
 [761446-45-1]

**6-(Piperidin-1-yl)pyridine-3-boronic acid pinacol ester** NEW

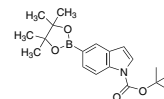
676519
 $C_{16}H_{25}BN_2O_2$
 MW: 288.19
 [852228-08-1]

**9H-Carbazole-2-boronic acid pinacol ester**

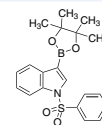
655724
 $C_{18}H_{20}BNO_2$
 MW: 293.17
 [871125-67-6]

**1-Boc-indole-5-boronic acid pinacol ester**

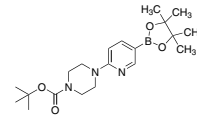
640387
 $C_{19}H_{26}BNO_4$
 MW: 343.23
 [777061-36-6]

**1-(Phenylsulfonyl)indole-3-boronic acid pinacol ester**

654280
 $C_{20}H_{22}BNO_4S$
 MW: 383.27
 [870717-93-4]

**6-(4-Boc-piperazin-1-yl)pyridine-3-boronic acid pinacol ester**

654337
 $C_{20}H_{32}BN_3O_4$
 MW: 389.30
 [496786-98-2]



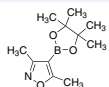
O-Heteroarylboronate esters

3,5-Dimethylisoxazole-4-boronic acid pinacol ester

643882C₁₁H₁₈BN₂O₃

MW: 223.08

[832114-00-8]



3-Furanboronic acid pinacol ester

575054C₁₀H₁₅BO₃

MW: 194.04

[248924-59-6]

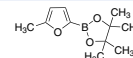


5-Methylfuran-2-boronic acid pinacol ester

NEW

680109C₁₁H₁₇BO₃

MW: 208.06



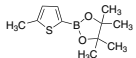
S-Heteroarylboronate esters

5-Methylthiophene-2-boronic acid pinacol ester

655074C₁₁H₁₇BO₂S

MW: 224.13

[476004-80-5]

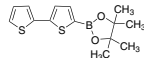


2,2'-Bithiophene-5-boronic acid pinacol ester

578878C₁₄H₁₇BO₂S₂

MW: 292.22

[479719-88-5]

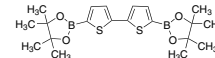


2,2'-Bithiophene-5,5'-diboronic acid bis(pinacol) ester

647020C₂₀H₂₈B₂O₄S₂

MW: 418.19

[239075-02-6]

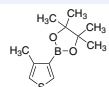


4-Methylthiophene-3-boronic acid pinacol ester

NEW

680001C₁₁H₁₇BO₂S

MW: 224.13

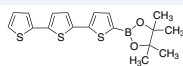


2,2':5',2''-Terthiophene-5-boronic acid pinacol ester

645141C₁₈H₁₉BO₂S₃

MW: 374.35

[849062-17-5]

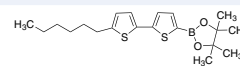


5'-Hexyl-2,2'-bithiophene-5-boronic acid pinacol ester

632961C₂₀H₂₉BO₂S₂

MW: 376.38

[579503-59-6]



Trifluoroborate Salts

Alkyltrifluoroborates

Potassium (bromomethyl)-trifluoroborate

NEW

681342CH₂BBrF₃K

MW: 200.84

[888711-44-2]



Potassium isopropyltrifluoroborate

667153C₃H₇BF₃K

MW: 149.99



Potassium methyltrifluoroborate

637890CH₃BF₃K

MW: 121.94

[13862-28-7]



Potassium butyltrifluoroborate

660094C₄H₉BF₃K

MW: 164.02

[444343-55-9]



Potassium sec-butyltrifluoroborate

667145C₄H₉BF₃K

MW: 164.02



Potassium cyclopropyltrifluoroborate

NEW

662984C₃H₅BF₃K

MW: 147.98



Potassium cyclopentyltrifluoroborate

666017C₅H₉BF₃K

MW: 176.03



Potassium allyltrifluoroborate

659274C₃H₅BF₃K

MW: 147.98



Potassium (3,3-dimethylbutyl)-trifluoroborate

NEW

667161C₆H₁₃BF₃K

MW: 192.07



Potassium benzyltrifluoroborate

563056C₇H₇BF₃K

MW: 198.03

[329976-73-0]



Potassium phenethyltrifluoroborate

563099C₈H₉BF₃K

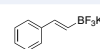
MW: 212.06

[329976-74-1]



Alkenyltrifluoroborates

Potassium vinyltrifluoroborate

655228C₂H₃BF₃K
MW: 133.95
[13682-77-4]Potassium *trans*-styryltrifluoroborate**576158**C₈H₇BF₃K
MW: 210.05
[201852-49-5]Potassium *trans*-3-methoxy-1-propenyltrifluoroborate

NEW

683590C₄H₇BF₃KO
MW: 178.00Potassium *trans*-1-decenyltrifluoroborate**637882**C₁₀H₁₉BF₃K
MW: 246.16
[479678-72-3]

Aryltrifluoroborates

Potassium 2,6-difluorophenyltrifluoroborate

597929C₆H₃BF₅K
MW: 219.99
[267006-25-7]

Potassium 2,4-difluorophenyltrifluoroborate

656992C₆H₃BF₅K
MW: 219.99
[871231-41-3]

Potassium 2-bromophenyltrifluoroborate

576107C₆H₄BBBrF₃K
MW: 262.90
[480445-38-3]

Potassium 4-bromophenyltrifluoroborate

571547C₆H₄BBBrF₃K
MW: 262.90
[374564-35-9]

Potassium 4-chlorophenyltrifluoroborate

657077C₆H₄BClF₃K
MW: 218.45
[661465-44-7]

Potassium 3-nitrophenyltrifluoroborate

571598C₆H₄BF₃KNO₂
MW: 229.01
[192863-40-4]

Potassium 3-fluorophenyltrifluoroborate

659770C₆H₄BF₄K
MW: 202.00
[267006-24-6]

Potassium 4-fluorophenyltrifluoroborate

592846C₆H₄BF₄K
MW: 202.00

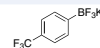
Potassium phenyltrifluoroborate

563951C₆H₅BF₃K
MW: 184.01
[153766-81-5]

Potassium 3-hydroxyphenyltrifluoroborate

659746C₆H₅BF₃KO
MW: 200.01
[871231-45-7]

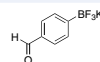
Potassium 4-(trifluoromethyl)phenyltrifluoroborate

576131C₇H₄BF₆K
MW: 252.01
[166328-08-1]

Potassium 3-formylphenyltrifluoroborate

657093C₇H₅BF₃KO
MW: 212.02
[871231-44-6]

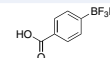
Potassium 4-formylphenyltrifluoroborate

576093C₇H₅BF₃KO
MW: 212.02
[374564-36-0]

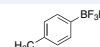
Potassium 3-carboxyphenyltrifluoroborate

659789C₇H₅BF₃KO₂
MW: 228.02
[850313-91-6]

Potassium 4-carboxyphenyltrifluoroborate

657069C₇H₅BF₃KO₂
MW: 228.02
[850623-38-0]

Potassium 3,4-(methylenedioxy)phenyltrifluoroborate

659754C₇H₅BF₃KO₂
MW: 228.02
[871231-46-8]Potassium *o*-tolyltrifluoroborate**576123**C₇H₇BF₃K
MW: 198.03
[274257-34-0]Potassium *p*-tolyltrifluoroborate**571555**C₇H₇BF₃K
MW: 198.03
[216434-82-1]

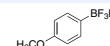
Potassium 2-methoxyphenyltrifluoroborate

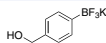
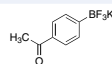
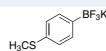
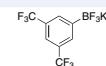
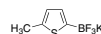
654930C₇H₇BF₃KO
MW: 214.03
[236388-46-8]

Potassium 3-methoxyphenyltrifluoroborate

576115C₇H₇BF₃KO
MW: 214.03
[438553-44-7]

Potassium 4-methoxyphenyltrifluoroborate

579068C₇H₇BF₃KO
MW: 214.03
[192863-36-8]

Potassium 4-(hydroxymethyl)phenyltrifluoroborate**659762**C₇H₇BF₃KO
MW: 214.03**Potassium 4-acetylphenyltrifluoroborate****657050**C₈H₇BF₃KO
MW: 226.05
[252726-24-2]**Potassium 2-naphthalenetrifluoroborate****657018**C₁₀H₇BF₃K
MW: 234.07
[668984-08-5]**Potassium 4-(methylthio)phenyltrifluoroborate****657115**C₇H₇BF₃KS
MW: 230.10
[871231-43-5]**Potassium 2,6-dimethylphenyltrifluoroborate****592943**C₈H₉BF₃K
MW: 212.06**Potassium 4-tert-butylphenyltrifluoroborate****654728**C₁₀H₁₃BF₃K
MW: 240.11
[423118-47-2]**Potassium 3,5-bis(trifluoromethyl)phenyltrifluoroborate****571563**C₈H₃BF₉K
MW: 320.00
[166328-09-2]**Potassium 2,6-dimethoxyphenyltrifluoroborate****657085**C₈H₉BF₃KO₂
MW: 244.06
[871231-42-4]**S-Heteroaryltrifluoroborates****Potassium 3-thiophenetrifluoroborate****571571**C₄H₃BF₃KS
MW: 190.04
[192863-37-9]**Potassium 5-methyl-2-thiophene trifluoroborate****654949**C₅H₅BF₃KS
MW: 204.06
[871231-40-2]

Argentina

SIGMA-ALDRICH DE ARGENTINA S.A.
Free Tel: 0810 888 7446
Tel: (+54) 11 4556 1472
Fax: (+54) 11 4552 1698

Australia

SIGMA-ALDRICH PTY LTD.
Free Tel: 1800 800 097
Free Fax: 1800 800 096
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Austria

SIGMA-ALDRICH HANDELS GmbH
Tel: (+43) 1 605 81 10
Fax: (+43) 1 605 81 20

Belgium

SIGMA-ALDRICH NV/SA.
Free Tel: 0800 14747
Free Fax: 0800 14745
Tel: (+32) 3 899 13 01
Fax: (+32) 3 899 13 11

Brazil

SIGMA-ALDRICH BRASIL LTDA.
Free Tel: 0800 701 7425
Tel: (+55) 11 3732 3100
Fax: (+55) 11 5522 9895

Canada

SIGMA-ALDRICH CANADA LTD.
Free Tel: 1800 565 1400
Free Fax: 1800 265 3858
Tel: (+1) 905 829 9500
Fax: (+1) 905 829 9292

China

SIGMA-ALDRICH (SHANGHAI)
TRADING CO. LTD.
Free Tel: 800 819 3336
Tel: (+86) 21 6141 5566
Fax: (+86) 21 6141 5567

Czech Republic

SIGMA-ALDRICH S.R.O.
Tel: (+420) 246 003 200
Fax: (+420) 246 003 291

Denmark

SIGMA-ALDRICH DENMARK A/S
Tel: (+45) 43 56 59 10
Fax: (+45) 43 56 59 05

Finland

SIGMA-ALDRICH FINLAND OY
Tel: (+358) 9 350 9250
Fax: (+358) 9 350 9255

France

SIGMA-ALDRICH CHIMIE S.à.r.l.
Free Tel: 0800 211 408
Free Fax: 0800 031 052
Tel: (+33) 474 82 28 00
Fax: (+33) 474 95 68 08

Germany

SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800 51 55 000
Free Fax: 0800 64 90 000
Tel: (+49) 89 6513 0
Fax: (+49) 89 6513 1160

Greece

SIGMA-ALDRICH (O.M.) LTD.
Tel: (+30) 210 994 8010
Fax: (+30) 210 994 3831

Hungary

SIGMA-ALDRICH Kft
Ingyenes zöld telefon: 06 80 355 355
Ingyenes zöld fax: 06 80 344 344
Tel: (+36) 1 235 9055
Fax: (+36) 1 235 9050

India

SIGMA-ALDRICH CHEMICALS
PRIVATE LIMITED
Telephone
Bangalore: (+91) 80 6621 9600
New Delhi: (+91) 11 4165 4255
Mumbai: (+91) 22 2570 2364
Hyderabad: (+91) 40 6684 5488
Fax
Bangalore: (+91) 80 6621 9650
New Delhi: (+91) 11 4165 4266
Mumbai: (+91) 22 2579 7589
Hyderabad: (+91) 40 6684 5466

Ireland

SIGMA-ALDRICH IRELAND LTD.
Free Tel: 1800 200 888
Free Fax: 1800 600 222
Tel: (+353) 1 404 1900
Fax: (+353) 1 404 1910

Israel

SIGMA-ALDRICH ISRAEL LTD.
Free Tel: 1 800 70 2222
Tel: (+972) 8 948 4100
Fax: (+972) 8 948 4200

Italy

SIGMA-ALDRICH S.r.l.
Numero Verde: 800 827018
Tel: (+39) 02 3341 7310
Fax: (+39) 02 3801 0737

Japan

SIGMA-ALDRICH JAPAN K.K.
Tokyo Tel: (+81) 3 5796 7300
Tokyo Fax: (+81) 3 5796 7315

Korea

SIGMA-ALDRICH KOREA
Free Tel: (+82) 80 023 7111
Free Fax: (+82) 80 023 8111
Tel: (+82) 31 329 9000
Fax: (+82) 31 329 9090

Malaysia

SIGMA-ALDRICH (M) SDN. BHD
Tel: (+60) 3 5635 3321
Fax: (+60) 3 5635 4116

Mexico

SIGMA-ALDRICH QUÍMICA, S.A. de C.V.
Free Tel: 01 800 007 5300
Free Fax: 01 800 712 9920
Tel: 52 722 276 1600
Fax: 52 722 276 1601

The Netherlands

SIGMA-ALDRICH CHEMIE BV
Free Tel: 0800 022 9088
Free Fax: 0800 022 9089
Tel: (+31) 78 620 5411
Fax: (+31) 78 620 5421

New Zealand

SIGMA-ALDRICH NEW ZEALAND LTD.
Free Tel: 0800 936 666
Free Fax: 0800 937 777
Tel: (+61) 2 9841 0555
Fax: (+61) 2 9841 0500

Norway

SIGMA-ALDRICH NORWAY AS
Tel: (+47) 23 17 60 60
Fax: (+47) 23 17 60 50

Poland

SIGMA-ALDRICH Sp. z o.o.
Tel: (+48) 61 829 01 00
Fax: (+48) 61 829 01 20

Portugal

SIGMA-ALDRICH QUÍMICA, S.A.
Free Tel: 800 202 180
Free Fax: 800 202 178
Tel: (+351) 21 924 2555
Fax: (+351) 21 924 2610

Russia

SIGMA-ALDRICH RUS, LLC
Tel: +7 (495) 621 6037
Fax: +7 (495) 621 5923

Singapore

SIGMA-ALDRICH PTE. LTD.
Tel: (+65) 6779 1200
Fax: (+65) 6779 1822

South Africa

SIGMA-ALDRICH
SOUTH AFRICA (PTY) LTD.
Free Tel: 0800 1100 75
Free Fax: 0800 1100 79
Tel: (+27) 11 979 1188
Fax: (+27) 11 979 1119

Spain

SIGMA-ALDRICH QUÍMICA, S.A.
Free Tel: 900 101 376
Free Fax: 900 102 028
Tel: (+34) 91 661 99 77
Fax: (+34) 91 661 96 42

Sweden

SIGMA-ALDRICH SWEDEN AB
Tel: (+46) 8 742 4200
Fax: (+46) 8 742 4243

Switzerland

SIGMA-ALDRICH CHEMIE GmbH
Free Tel: 0800 80 00 80
Free Fax: 0800 80 00 81
Tel: (+41) 81 755 2828
Fax: (+41) 81 755 2815

United Kingdom

SIGMA-ALDRICH COMPANY LTD.
Free Tel: 0800 717 181
Free Fax: 0800 378 785
Tel: (+44) 1747 833 000
Fax: (+44) 1747 833 313
SAFC (UK) Free Tel: 0800 71 71 17

United States

SIGMA-ALDRICH
P.O. Box 14508
St. Louis, Missouri 63178
Toll-Free: 800 325 3010
Toll-Free Fax: 800 325 5052
Call Collect: (+1) 314 771 5750
Tel: (+1) 314 771 5765
Fax: (+1) 314 771 5757

Internet

sigma-aldrich.com

World Headquarters

3050 Spruce St., St. Louis, MO 63103
(314) 771-5765
sigma-aldrich.com

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052

Technical Service (800) 325-5832 • sigma-aldrich.com/techservice

Development/Bulk Manufacturing Inquiries SAFC™ (800) 244-1173

The SIGMA-ALDRICH Group

©2007 Sigma-Aldrich Co. All rights reserved.

SIGMA, SAFC, SAFC, SIGMA-ALDRICH, ISOTEC, ALDRICH, FLUKA, and SUPELCO are trademarks belonging to Sigma-Aldrich Co. and its affiliate Sigma-Aldrich Biotechnology, L.P. Riedel-de Haën® trademark under license from Riedel-de Haën GmbH. Sigma brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.



SIGMA-ALDRICH

*Accelerating Customers' Success
through Leadership in Life Science,
High Technology and Service*

JRD
02750-504740
06-2007