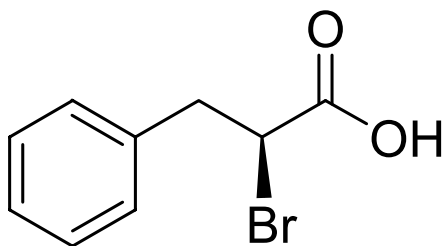


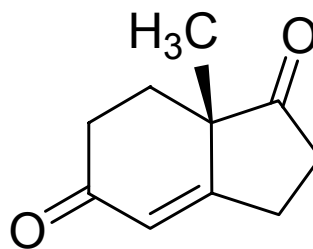
FLUKA 00031 (35016-63-8)

(S)-2-Bromo-3-phenylpropionic acid, technical, >= 95.0 % T



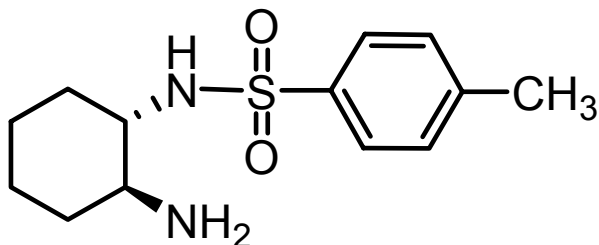
ALDRICH 664073 (17553-86-5)

(S)-(+)-2,3,7,7a-Tetrahydro-7a-methyl-1H-indene-1,5(6H)-dione, 97 %



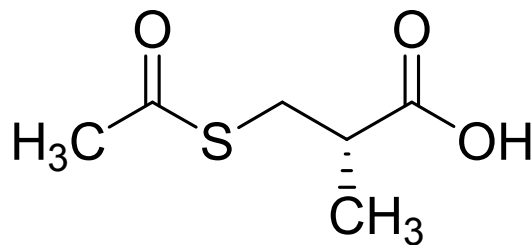
ALDRICH 677310

(1S,2S)-(+)-N-p-Tosyl-1,2-cyclohexanediamine, 97 %



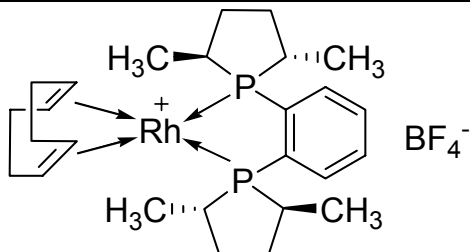
ALDRICH 472522 (76497-39-7)

(S)-(-)-3-Acetylthio-2-methylpropanoic acid, 97 %



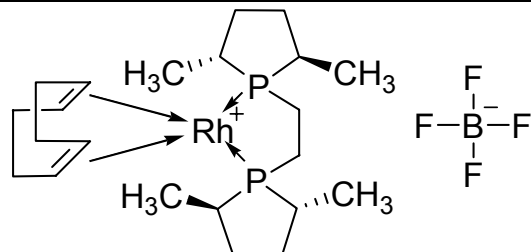
ALDRICH 675539 (205064-10-4)

1,2-Bis[(2S,5S)-2,5-dimethylphospholano]benzene(cyclooctadiene)rhodium(I) tetrafluoroborate



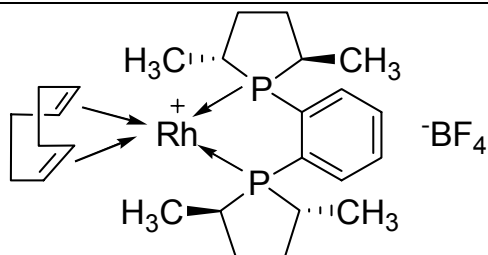
ALDRICH 675555 (305818-67-1)

1,2-Bis[(2R,5R)-2,5-dimethylphospholano]ethane(cyclooctadiene)rhodium(I) tetrafluoroborate



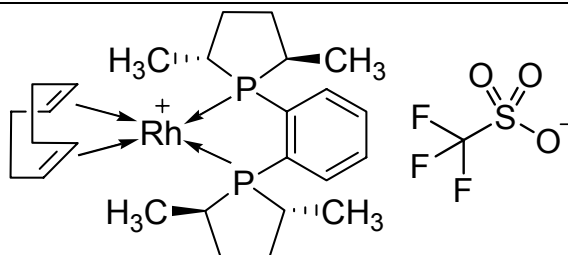
ALDRICH 675520 (210057-23-1)

1,2-Bis[(2R,5R)-2,5-dimethylphospholano]benzene(cyclooctadiene)rhodium(I) tetrafluoroborate



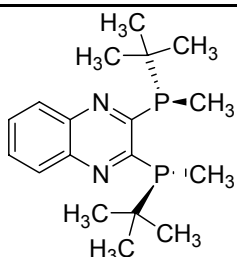
ALDRICH 675571 (187682-63-9)

1,2-Bis[(2R,5R)-2,5-dimethylphospholano]benzene(cyclooctadiene)rhodium(I) trifluoromethanesulfonate, >= 97 %



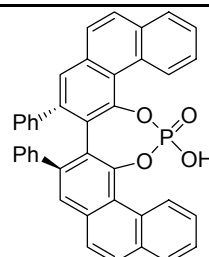
ALDRICH 676403

(R,R)-(-)-2,3-Bis(tert-butylmethylphosphino)quinoxaline, 97 %



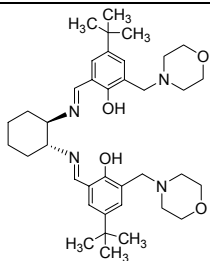
ALDRICH 675512 (871130-18-6)

(R)-(-)-VAPOL hydrogenphosphate



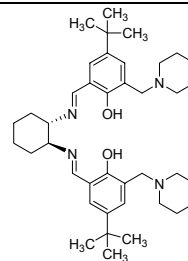
ALDRICH 673331 (323193-85-7)

2,2'-[[[(1R,2R)-(-)-1,2-Cyclohexanediylbis[(E)-(nitrilomethylidene)]]bis[4-(tert-butyl)-6-(4-morpholinylmethyl)phenol], 97 %



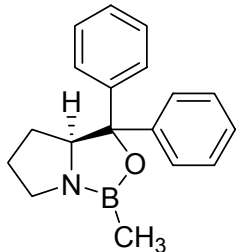
ALDRICH 673366 (478282-27-8)

2,2'-[[[(1S,2S)-1,2-Cyclohexanediylbis[(E)-(nitrilomethylidene)]]bis[4-(tert-butyl)-6-(4-piperidinylmethyl)phenol], 97 %



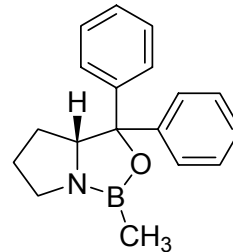
ALDRICH 674648 (112022-81-8)

(S)-(-)-2-Methyl-CBS-oxazaborolidine solution, 1M THF



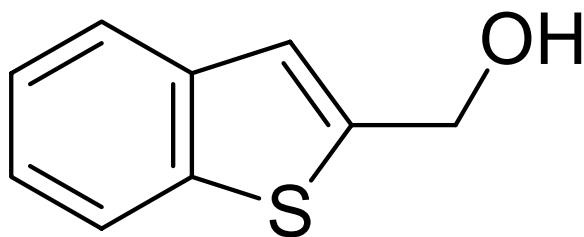
ALDRICH 674656 (112022-83-0)

(R)-(+)-2-Methyl-CBS-oxazaborolidine solution, 1M THF



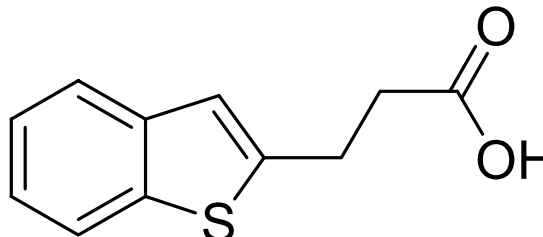
ALDRICH 679542 (17890-56-1)

Benzothiophene-2-methanol, 97 %



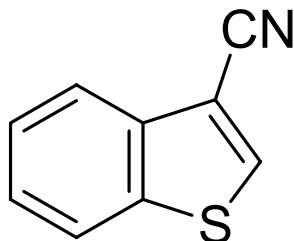
ALDRICH 679569 (42768-60-5)

Benzo[b]thiophene-2-propionic acid, 97 %



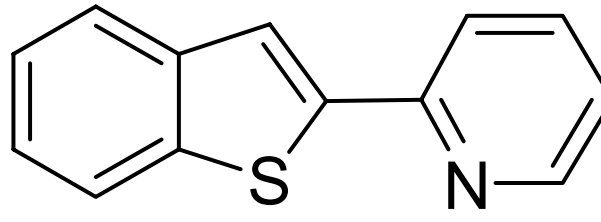
ALDRICH 665975 (24434-84-2)

Benzo[b]thiophene-3-carbonitrile, 97 %



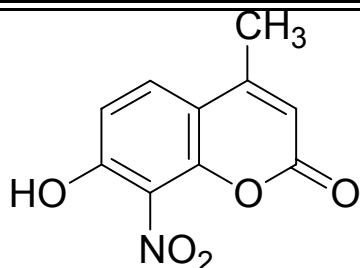
ALDRICH 667684 (38210-35-4)

2-(2-Pyridyl)benzothiophene, 97 %



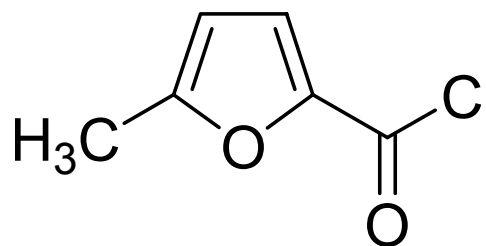
ALDRICH 640646 (19037-69-5)

7-Hydroxy-4-methyl-8-nitrocoumarin, 96 %



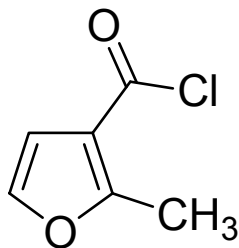
ALDRICH 678570 (14003-11-3)

5-Methylfuran-2-carbonyl chloride, 97 %



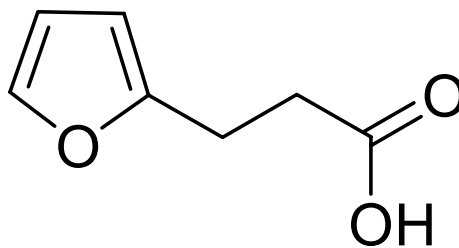
ALDRICH 678597 (5555-00-0)

2-Methylfuran-3-carbonyl chloride, 97 %



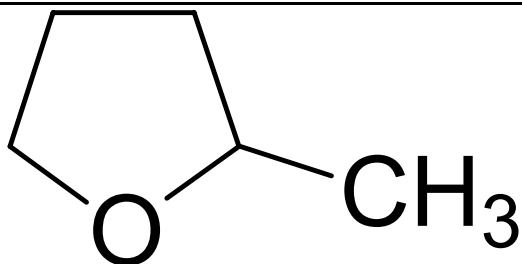
ALDRICH 667730 (935-13-7)

3-(2-Furyl)propionic acid, 97 %



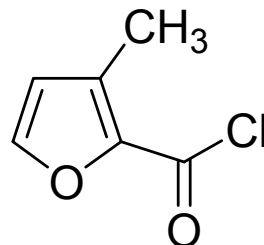
ALDRICH 673277 (96-47-9)

2-Methyltetrahydrofuran, >= 99 %, anhydrous, Inhibitor-free



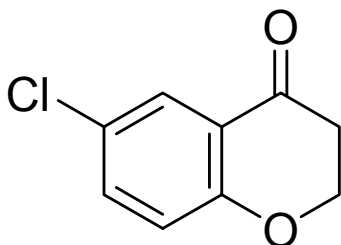
ALDRICH 678589 (22601-06-5)

3-Methylfuran-2-carbonyl chloride, 97 %



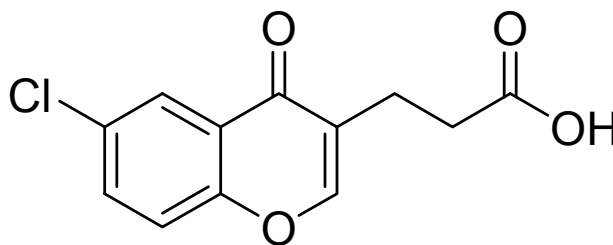
ALDRICH 640476 (37674-72-9)

6-Chloro-4-chromanone, 97 %



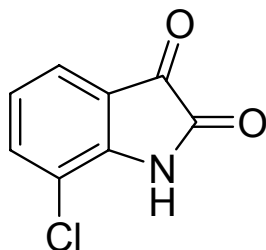
ALDRICH 658014

6-Chlorochromone-3-propionic acid, 96 %



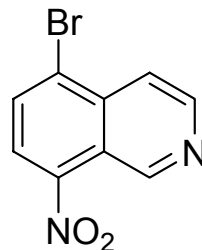
ALDRICH 675040 (7477-63-6)

7-Chloroisatin, 97 %



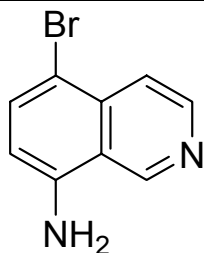
ALDRICH 675148 (63927-23-1)

5-Bromo-8-nitroisquinoline, 90 %



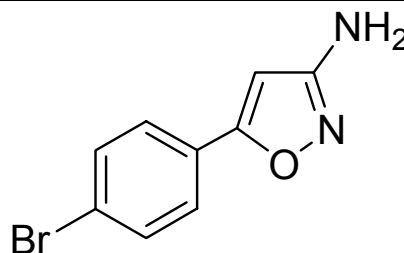
ALDRICH 675954 (90721-35-0)

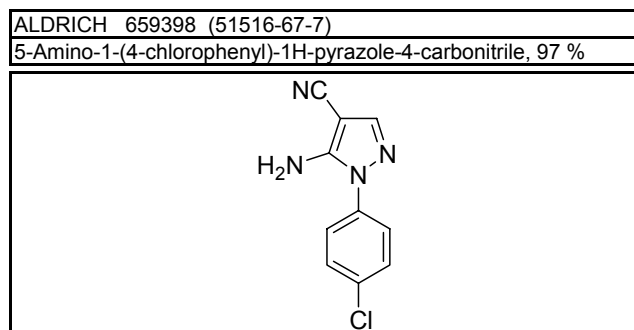
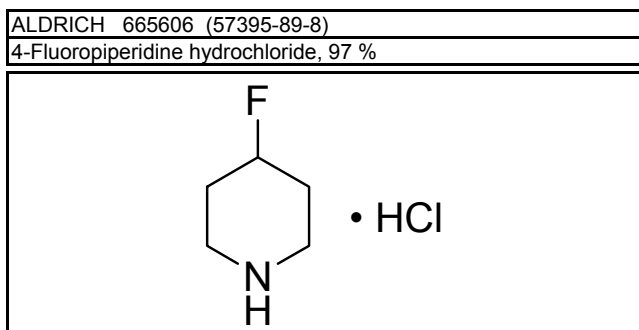
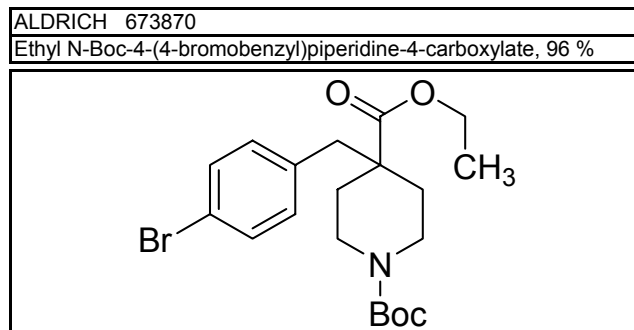
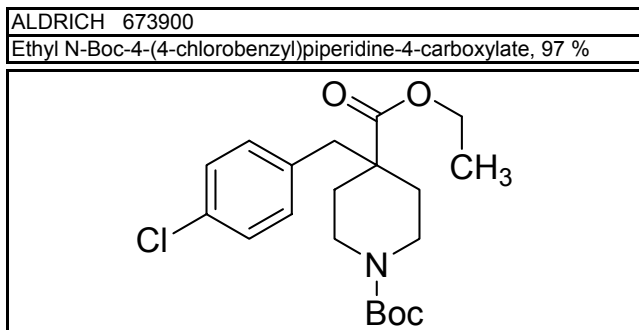
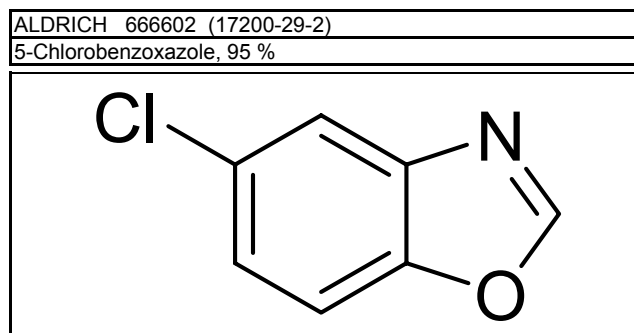
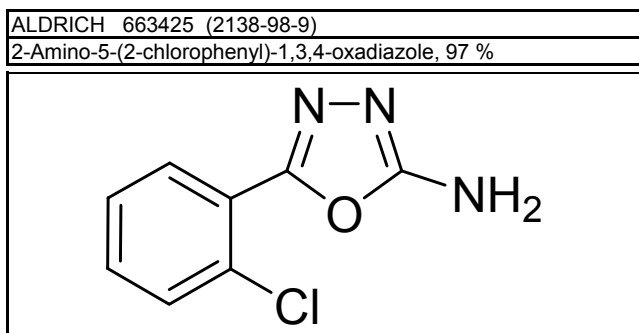
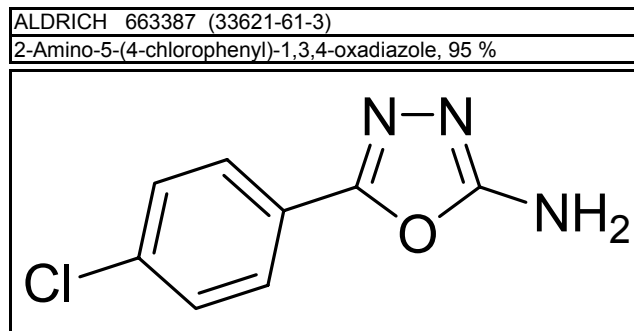
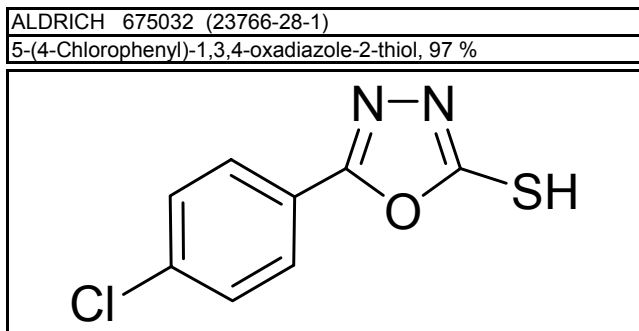
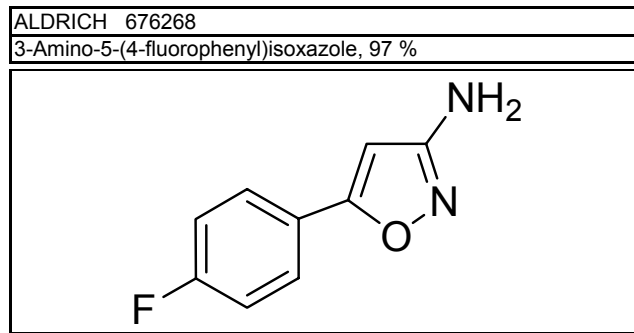
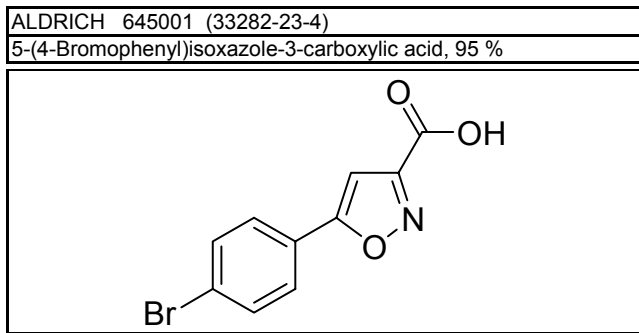
8-Amino-5-bromoisquinoline, 90 %



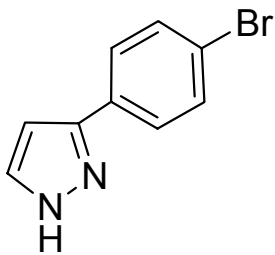
ALDRICH 676276

3-Amino-5-(4-bromophenyl)isoxazole, 97 %

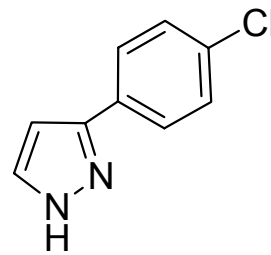




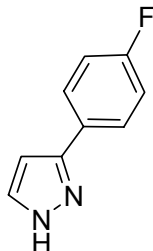
ALDRICH 676292 (73387-46-9)
3-(4-Bromophenyl)-1H-pyrazole, 97 %



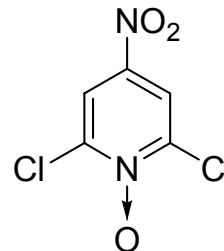
ALDRICH 676284 (59843-58-2)
3-(4-Chlorophenyl)-1H-pyrazole, 97 %



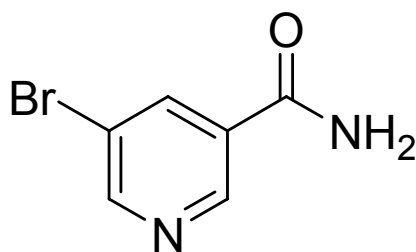
ALDRICH 677221 (154258-82-9)
3-(4-Fluorophenyl)-1H-pyrazole, 97 %



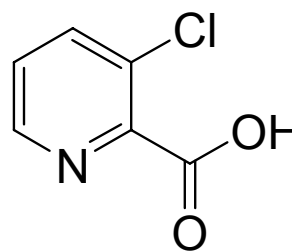
ALDRICH 673161 (2587-01-1)
2,6-Dichloro-4-nitropyridine-N-oxide, 97 %



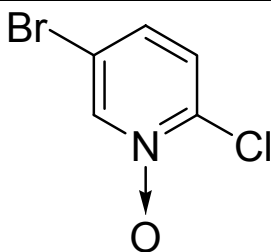
ALDRICH 667137 (28733-43-9)
5-Bromo-3-pyridinecarboxamide, 97 %



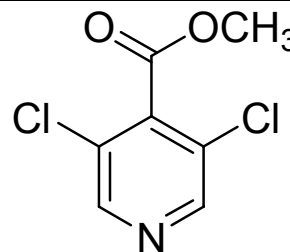
ALDRICH 680052 (57266-69-0)
3-Chloropyridine-2-carboxylic acid, 97 %



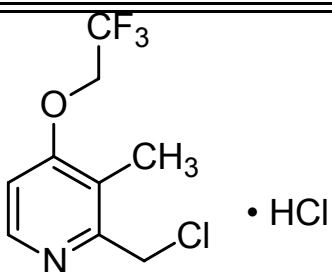
ALDRICH 677833
5-Bromo-2-chloropyridine N-oxide, 97 %



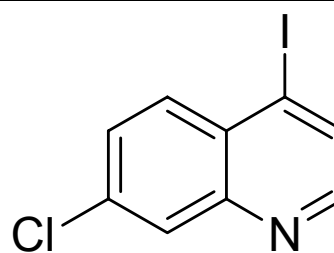
ALDRICH 679577 (773874-72-9)
Methyl 3,5-dichloropyridine-4-carboxylate, 97 %



ALDRICH 631388 (127337-60-4)
2-(Chloromethyl)-3-methyl-4-(2,2,2-trifluoroethoxy)pyridine hydrochloride, 97 %

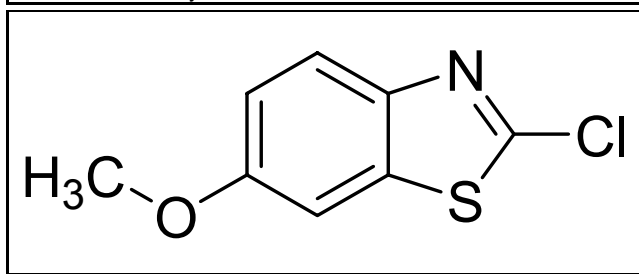


ALDRICH 652903 (98591-57-2)
7-Chloro-4-iodoquinoline, purum, 97 %



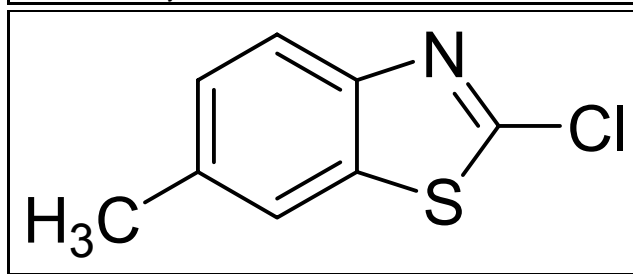
ALDRICH 667331 (2605-14-3)

2-Chloro-6-methoxybenzothiazole, 97 %



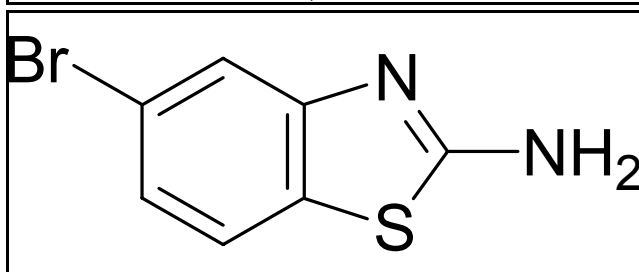
ALDRICH 667374 (3507-26-4)

2-Chloro-6-methylbenzothiazole, 96 %



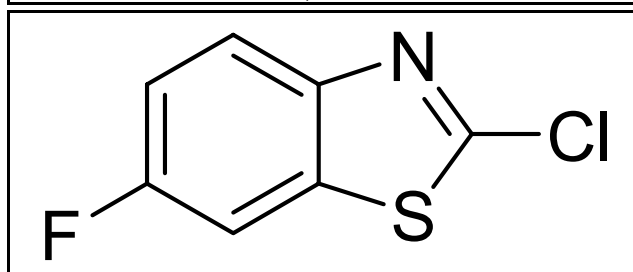
ALDRICH 647683 (20358-03-6)

2-Amino-5-bromobenzothiazole, 97 %



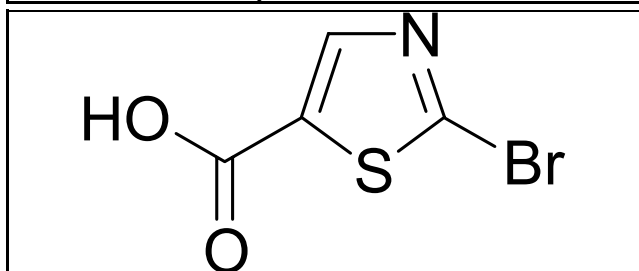
ALDRICH 667358 (399-74-6)

2-Chloro-6-fluorobenzothiazole, 97 %



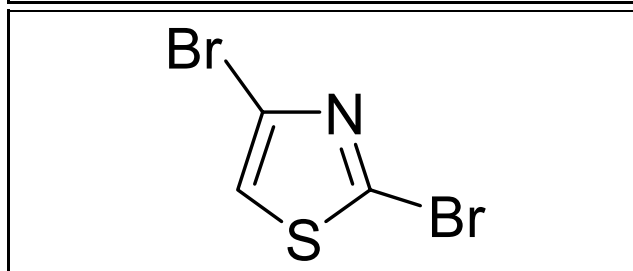
ALDRICH 642487 (54045-76-0)

2-Bromothiazole-5-carboxylic acid, 97 %



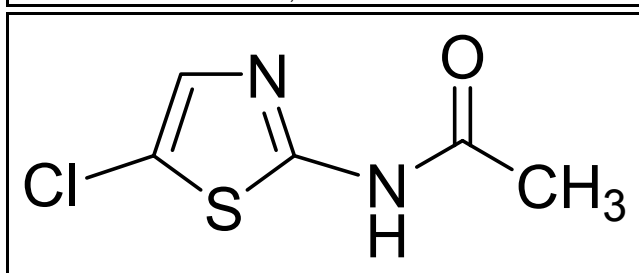
ALDRICH 677914 (4175-77-3)

2,4-Dibromothiazole, 97 %



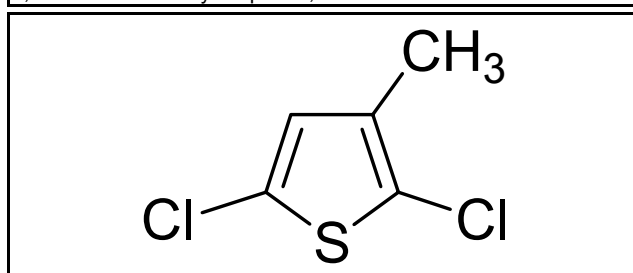
ALDRICH 673129 (20256-39-7)

2-Acetamido-5-chlorothiazole, 96 %



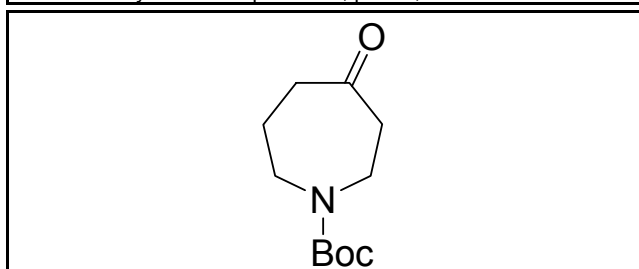
ALDRICH 668362 (17249-90-0)

2,5-Dichloro-3-methylthiophene, 96 %



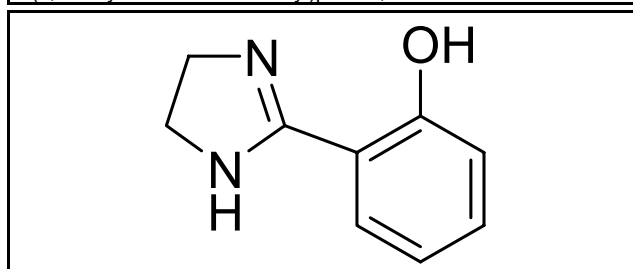
FLUKA 00833 (188975-88-4)

N-Boc-hexahydro-1H-azepin-4-one, purum, >= 95.0 % HPLC



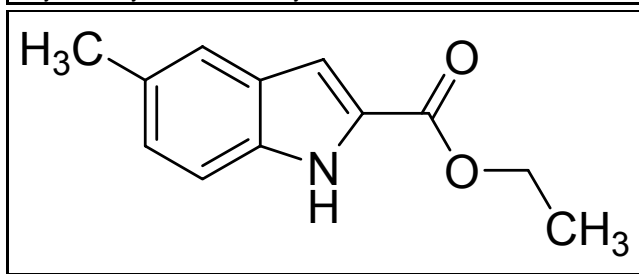
ALDRICH 654698 (1565-39-5)

2-(4,5-Dihydro-1H-imidazol-2-yl)phenol, 97 %



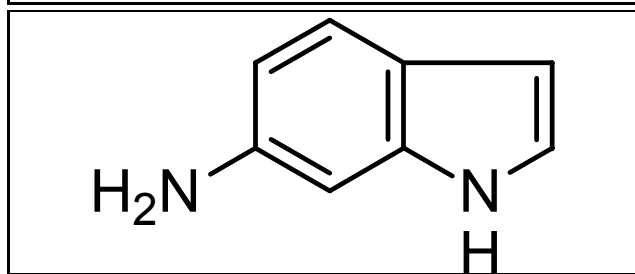
ALDRICH 663190 (16382-15-3)

Ethyl 5-methylindole-2-carboxylate, 97 %



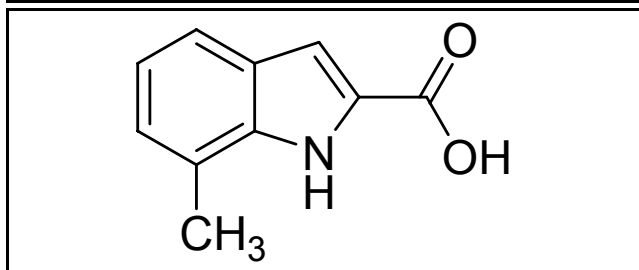
ALDRICH 630721 (5318-27-4)

6-Aminoindole, 97 %



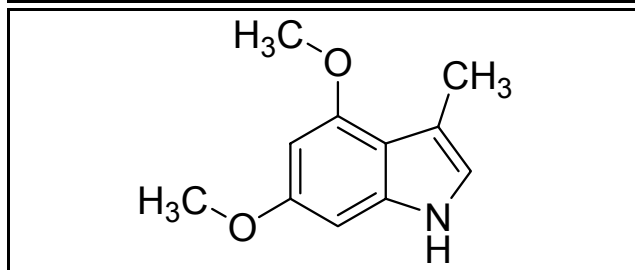
ALDRICH 668680 (18474-60-7)

7-Methylindole-2-carboxylic acid, 95 %



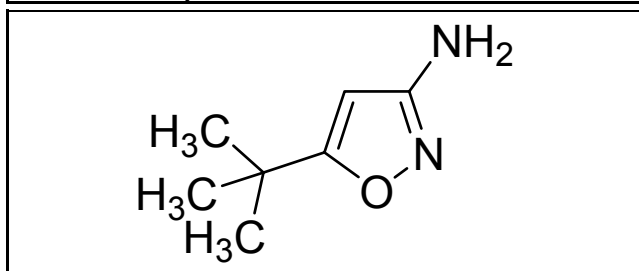
ALDRICH 658588 (74973-30-1)

4,6-Dimethoxy-3-methylindole, 97 %



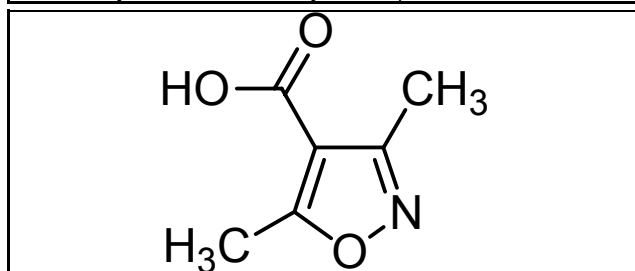
ALDRICH 666580 (55809-36-4)

3-Amino-5-tert-butylisoxazole, 97 %



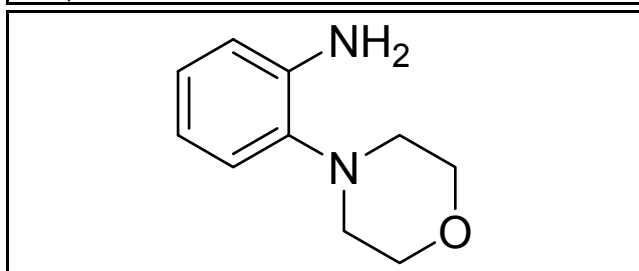
ALDRICH 652873 (2510-36-3)

3,5-Dimethylisoxazole-4-carboxylic acid, puriss., 98 %



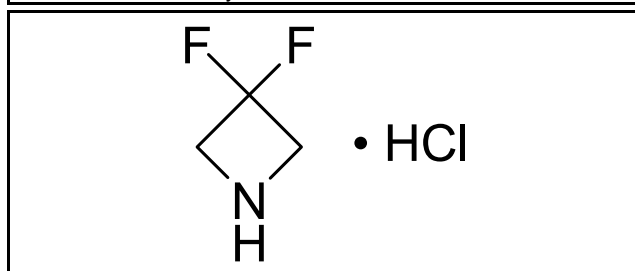
ALDRICH 663204 (5585-33-1)

2-Morpholinoaniline, 97 %



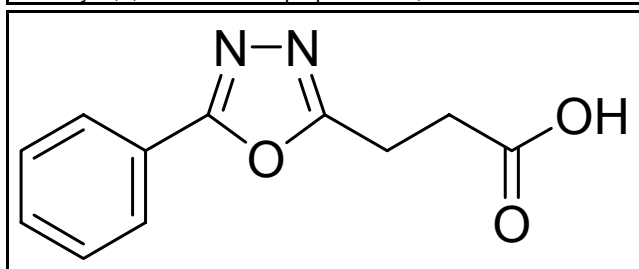
ALDRICH 665584 (288315-03-7)

3,3-Difluoroazetidine hydrochloride, 97 %



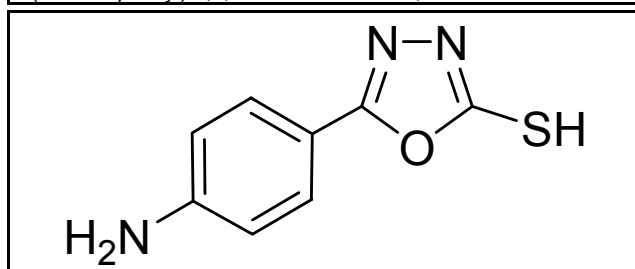
ALDRICH 663409 (23464-98-4)

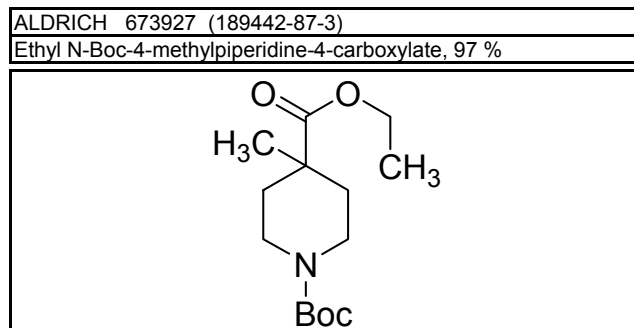
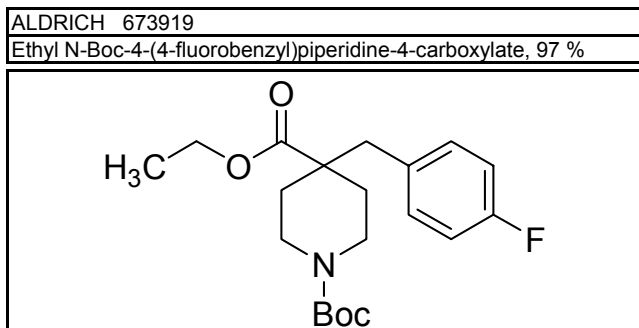
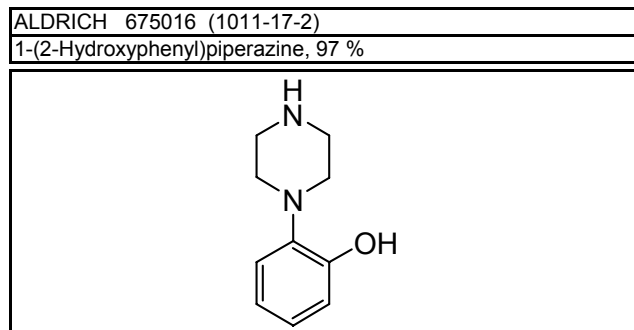
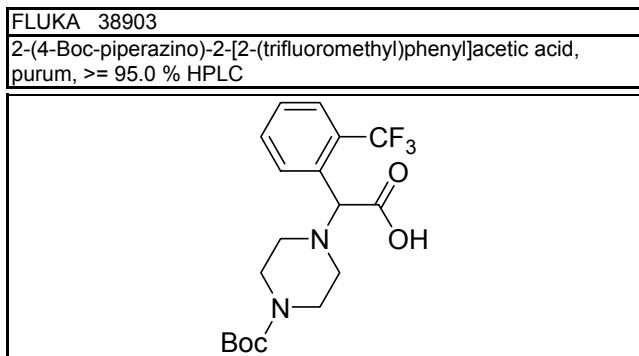
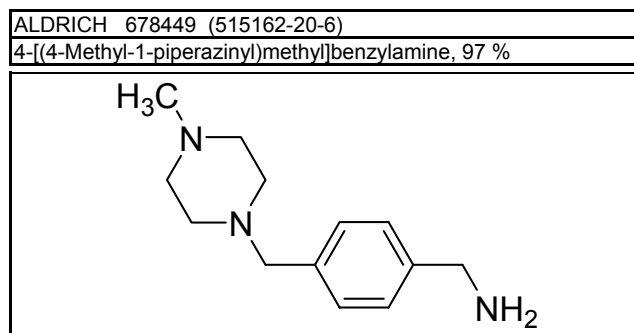
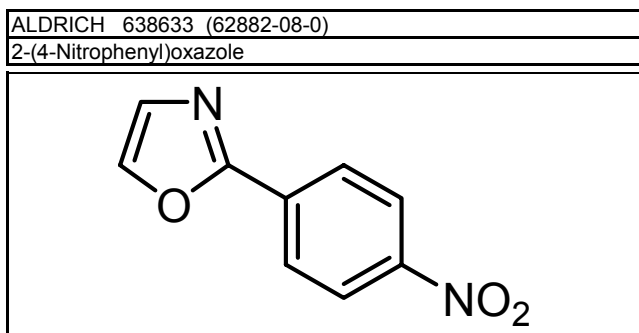
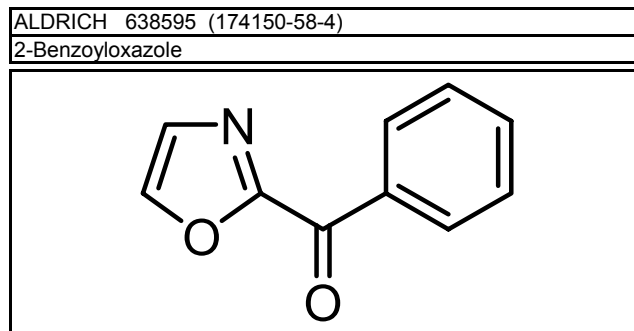
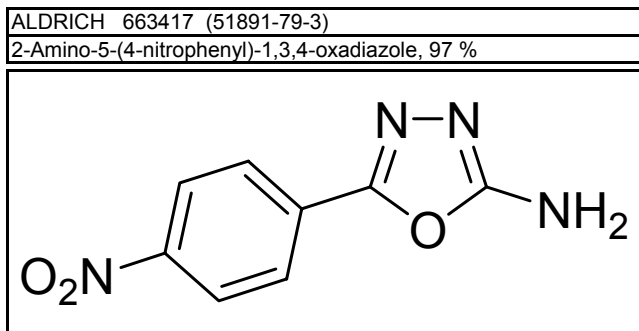
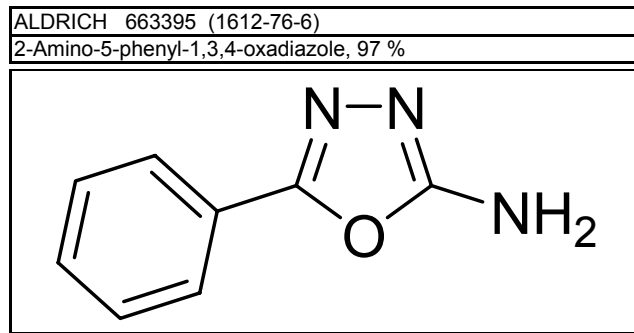
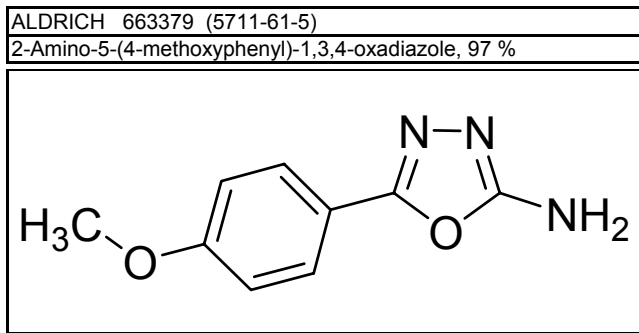
5-Phenyl-1,3,4-oxadiazole-2-propionic acid, 97 %



ALDRICH 675024 (32058-82-5)

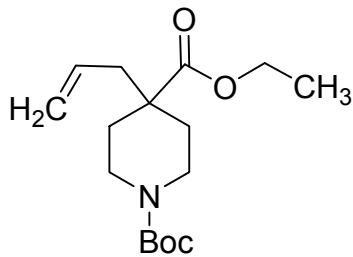
5-(4-Aminophenyl)-1,3,4-oxadiazole-2-thiol, 97 %





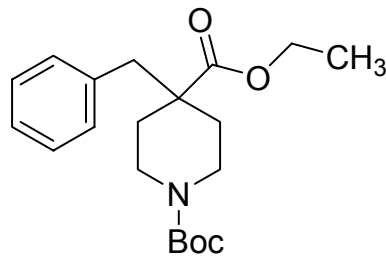
ALDRICH 673730

Ethyl N-Boc-4-allylpiperidine-4-carboxylate, 90 %



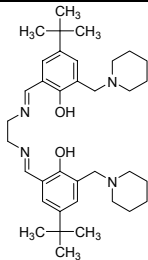
ALDRICH 673749

Ethyl N-Boc-4-benzylpiperidine-4-carboxylate, 96 %



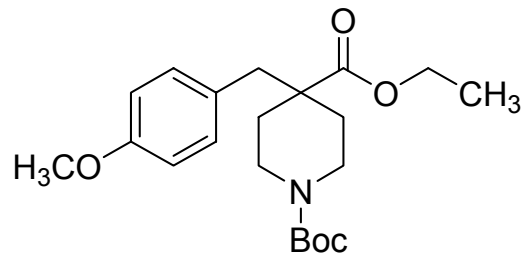
ALDRICH 674486

2,2'-[1,2-Ethanediy]bis[(E)-(nitrilomethylidene)] bis[4-(tert-butyl)-6-(1-piperidinylmethyl)phenol], 97 %



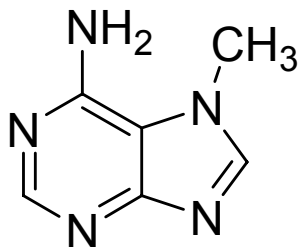
ALDRICH 673897

Ethyl N-Boc-4-(4-methoxybenzyl)piperidine-4-carboxylate, 97 %



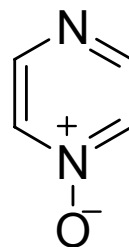
ALDRICH 666548 (935-69-3)

7-Methyladenine, 97 %



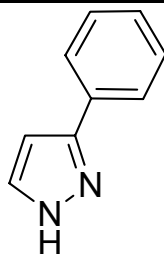
ALDRICH 678260 (2423-65-6)

Pyrazine N-oxide, 97 %



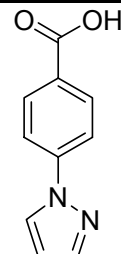
ALDRICH 676349

3-Phenyl-1H-pyrazole, 97 %



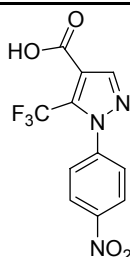
ALDRICH 668672 (16209-00-0)

4-(1H-Pyrazol-1-yl)benzoic acid, 90 %



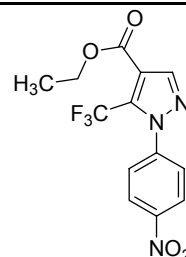
ALDRICH 659347 (142818-03-9)

1-(4-Nitrophenyl)-5-(trifluoromethyl)-1H-pyrazole-4-carboxylic acid, 97 %



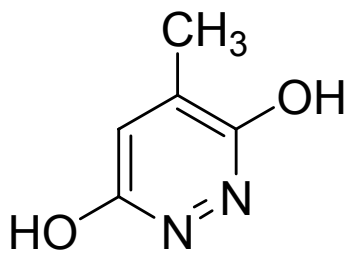
ALDRICH 659355 (175137-35-6)

Ethyl 1-(4-nitrophenyl)-5-(trifluoromethyl)-1H-pyrazole-4-carboxylate, 97 %



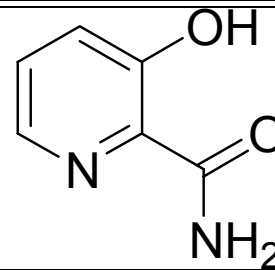
ALDRICH 668141 (5754-18-7)

3,6-Dihydroxy-4-methylpyridazine, 97 %



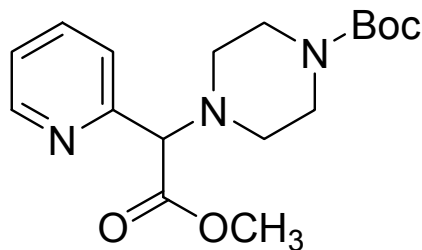
ALDRICH 646423 (933-90-4)

3-Hydroxypicolinamide, 98 %



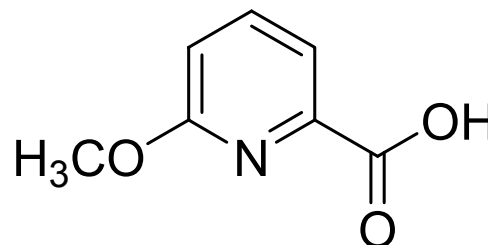
FLUKA 19578

Methyl 2-(4-Boc-piperazino)-2-(2-pyridyl)acetate, purum, >= 95.0 % HPLC



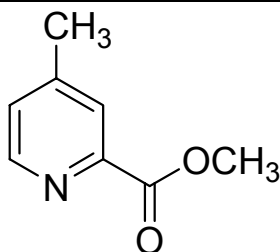
ALDRICH 679607

6-Methoxypyridine-2-carboxylic acid, 97 %



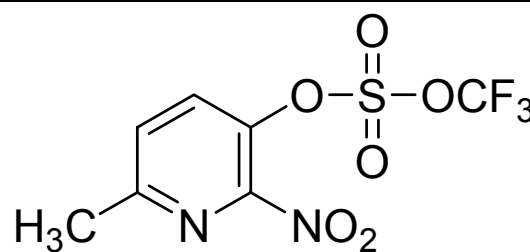
ALDRICH 678368

Methyl 4-methyl-2-pyridinecarboxylate, 97 %



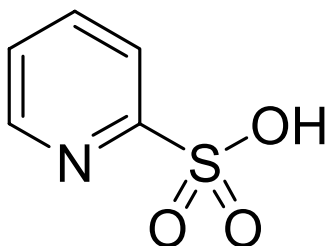
ALDRICH 649937 (163083-48-5)

6-Methyl-2-nitro-3-pyridyl trifluoromethanesulfonate, 97 %



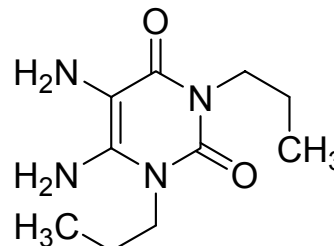
ALDRICH 661759 (15103-48-7)

2-Pyridinesulfonic acid, 97 %



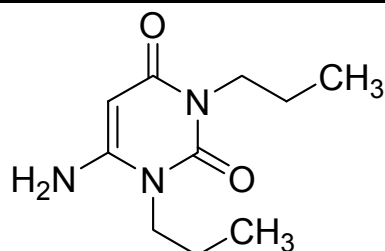
ALDRICH 665452

5,6-Diamino-1,3-dipropyluracil, 97 %



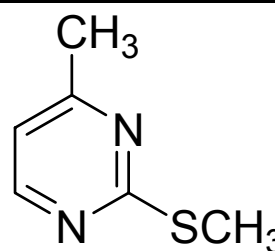
ALDRICH 665444 (41862-14-0)

6-Amino-1,3-dipropyluracil, 97 %

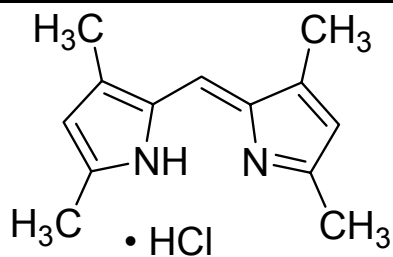


ALDRICH 674885 (14001-63-9)

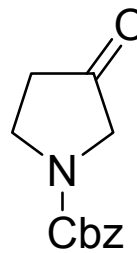
4-Methyl-2-(methylthio)pyrimidine, 97 %



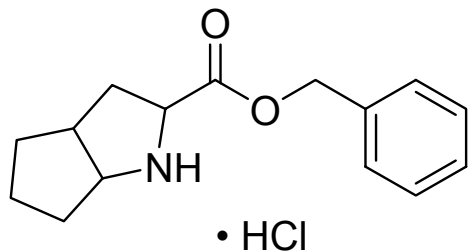
ALDRICH 678058 (67036-97-9)
2-[(3,5-Dimethyl-2H-pyrrol-2-ylidene)methyl]-3,5-dimethyl-1H-pyrrole monohydrochloride, 97 %



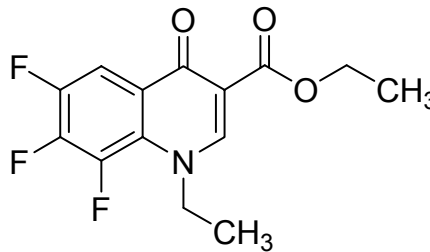
ALDRICH 661201 (130312-02-6)
1-Z-3-Pyrrolidinone, 96 %



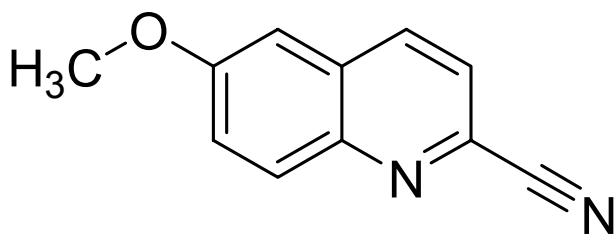
ALDRICH 652547 (93779-29-4)
Benzyl 2-azabicyclo[3.3.0]octane-3-carboxylate hydrochloride, 97 %



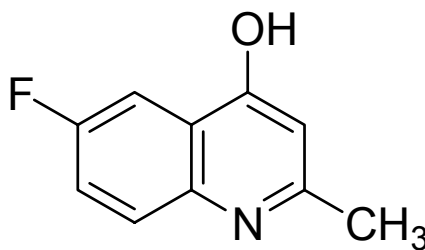
ALDRICH 677701 (100501-62-0)
Ethyl 1-ethyl-6,7,8-trifluoro-4-oxo-1,4-dihydroquinoline-3-carboxylate, 97 %



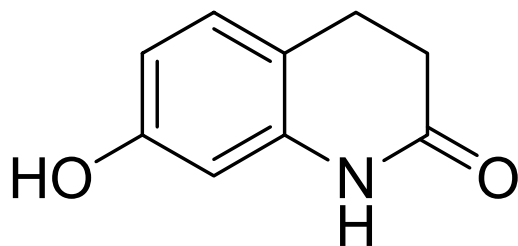
ALDRICH 673528 (5467-79-8)
6-Methoxy-2-quinolinecarbonitrile, 97 %



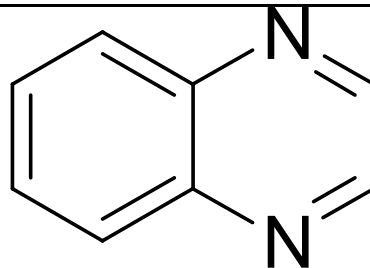
ALDRICH 586501
6-Fluoro-4-hydroxy-2-methylquinoline, 97 %



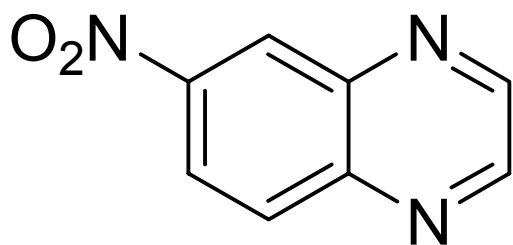
ALDRICH 675369 (22246-18-0)
7-Hydroxy-3,4-dihydro-2(1H)-quinolinone, 97 %



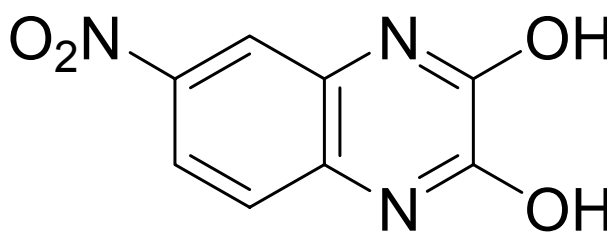
ALDRICH Q1603 (91-19-0)
Quinoxaline, 99 %



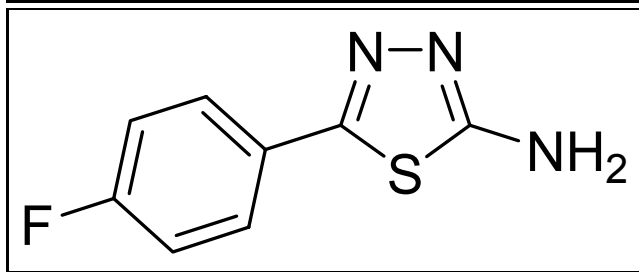
ALDRICH 678384 (6639-87-8)
6-Nitroquinoxaline, 97 %



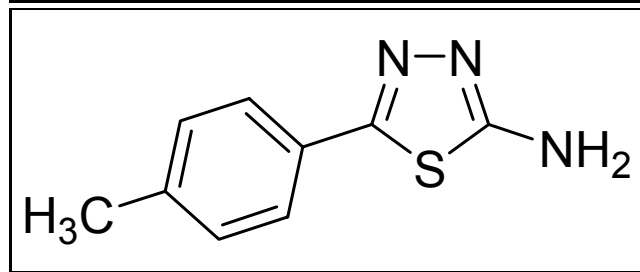
ALDRICH 657239 (2379-56-8)
6-Nitro-2,3-dihydroxyquinoxaline, 96 %



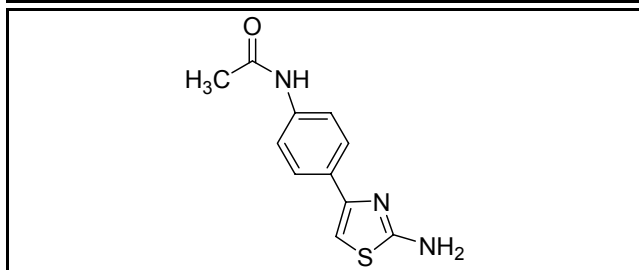
ALDRICH 663557 (942-70-1)
2-Amino-5-(4-fluorophenyl)-1,3,4-thiadiazole, 97 %



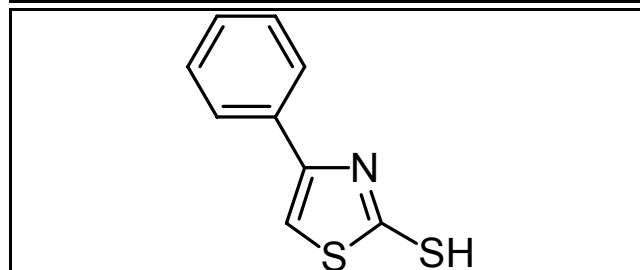
ALDRICH 663565 (26907-54-0)
2-Amino-5-(4-methylphenyl)-1,3,4-thiadiazole, 97 %



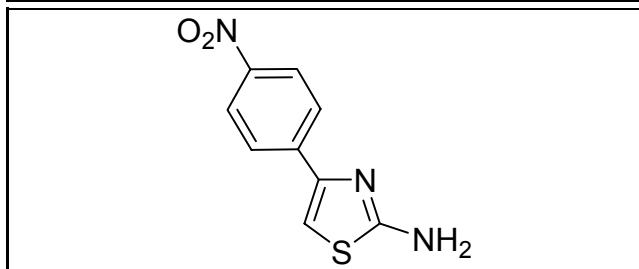
ALDRICH 653527
4-(4-Acetamidophenyl)-2-aminothiazole, 97 %



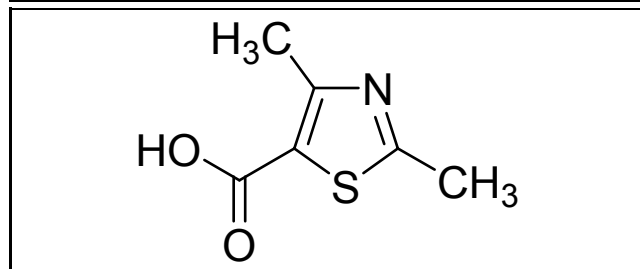
ALDRICH 632570 (2103-88-0)
4-Phenylthiazole-2-thiol, technical grade, 90 %



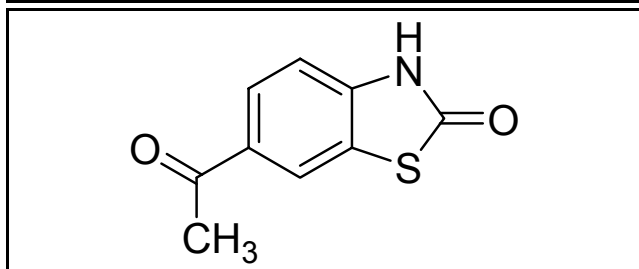
ALDRICH 658499 (2104-09-8)
2-Amino-4-(4-nitrophenyl)thiazole, 96 %



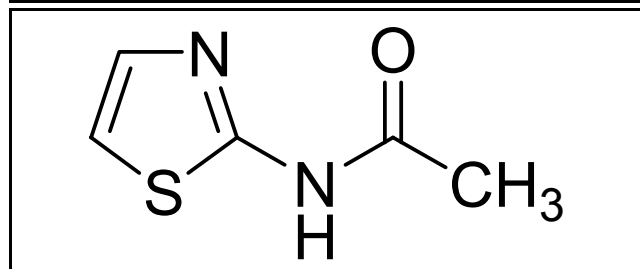
ALDRICH 642509 (53137-27-2)
2,4-Dimethylthiazole-5-carboxylic acid, 97 %



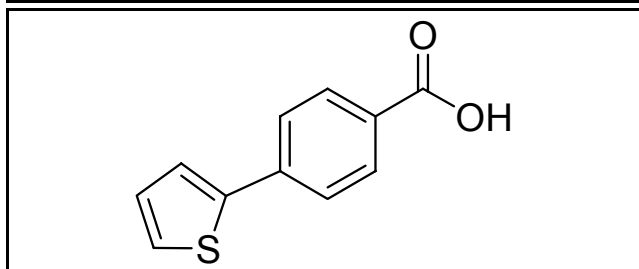
ALDRICH 663182
6-Acetyl-2(3H)-benzothiazolone, 97 %



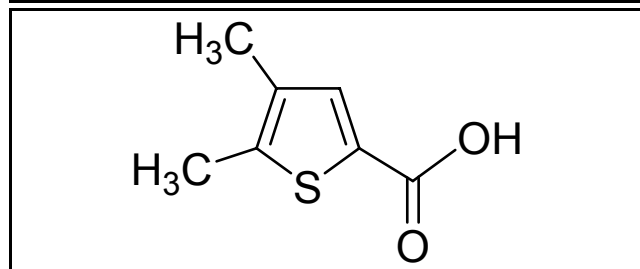
ALDRICH 652954 (2719-23-5)
2-Acetamidothiazole, purum, 98 %



ALDRICH 654566
4-(Thiophen-2-yl)benzoic acid, 97 %

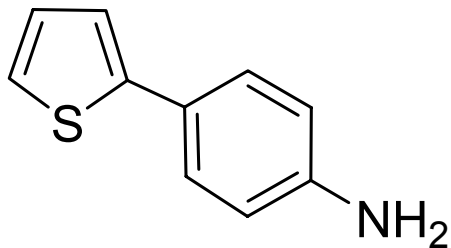


ALDRICH 668346 (40808-24-0)
4,5-Dimethylthiophene-2-carboxylic acid, 97 %



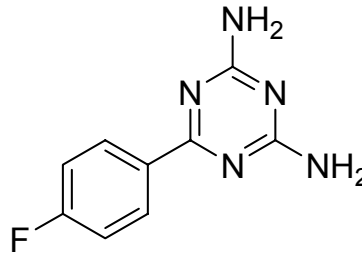
ALDRICH 654558 (70010-48-9)

4-(Thiophen-2-yl)aniline, 97 %



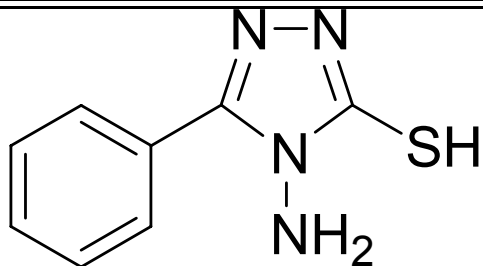
ALDRICH 597570 (30530-44-0)

2,4-Diamino-6-(4-fluorophenyl)-1,3,5-triazine, 95 %



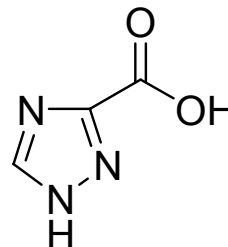
ALDRICH 675008 (22706-11-2)

4-Amino-5-phenyl-4H-1,2,4-triazole-3-thiol, 97 %



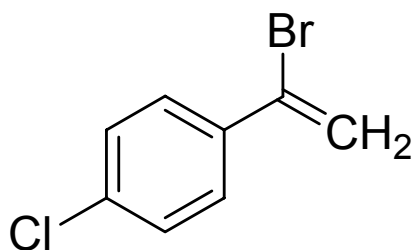
ALDRICH 658472 (4928-87-4)

1,2,4-Triazole-3-carboxylic acid, 97 %



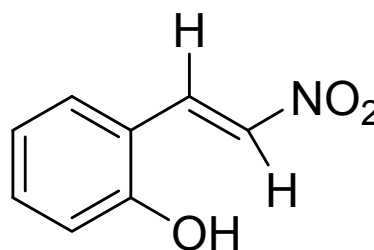
ALDRICH 675113 (89619-10-3)

1-(1-Bromovinyl)-4-chlorobenzene, 90 %



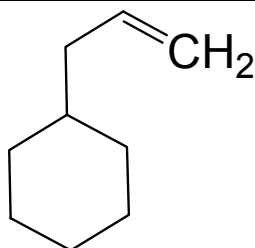
ALDRICH 639680 (3156-43-2)

trans-2-Hydroxy-beta-nitrostyrene



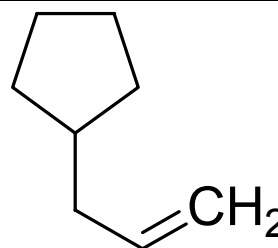
ALDRICH 662844 (2114-42-3)

Allylcyclohexane, 96 %



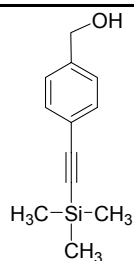
ALDRICH 662852 (3524-75-2)

Allylcyclopentane, 97 %



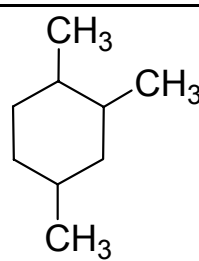
ALDRICH 668273 (275386-60-2)

4-(Trimethylsilylethynyl)benzyl alcohol, 97 %

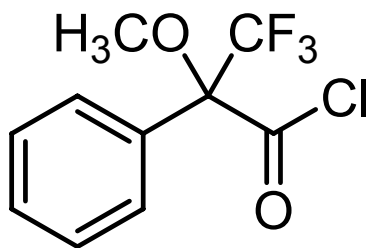


ALDRICH 677345 (2234-75-5)

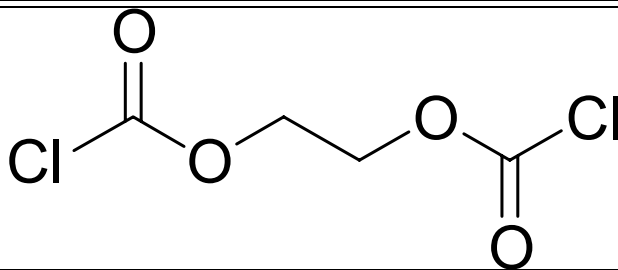
1,2,4-Trimethylcyclohexane mixture of isomers



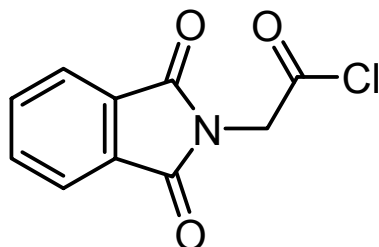
ALDRICH 667773 (64312-89-6)
alpha-Methoxy-alpha-trifluoromethylphenylacetyl chloride, 97 %



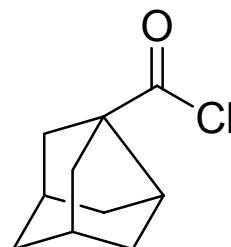
ALDRICH 653071 (124-05-0)
Ethylenebis(chloroformate), puriss., 98 %



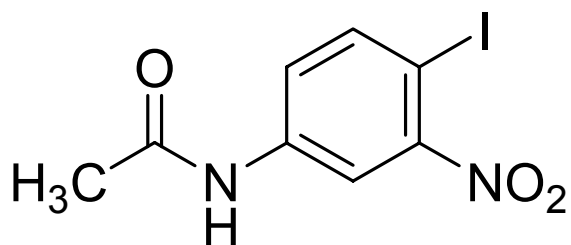
ALDRICH 636401 (6780-38-7)
Phthalylglycyl chloride



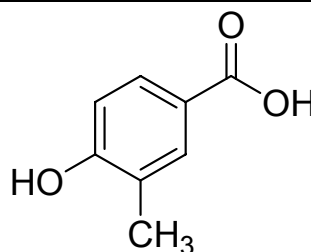
ALDRICH 678678 (67064-09-9)
3-Noradamantanecarbonyl chloride, 97 %



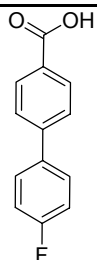
ALDRICH 667692
N-(4-Iodo-3-nitrophenyl)acetamide, 97 %



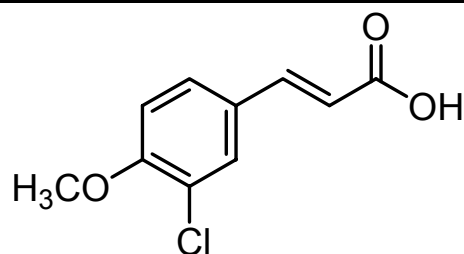
ALDRICH 659282 (499-76-3)
4-Hydroxy-3-methylbenzoic acid, 97 %



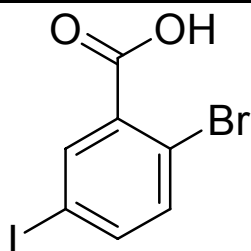
ALDRICH 651710 (5731-10-2)
4-(4-Fluorophenyl)benzoic acid, 97 %



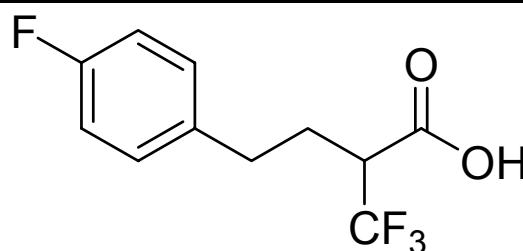
ALDRICH 638765 (58236-76-3)
3-Chloro-4-methoxycinnamic acid



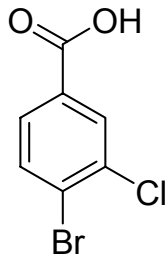
ALDRICH 665339 (25252-00-0)
2-Bromo-5-iodobenzoic acid, 97 %



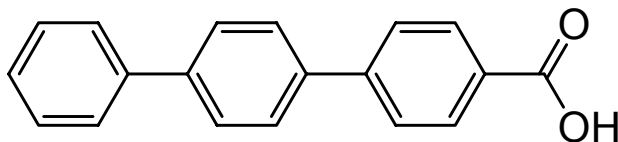
ALDRICH 674826
4-(4-Fluorophenyl)-2-(trifluoromethyl)butanoic acid



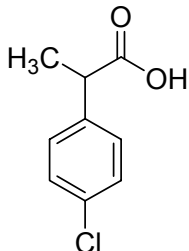
ALDRICH 665312 (25118-59-6)
4-Bromo-3-chlorobenzoic acid, 97 %



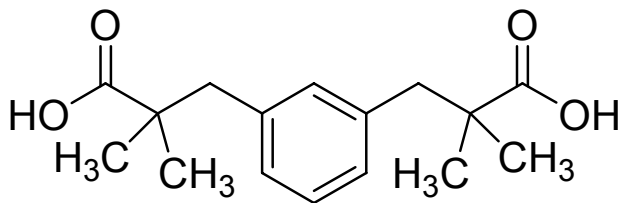
ALDRICH 674168 (5731-15-7)
p-Terphenyl-4-carboxylic acid, 97 %



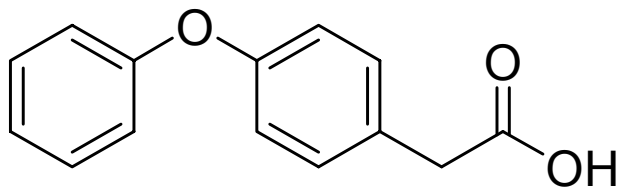
ALDRICH 650714 (938-95-4)
DL-4-Chloro-alpha-methylphenylacetic acid, 96 %



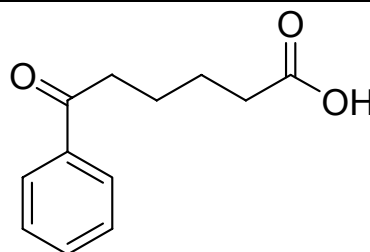
ALDRICH 666122
alpha,alpha,alpha',alpha'-Tetramethyl-1,3-benzenedipropionic acid, 97 %



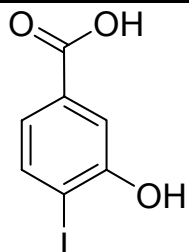
ALDRICH 665053 (6328-74-1)
4-Phenoxyphenylacetic acid, 97 %



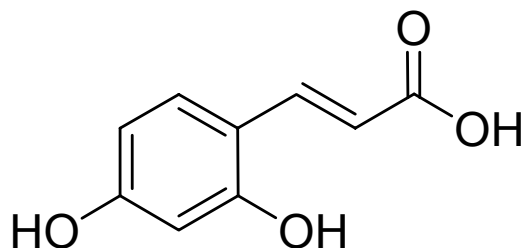
ALDRICH 663166 (4144-62-1)
5-Benzoylpentanoic acid, 97 %



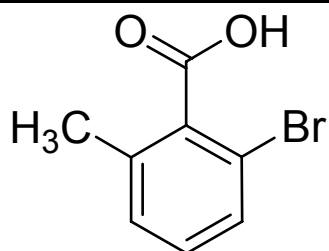
ALDRICH 664057 (58123-77-6)
3-Hydroxy-4-iodobenzoic acid, 97 %



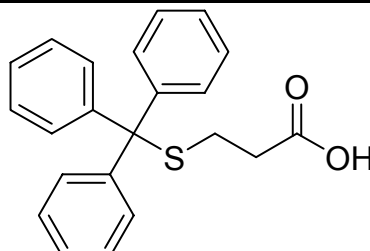
ALDRICH 663158
2,4-Dihydroxycinnamic acid, 97 %



ALDRICH 665673 (90259-31-7)
2-Bromo-6-methylbenzoic acid, 97 %

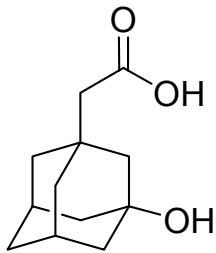


FLUKA 53695 (27144-18-9)
3-(Tritylthio)propionic acid, purum, >= 97.0 % T



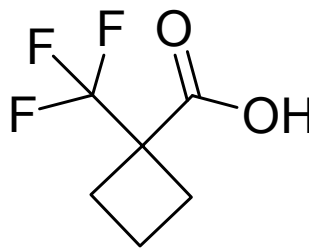
ALDRICH 678651 (17768-36-4)

3-Hydroxyadamantane-1-acetic acid, 97 %



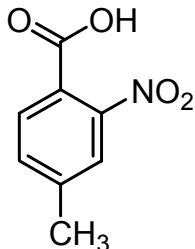
ALDRICH 665592 (277756-45-3)

1-(Trifluoromethyl)cyclobutanecarboxylic acid, 97 %



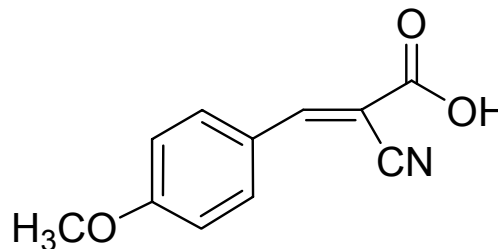
ALDRICH 638293 (27329-27-7)

4-Methyl-2-nitrobenzoic acid, 97 %



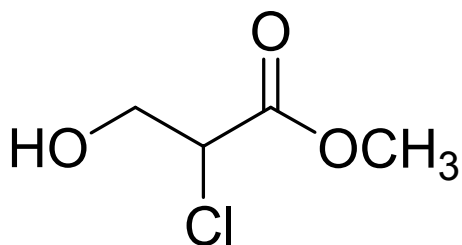
ALDRICH 586536 (1519-55-7)

alpha-Cyano-4-methoxycinnamic acid, 97 %



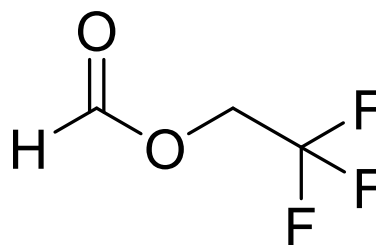
ALDRICH 678724 (98070-39-4)

Methyl 2-chloro-3-hydroxypropionate, 97 %



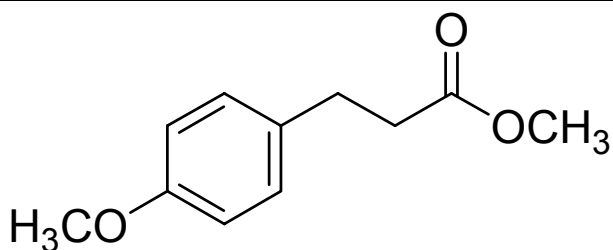
ALDRICH 669083 (32042-38-9)

2,2,2-Trifluoroethyl formate, 95 %



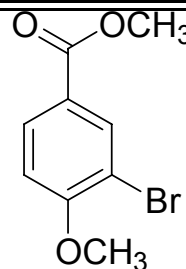
ALDRICH 666424 (15823-04-8)

Methyl 3-(4-methoxyphenyl)propionate, 97 %



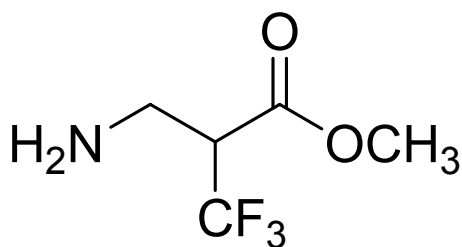
ALDRICH 652830 (35450-37-4)

Methyl 3-bromo-4-methoxybenzoate, puriss., 98 %



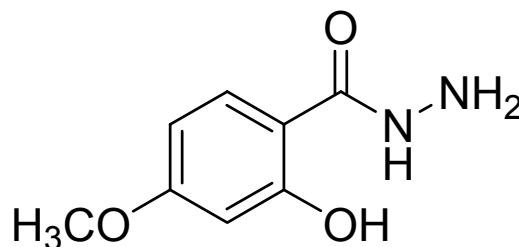
ALDRICH 666270

Methyl 3-amino-2-(trifluoromethyl)propionate, 90 %



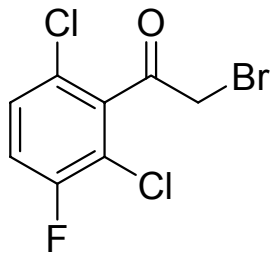
ALDRICH 538574

2-Hydroxy-4-methoxybenzoic acid hydrazide, 97 %



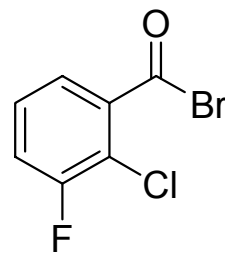
ALDRICH 669415

2-Bromo-2',6'-dichloro-3'-fluoroacetophenone, 95 %



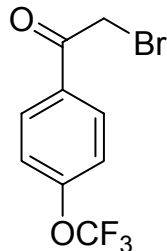
ALDRICH 669423 (63529-30-6)

2-Bromo-3'-chloro-4'-fluoroacetophenone, 96 %



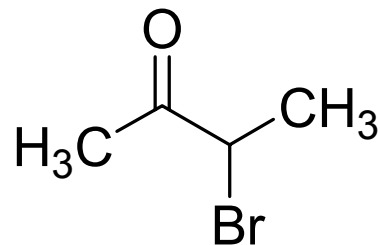
ALDRICH 669474 (103962-10-3)

2-Bromo-4'-(trifluoromethoxy)acetophenone, 97 %



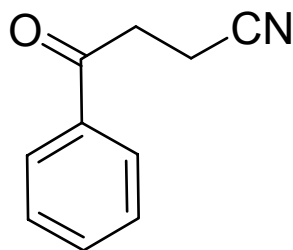
ALDRICH 662747 (814-75-5)

3-Bromo-2-butanone, 97 %



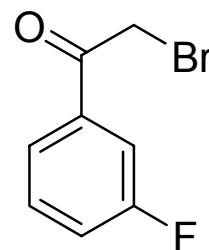
ALDRICH 667269 (5343-98-6)

4-Oxo-4-phenylbutanenitrile, 97 %



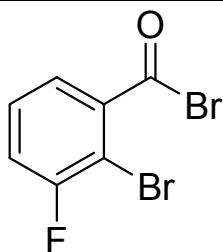
ALDRICH 669482 (53631-18-8)

2-Bromo-3'-fluoroacetophenone, 97 %



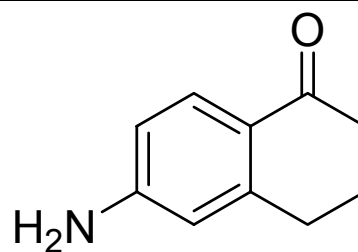
ALDRICH 669458 (435273-49-7)

2,3'-Dibromo-4'-fluoroacetophenone, 97 %



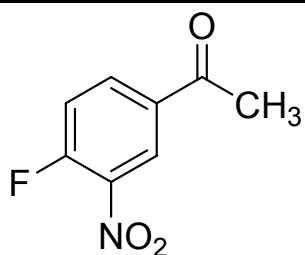
ALDRICH 678422 (3470-53-9)

6-Amino-3,4-dihydro-1(2H)-naphthalenone, 96 %



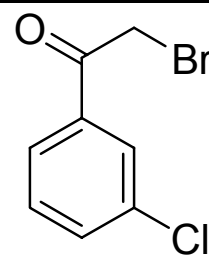
ALDRICH 676357 (400-93-1)

4'-Fluoro-3'-nitroacetophenone, 97 %



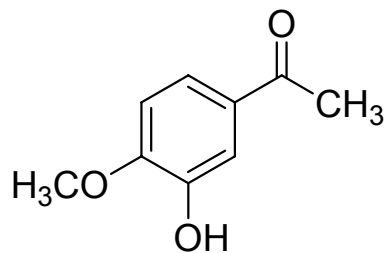
ALDRICH 672823 (41011-01-2)

2-Bromo-3'-chloroacetophenone, 97 %



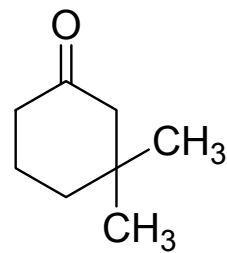
ALDRICH 679704 (6100-74-9)

3'-Hydroxy-4'-methoxyacetophenone, 97 %



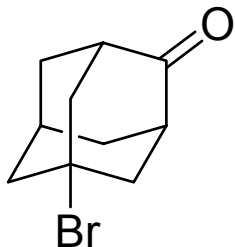
ALDRICH 674893 (2979-19-3)

3,3-Dimethylcyclohexanone, 90 %



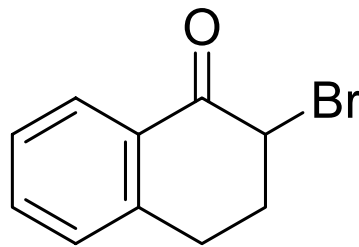
ALDRICH 678686 (20098-20-8)

5-Bromo-2-adamantanone, 97 %



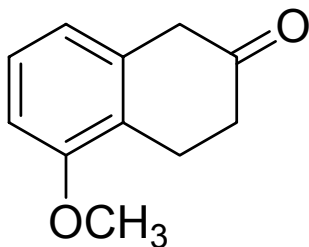
ALDRICH 667188 (13672-07-6)

2-Bromo-1-tetralone, 97 %



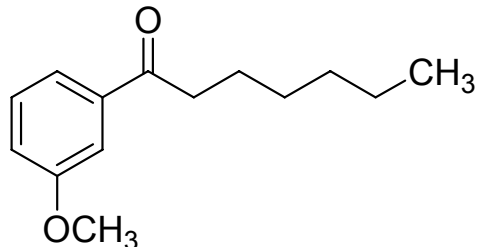
ALDRICH 579335 (32940-15-1)

5-Methoxy-2-tetralone, 97 %



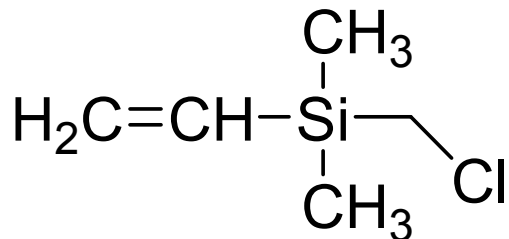
ALDRICH 649899 (100863-37-4)

1-(3-Methoxyphenyl)-1-heptanone



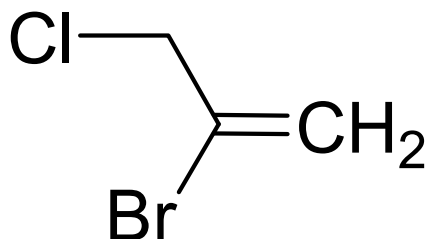
ALDRICH 649112 (16709-86-7)

(Chloromethyl)dimethylvinylsilane, purum



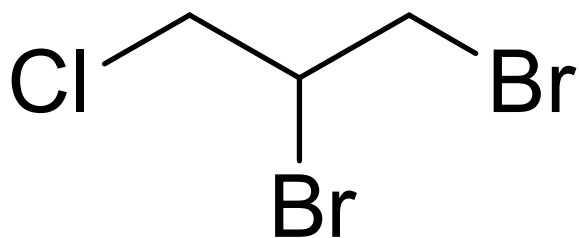
ALDRICH 679844 (16400-63-8)

2-Bromo-3-chloropropene, 97 %



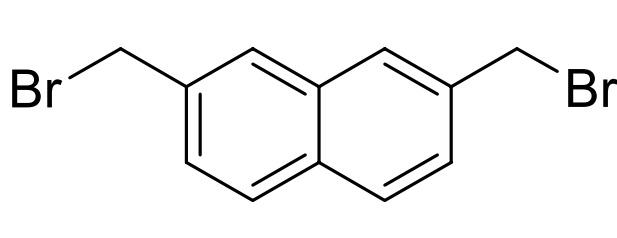
ALDRICH 676713 (96-12-8)

1,2-Dibromo-3-chloropropane, 97 %



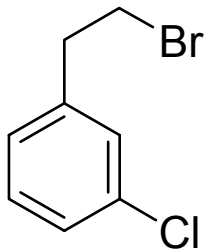
FLUKA 53658 (38309-89-6)

2,7-Bis(bromomethyl)naphthalene, purum, >= 97.0 % HPLC



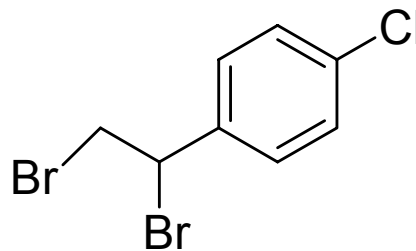
ALDRICH 655058 (16799-05-6)

3-Chlorophenethyl bromide



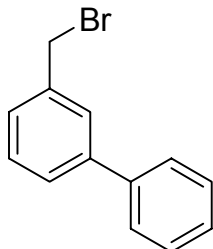
ALDRICH 674982 (23135-16-2)

1-Chloro-4-(1,2-dibromoethyl) benzene, 97 %



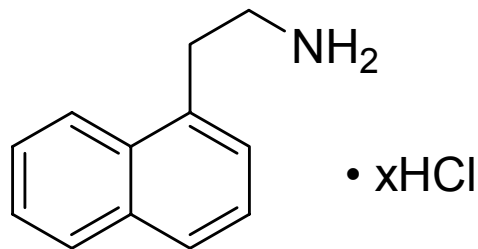
ALDRICH 538388 (14704-31-5)

3-Phenylbenzyl bromide, 97 %



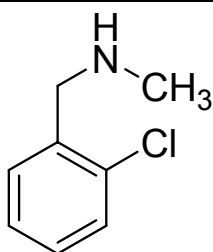
ALDRICH 661694 (942-05-2)

2-(1-Naphthyl)ethylamine hydrochloride, 95 %



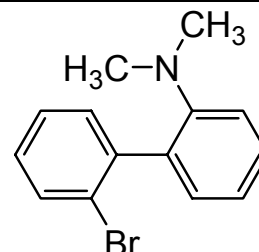
ALDRICH 126853 (94-64-4)

2-Chloro-N-methylbenzylamine, 97 %



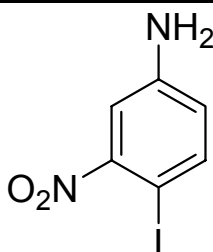
ALDRICH 642240 (213697-67-7)

2-(Dimethylamino)-2'-bromo-1,1'-biphenyl, 97 %



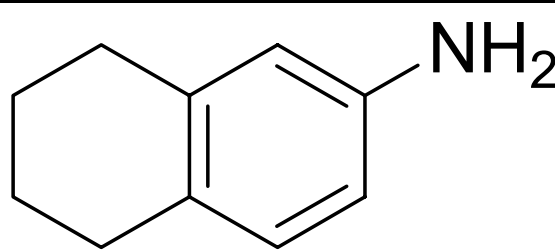
ALDRICH 667927 (105752-04-3)

4-Iodo-3-nitroaniline, 97 %



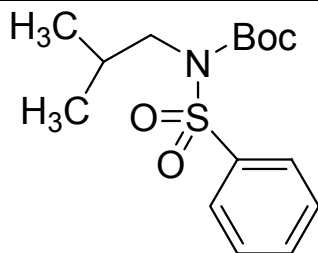
ALDRICH 678864 (2217-43-8)

5,6,7,8-Tetrahydro-2-naphthylamine, 96 %



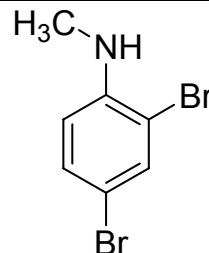
ALDRICH 679550

N-Boc-alpha-(phenylsulfonyl)isobutylamine, 97 %

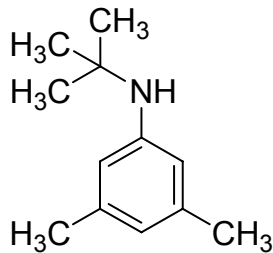


ALDRICH 665886 (73557-58-1)

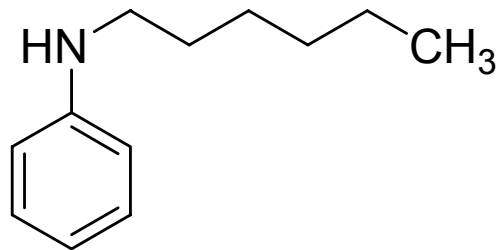
2,4-Dibromo-N-methylaniline, 97 %



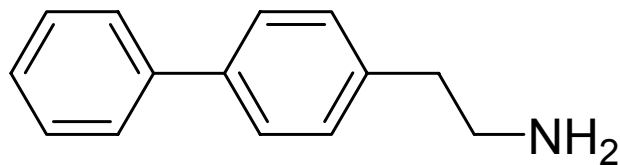
ALDRICH 677981 (110993-40-3)
N-tert-Butyl-3,5-dimethylaniline, 97 %



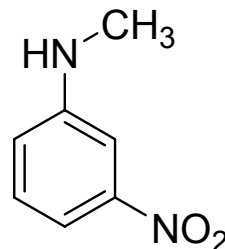
ALDRICH 673285 (4746-32-1)
N-Hexylaniline, 98 %



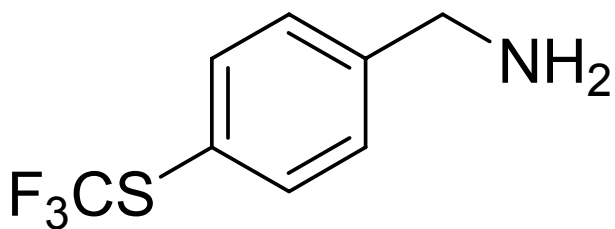
ALDRICH 560227 (17027-51-9)
2-(4-Biphenyl)ethylaniline, 98 %



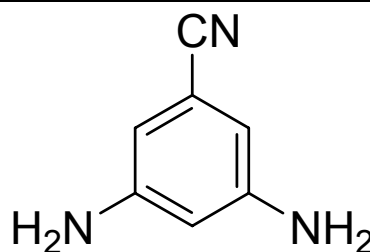
ALDRICH 665894 (619-26-1)
3-Nitro-N-methylaniline, 97 %



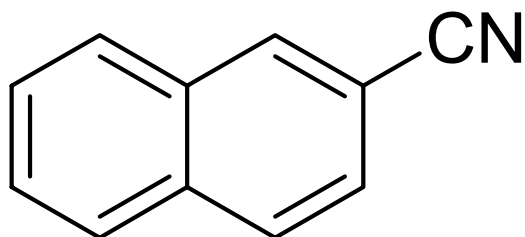
ALDRICH 555460 (128273-56-3)
4-(Trifluoromethylthio)benzylamine



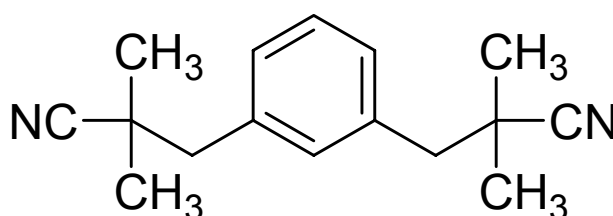
ALDRICH 669202 (33786-93-5)
3,5-Diaminobenzonitrile, 97 %



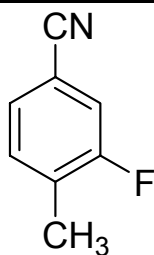
ALDRICH 673137 (613-46-7)
Naphthalene-2-carbonitrile, 97 %



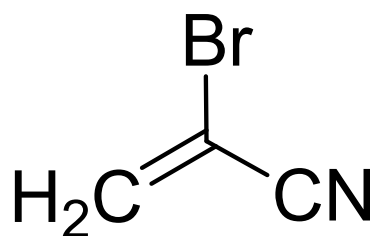
ALDRICH 666114
alpha,alpha,alpha'-alpha'-Tetramethyl-1,3-benzenedipropionitrile, 97 %



ALDRICH 664030 (170572-49-3)
3-Fluoro-4-methylbenzonitrile, 97 %

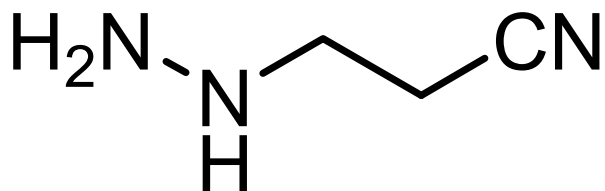


ALDRICH 674532 (920-34-3)
2-Bromoacrylonitrile, 96 %



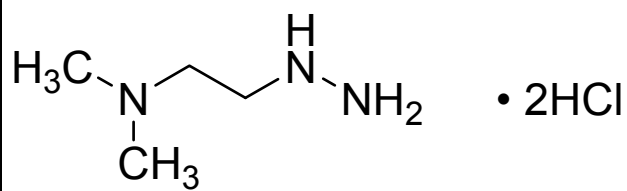
ALDRICH 636193 (353-07-1)

Cyanoethylhydrazine, 97 %



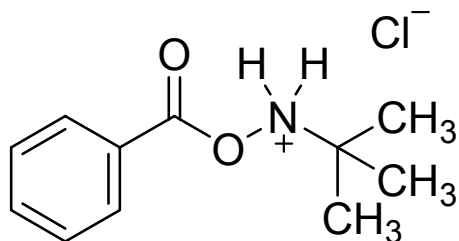
ALDRICH 674141

Dimethylaminoethylhydrazine dihydrochloride, 97 %



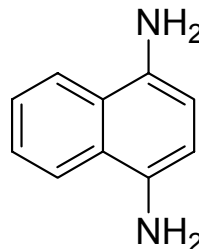
ALDRICH 669407

N-tert-Butyl-O-benzoylhydroxylamine hydrochloride, 97 %



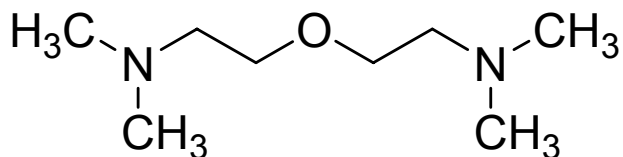
ALDRICH 667617 (2243-61-0)

1,4-Diaminonaphthalene, 97 %



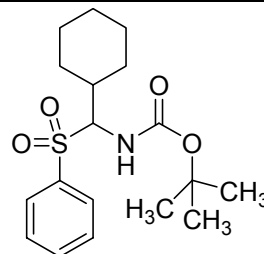
ALDRICH 667609 (3033-62-3)

Bis[2-(N,N-dimethylamino)ethyl] ether, 97 %



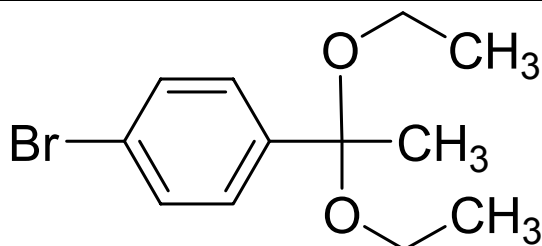
ALDRICH 669091 (253426-70-9)

N-Boc-alpha-(phenylsulfonyl)cyclohexylmethylamine, 97 %



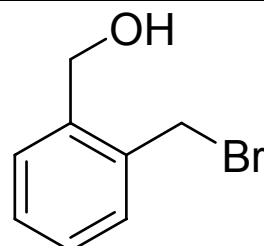
ALDRICH 568597 (61390-40-7)

4-Bromoacetophenone diethyl acetal, 95 %



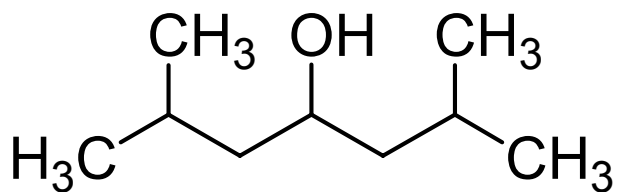
ALDRICH 667706 (74785-02-7)

2-(Bromomethyl)benzyl alcohol, 95 %



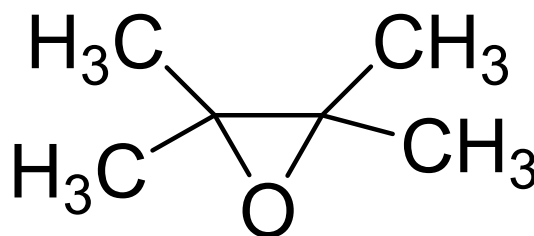
ALDRICH 541486 (108-82-7)

2,6-Dimethyl-4-heptanol, 80 %

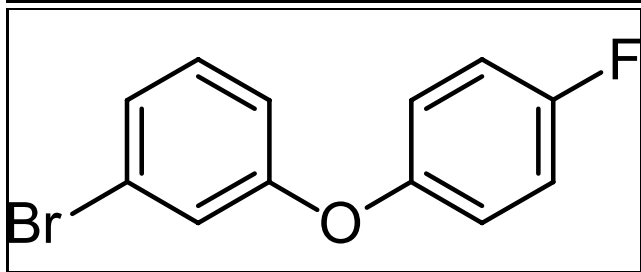


ALDRICH 662100 (5076-20-0)

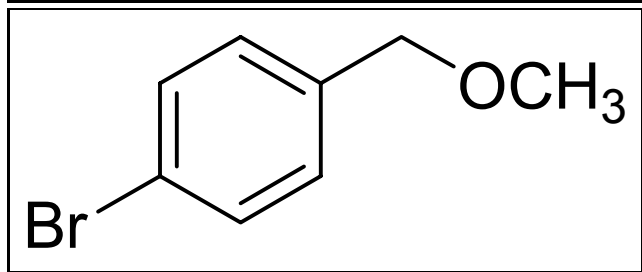
2,3-Dimethyl-2,3-epoxybutane, 99 %



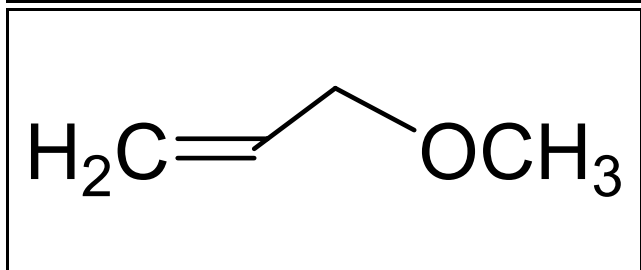
ALDRICH 636606 (50904-38-6)
3-Bromo-4'-fluorodiphenyl ether, 97 %



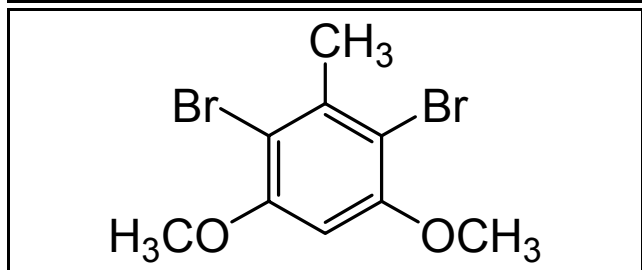
ALDRICH 680419 (1515-88-4)
4-Bromobenzyl methyl ether, 97 %



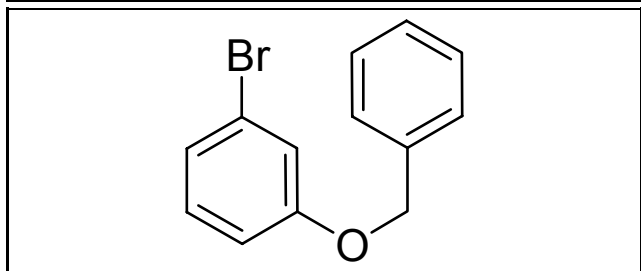
ALDRICH 678201 (627-40-7)
Allyl methyl ether, 97 %



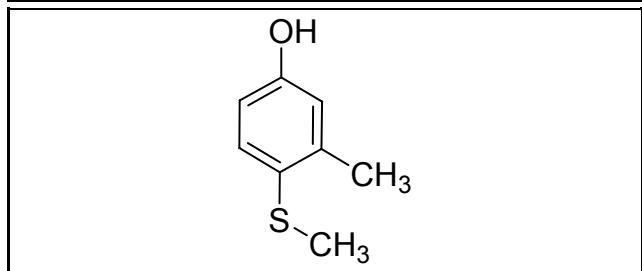
ALDRICH 677752 (13365-08-7)
2,6-Dibromo-3,5-dimethoxytoluene, 97 %



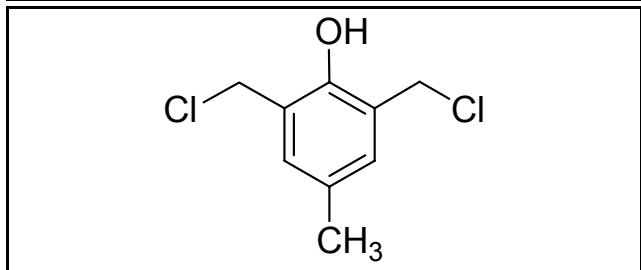
ALDRICH 668419 (53087-13-1)
3-Benzyloxybromobenzene, 96 %



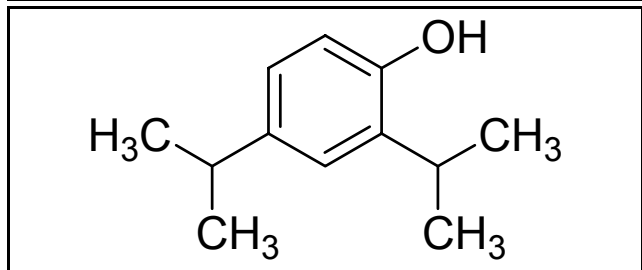
ALDRICH 673145 (3120-74-9)
3-Methyl-4-(methylthio)phenol, 97 %



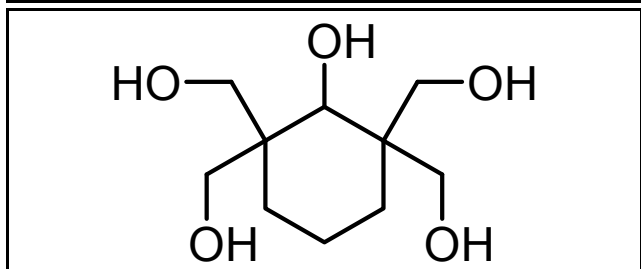
ALDRICH 668613 (5862-32-8)
2,6-Bis(chloromethyl)-4-methylphenol, 97 %



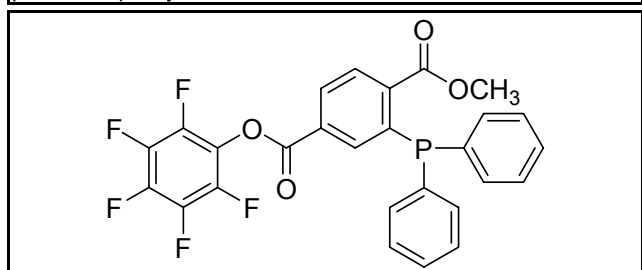
ALDRICH 557811 (2934-05-6)
2,4-Diisopropylphenol, 98 %



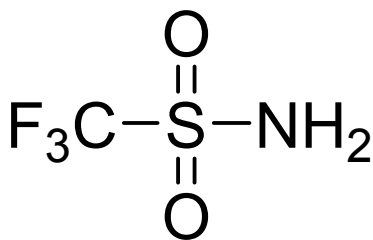
ALDRICH 632082 (5416-55-7)
2,2,6,6-Tetrakis(hydroxymethyl)cyclohexanol, 97 %



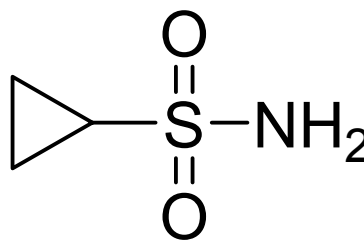
ALDRICH 679011
2-(Diphenylphosphino)terephthalic acid, 1-methyl 4-pentafluorophenyldiester, 97 %



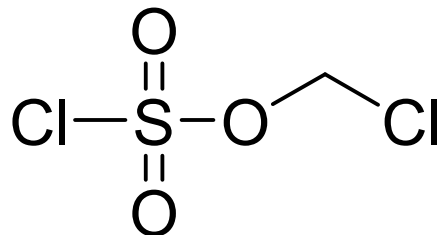
ALDRICH 638455 (421-85-2)
Trifluoromethanesulfonamide, 95 %



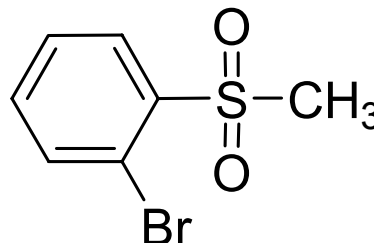
ALDRICH 674060 (154350-28-4)
Cyclopropanesulfonamide



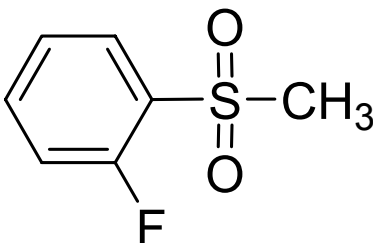
ALDRICH 656682 (49715-04-0)
Chloromethyl chlorosulfate, 97 %



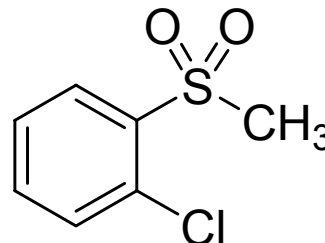
ALDRICH 669334 (33951-33-6)
2-Bromophenyl methyl sulfone, 97 %



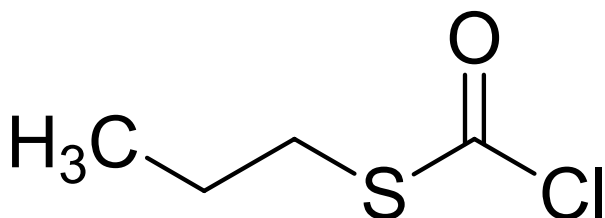
ALDRICH 669385 (654-47-7)
2-Fluorophenyl methyl sulfone, 97 %



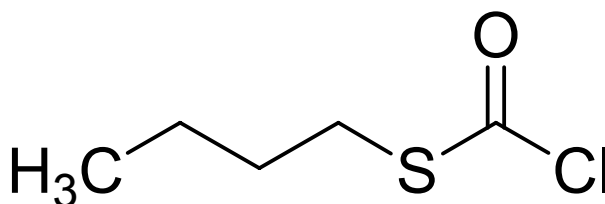
ALDRICH 669342 (17482-05-2)
2-Chlorophenyl methyl sulfone, 97 %



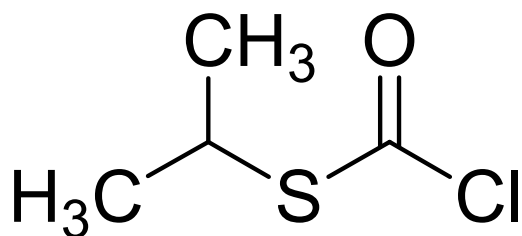
ALDRICH 668788 (13889-92-4)
S-Propyl chlorothioformate, 96 %



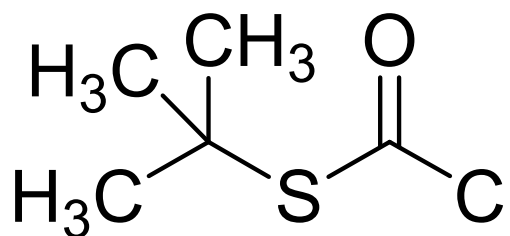
ALDRICH 672912 (13889-94-6)
S-Butyl chlorothioformate, 96 %



ALDRICH 674184 (13889-93-5)
S-Isopropyl chlorothioformate, 96 %

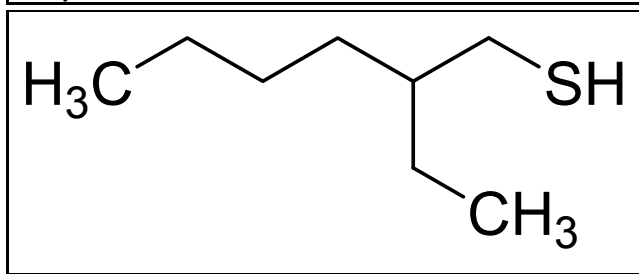


ALDRICH 672939 (13889-95-7)
S-tert-Butyl chlorothioformate, 96 %



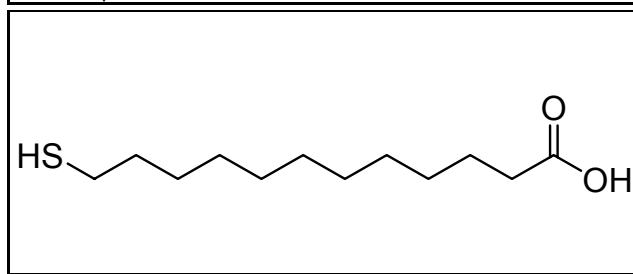
ALDRICH 669148 (7341-17-5)

2-Ethylhexanethiol, 97 %



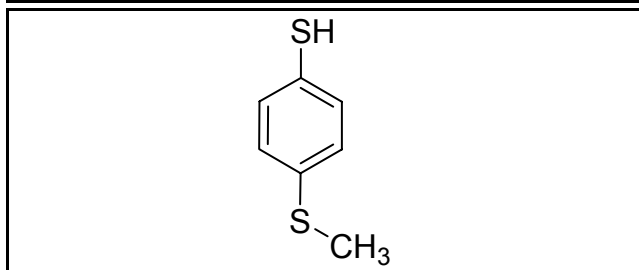
ALDRICH 675067 (82001-53-4)

12-Mercaptododecanoic acid, 96 %



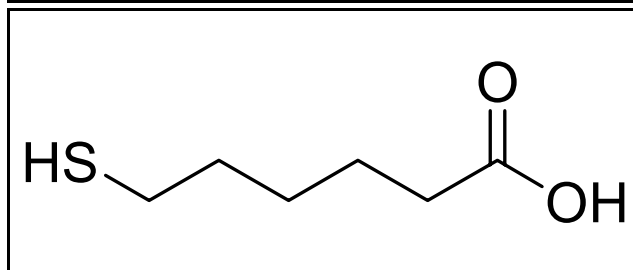
ALDRICH 662496 (1122-97-0)

4-(Methylsulfanyl)thiophenol, 97 %



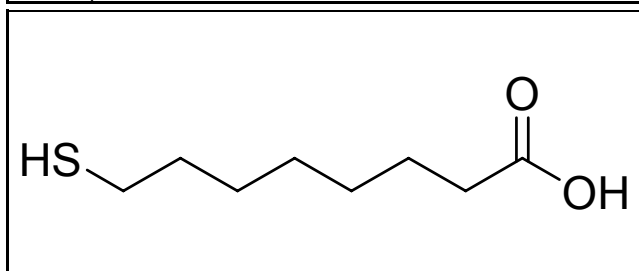
ALDRICH 674974 (17689-17-7)

6-Mercaptohexanoic acid, 90 %



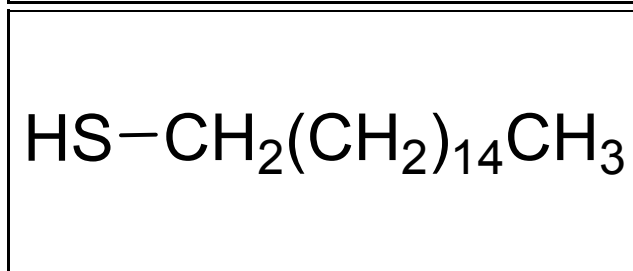
ALDRICH 675075 (74328-61-3)

8-Mercptooctanoic acid, 96 %



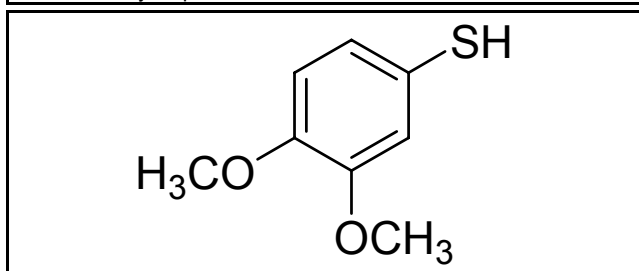
ALDRICH 674516 (2917-26-2)

1-Hexadecanethiol, 99 %



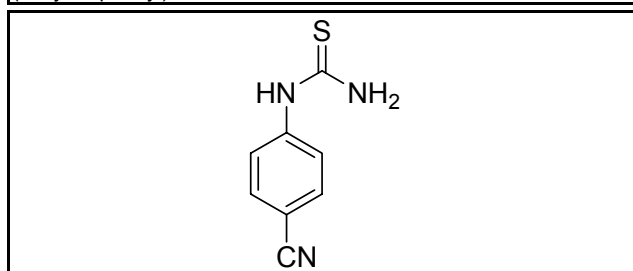
ALDRICH 662518 (700-96-9)

3,4-Dimethoxythiophenol, 97 %



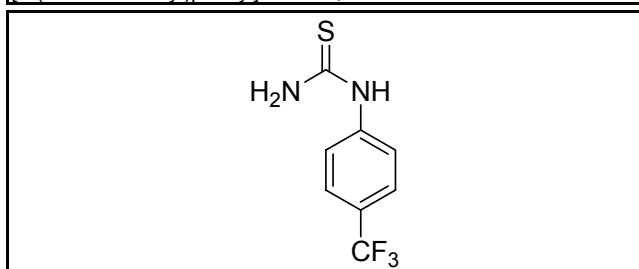
ALDRICH 654744 (3460-55-7)

(4-Cyanophenyl)thiourea, 97 %



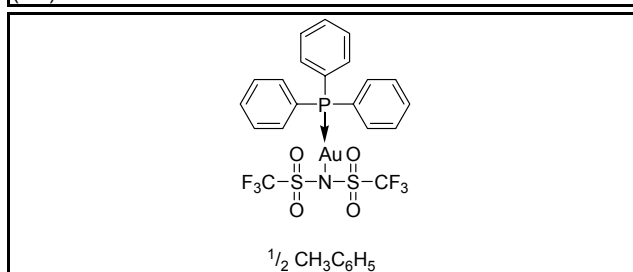
ALDRICH 654760 (1736-72-7)

[4-(Trifluoromethyl)phenyl]thiourea, 97 %



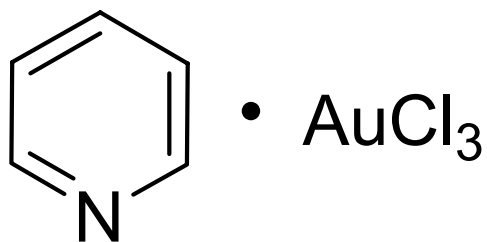
ALDRICH 677922

[Bis(trifluoromethanesulfonyl)imidate](triphenylphosphine)gold(I)
(2:1) toluene adduct



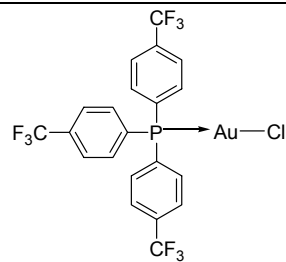
ALDRICH 677876 (14911-01-4)

Trichloro(pyridine)gold(III), 97 %



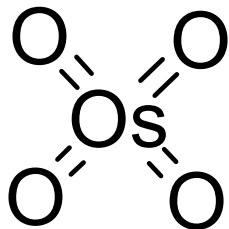
ALDRICH 665177 (385815-83-8)

Chloro(tris(para-trifluoromethylphenyl)phosphine)gold(I), 99 %



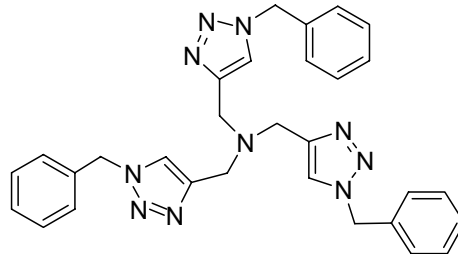
ALDRICH 658685 (20816-12-0)

Os EnCat(R), loading 0.3mmol/g Os



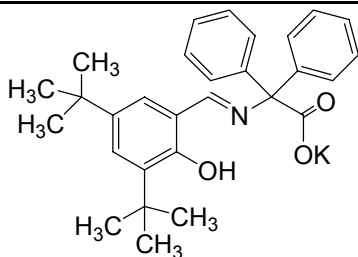
ALDRICH 678937

Tris[(1-benzyl-1H-1,2,3-triazol-4-yl)methyl]amine, 97 %



ALDRICH 676551

Potassium 2-(3,5-di-tert-butyl-2-hydroxybenzylideneamino)-2,2-diphenylacetate, 95 %



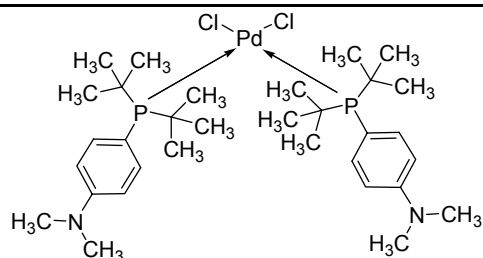
ALDRICH 674133

Palladium nanoparticles entrapped in aluminum hydroxide matrix

Pd/AlO(OH)

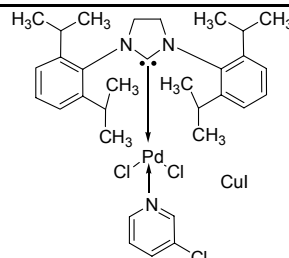
ALDRICH 678740 (887919-35-9)

(AMPHOS)2PdCl2



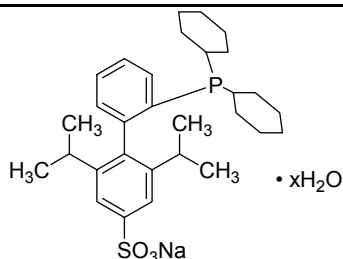
ALDRICH 677256

PEPPSI(R)-SONO



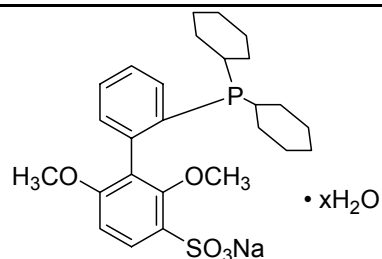
ALDRICH 677272

Sodium 2'-(dicyclohexylphosphino)-2,6-diisopropyl-1,1'-biphenyl-4-sulfonate hydrate



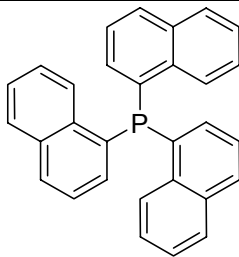
ALDRICH 677280

Sodium 2'-(dicyclohexylphosphino)-2,6-dimethoxy-1,1'-biphenyl-3-sulfonate hydrate



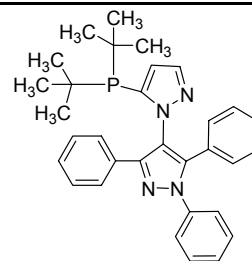
ALDRICH 667218 (3411-48-1)

Tri-1-naphthylphosphine, 97 %



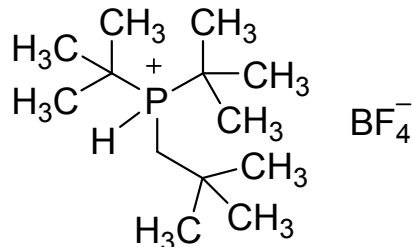
ALDRICH 676632

5-(Di-tert-butylphosphino)-1', 3', 5'-triphenyl-1'H-[1,4']bipyrazole, 97 %



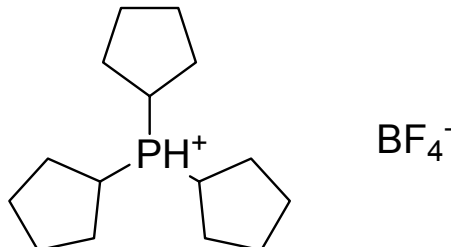
ALDRICH 676179

Di-tert-butylneopentylphosphonium tetrafluoroborate, 97 %



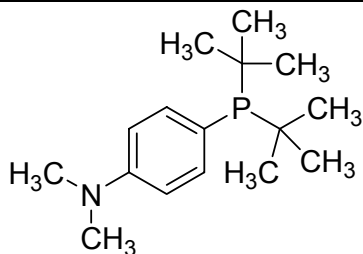
ALDRICH 673951 (610756-04-2)

Tricyclopentylphosphine tetrafluoroborate, 95 %



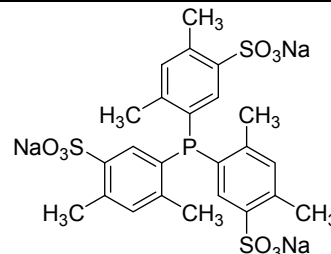
ALDRICH 677264

(4-(N,N-Dimethylamino)phenyl)di-tert-butyl phosphine, 95 %



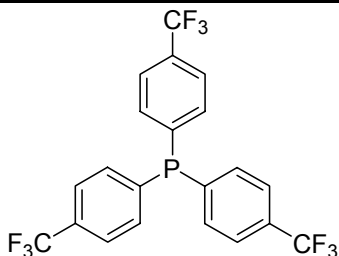
ALDRICH 667382 (443150-11-6)

Tris(2, 4-dimethyl-5-sulfophenyl)phosphine trisodium salt, 95 %



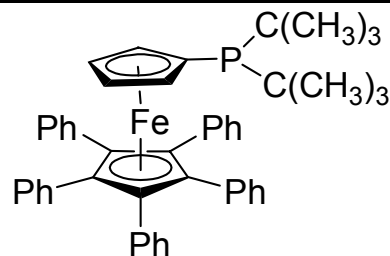
ALDRICH 666629 (13406-29-6)

Tris(4-trifluoromethylphenyl)phosphine, 97 %



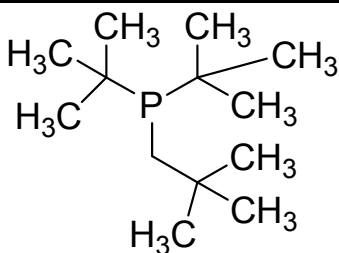
ALDRICH 675784 (312959-24-3)

1,2,3,4,5-Pentaphenyl-1'-(di-tert-butylphosphino)ferrocene



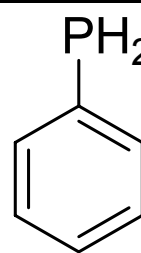
ALDRICH 676187

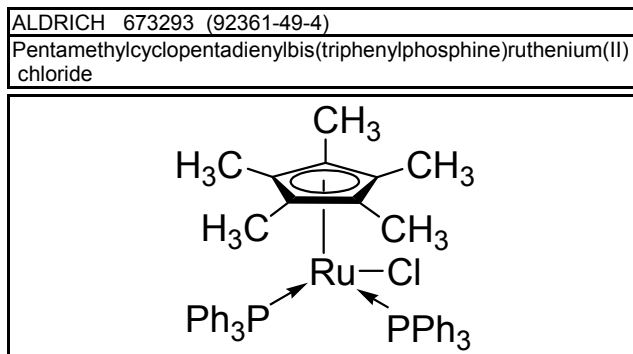
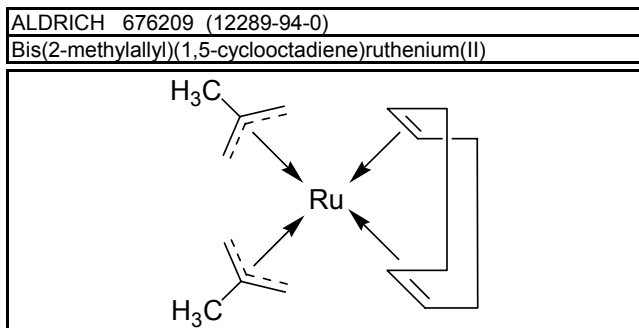
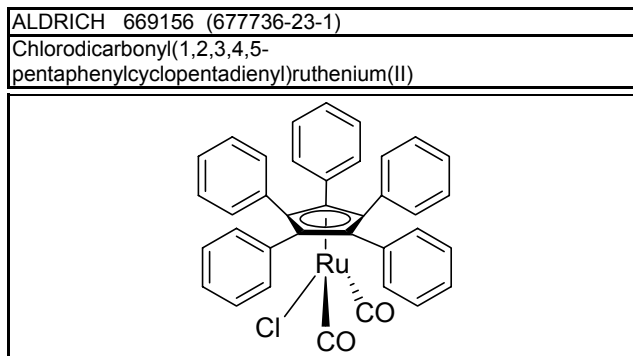
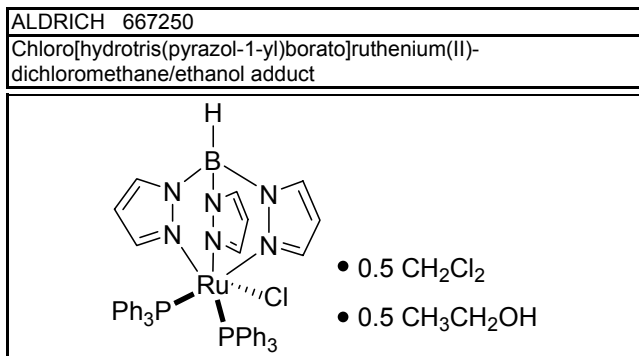
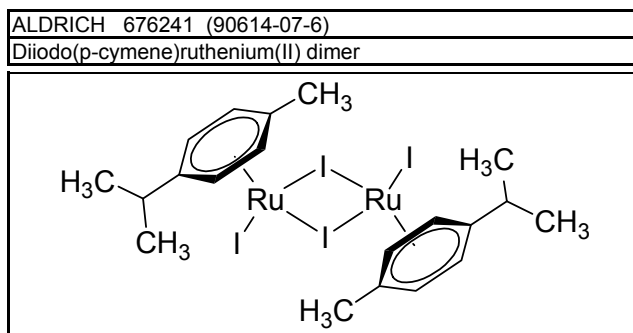
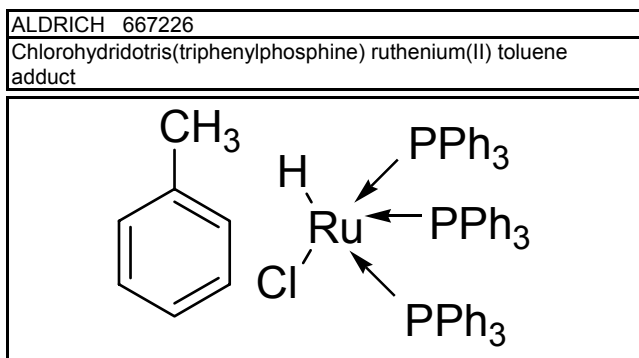
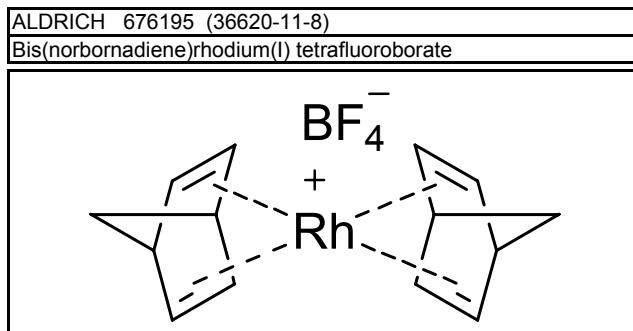
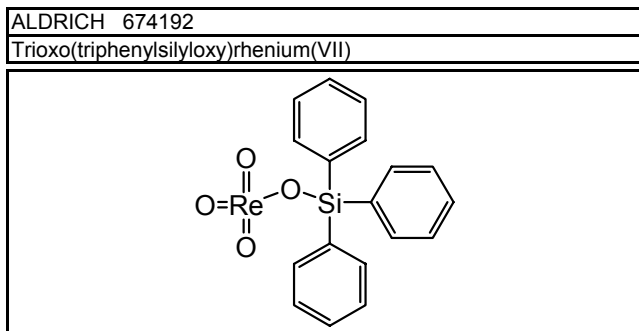
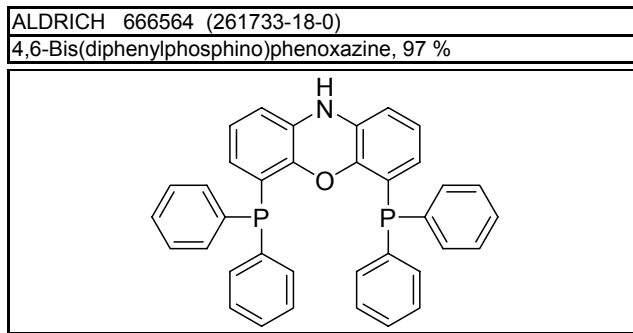
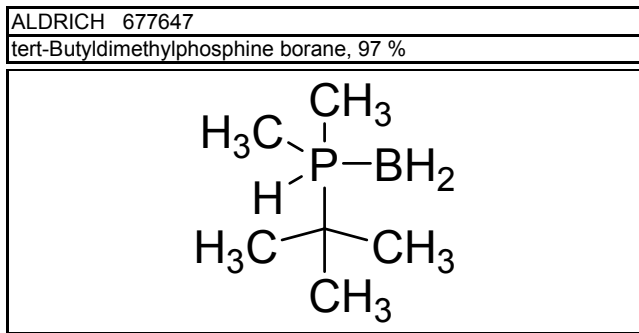
Di-tert-butylneopentylphosphine, 95 %



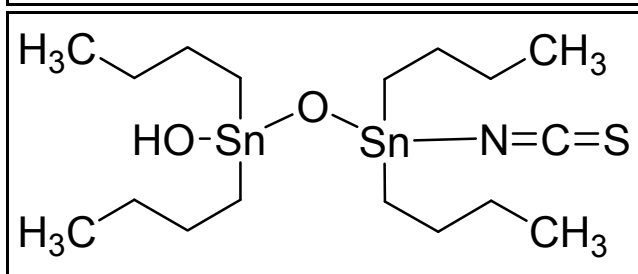
ALDRICH 674389 (638-21-1)

Phenylphosphine, ca. 10% weight in hexanes

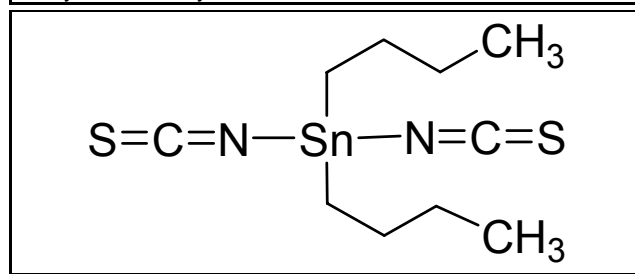




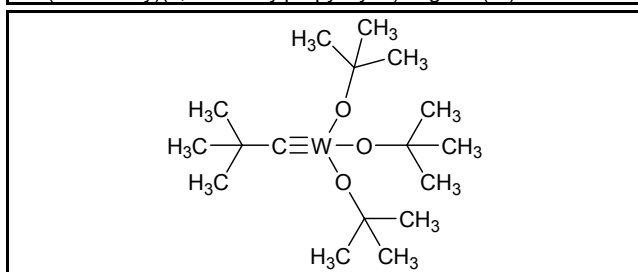
ALDRICH 674214 (27515-11-3)
1-Hydroxy-3-(isothiocyanato)-1,1,3,3-tetrabutylidistannoxane, 97 %



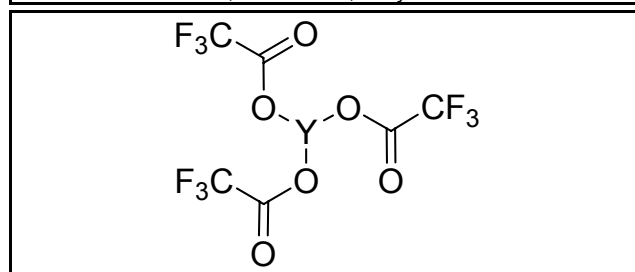
ALDRICH 669075 (15719-34-3)
Dibutyltin diisothiocyanate, 97 %



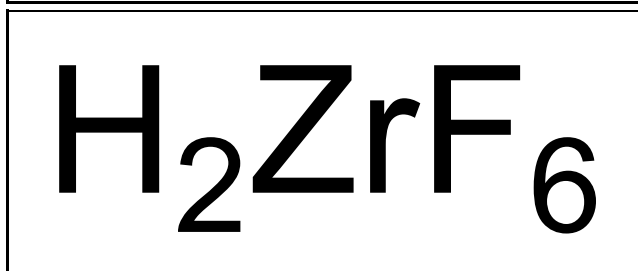
ALDRICH 666440 (78234-36-3)
Tris(tert-butoxy)(2,2-dimethylpropylidene)tungsten(VI)



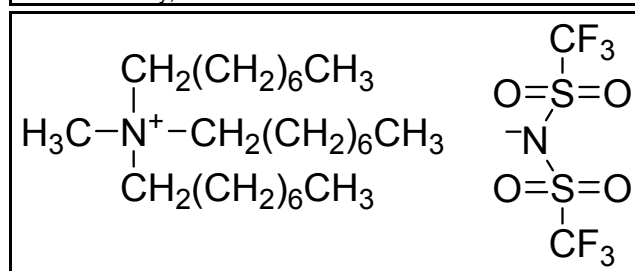
ALDRICH 659886 (304851-95-4)
Yttrium trifluoroacetate, >= 99.99 %, anhydrous



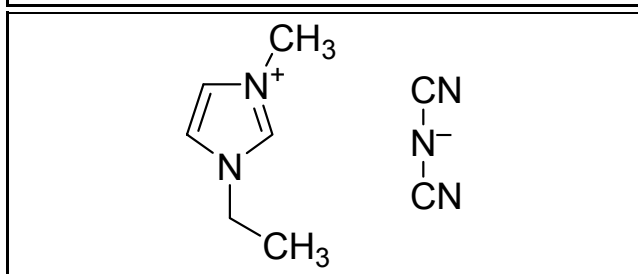
ALDRICH 679755 (12021-95-3)
Hexafluorozirconic acid solution, 50wt. % H2O



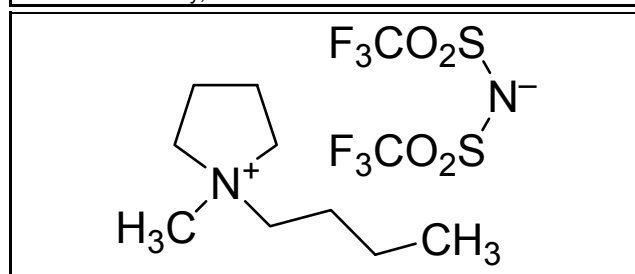
FLUKA 00797 (375395-33-8)
Methyl-trioctylammonium bis(trifluoromethylsulfonyl)imide, for electrochemistry, >= 99 % NMR



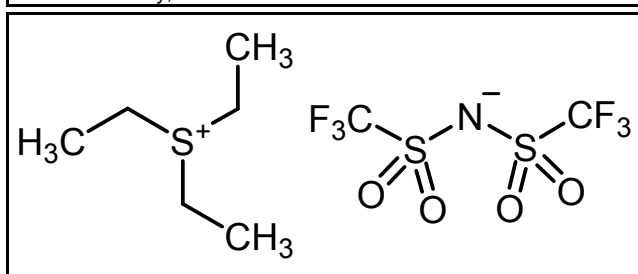
FLUKA 00796 (370865-89-7)
1-Ethyl-3-methylimidazolium dicyanamide, for electrochemistry, >= 99 % HPLC



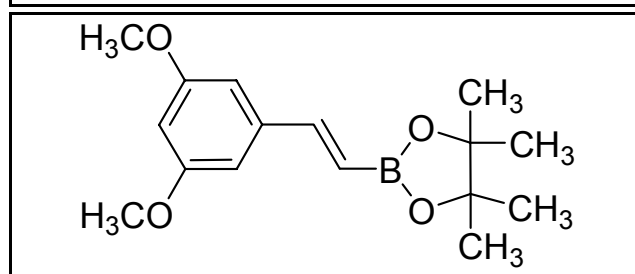
FLUKA 40963 (223437-11-4)
1-Butyl-1-methylpyrrolidinium bis(trifluoromethylsulfonyl)imide, for electrochemistry, >= 99.0 % T



FLUKA 08748 (321746-49-0)
Triethylsulfonium bis(trifluoromethylsulfonyl)imide, for electrochemistry, >= 99.0 % NMR

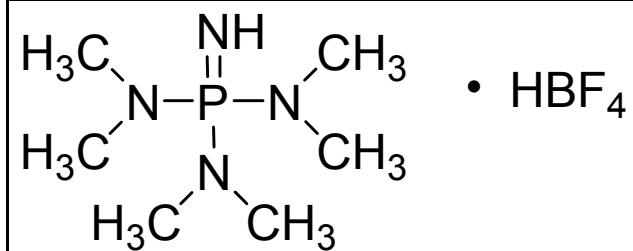


ALDRICH 676160
trans-2-(3,5-Dimethoxyphenyl)vinylboronic acid pinacol ester, 97 %



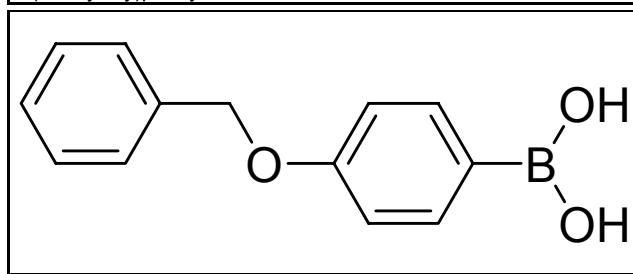
FLUKA 56815 (181470-75-7)

Iminotris(dimethylamino)phosphonium tetrafluoroborate salt,
purum, >= 95.0 % 31P-NMR



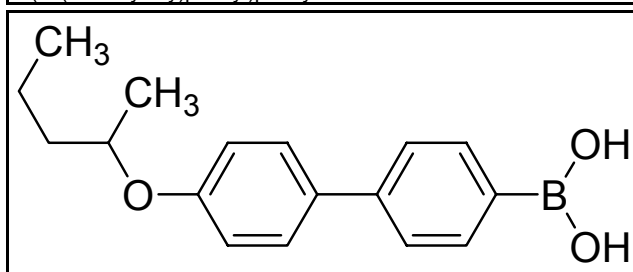
ALDRICH 666920 (146631-00-7)

4-(Benzyloxy)phenylboronic acid



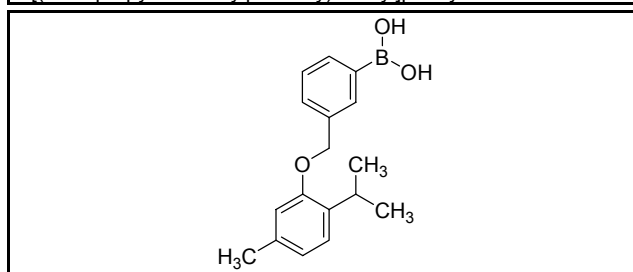
ALDRICH 666815

4-(4'-(2-Pentyl-2-methoxy)phenyl)phenylboronic acid



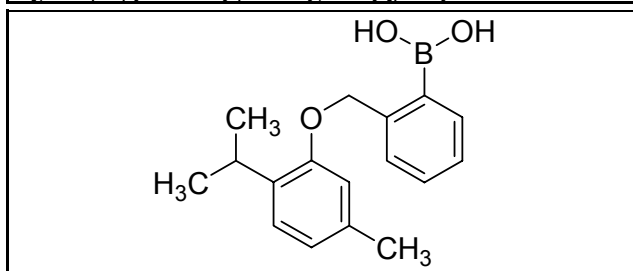
ALDRICH 661848

3-[(2'-Isopropyl-5'-methylphenoxy)methyl]phenylboronic acid



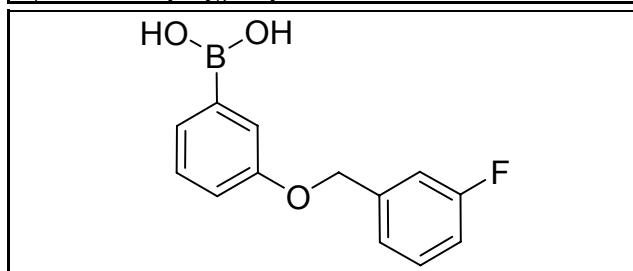
ALDRICH 661856

2-[(2'-Isopropyl-5'-methylphenoxy)methyl]phenylboronic acid



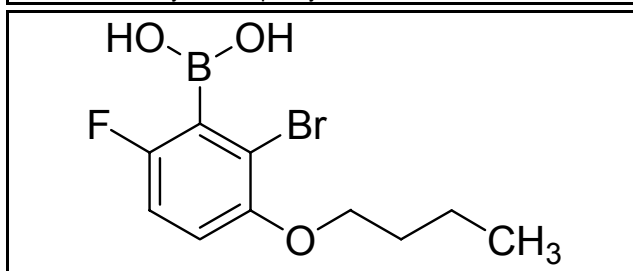
ALDRICH 661872

3-(3'-Fluorobenzoyloxy)phenylboronic acid



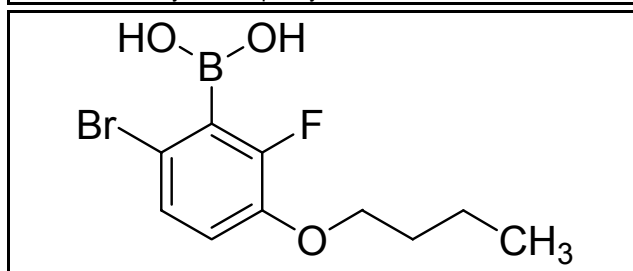
ALDRICH 661945

2-Bromo-3-butoxy-6-fluorophenylboronic acid



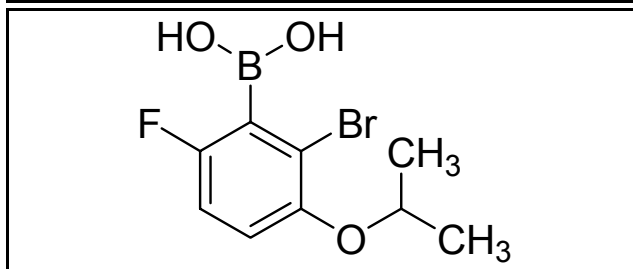
ALDRICH 661953

6-Bromo-3-butoxy-2-fluorophenylboronic acid



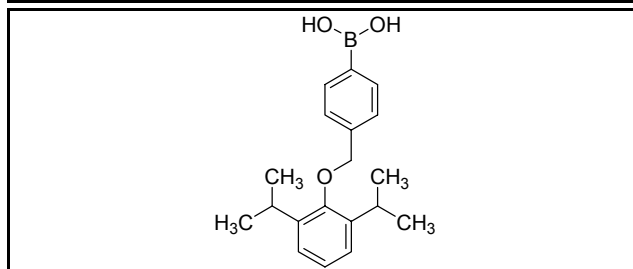
ALDRICH 661996

2-Bromo-6-fluoro-3-isopropoxyphenylboronic acid



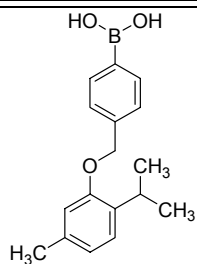
ALDRICH 662038

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



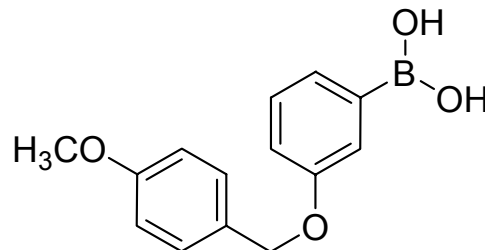
ALDRICH 662097

4-[(2-Isopropyl-5-methylphenoxy)methyl]phenylboronic acid



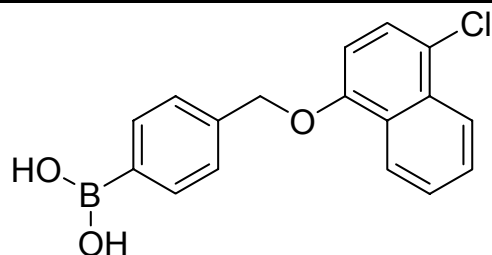
ALDRICH 666750

3-(4'-Methoxybenzyloxy)phenylboronic acid



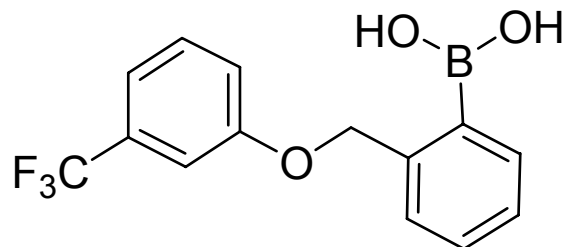
ALDRICH 645656 (870778-84-0)

4-[(4'-Chloro-1-naphthoxy)methyl]phenylboronic acid



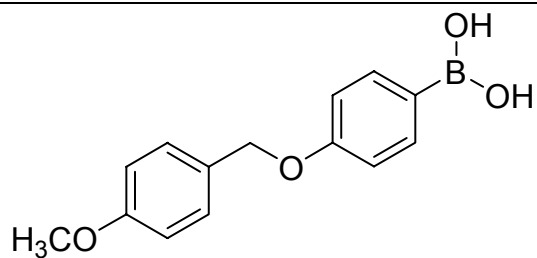
ALDRICH 666866

2-[(3'-(Trifluoromethyl)phenoxy)methyl]phenylboronic acid



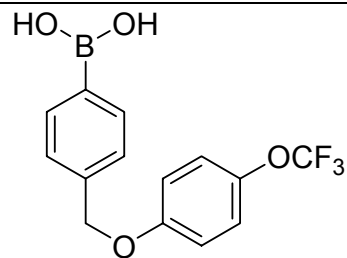
ALDRICH 666769

4-(4'-Methoxybenzyloxy)phenylboronic acid



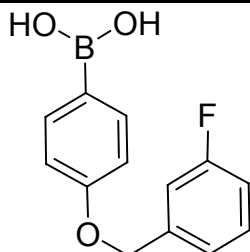
ALDRICH 651303 (870778-97-5)

4-[4'-(Trifluoromethoxy)phenoxy]methyl]phenylboronic acid



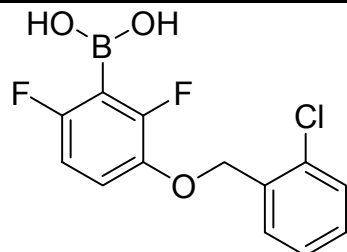
ALDRICH 661864

4-(3'-Fluorobenzyloxy)phenylboronic acid



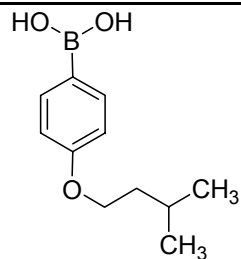
ALDRICH 651362 (870778-99-7)

2,6-Difluoro-3-(2'-chlorobenzyloxy)phenylboronic acid



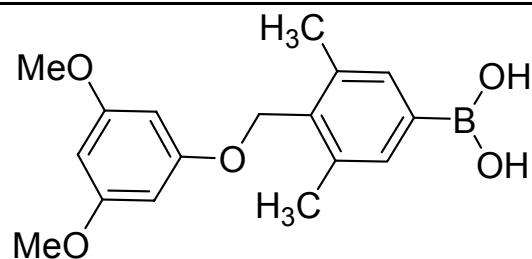
ALDRICH 662070

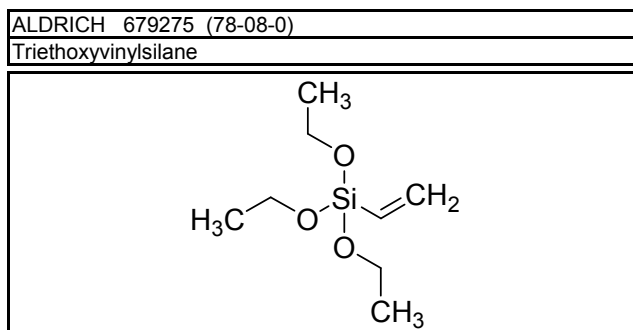
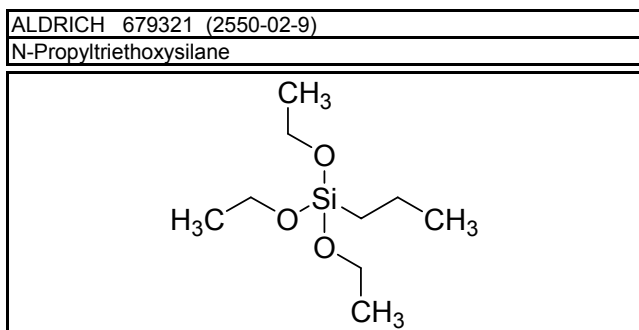
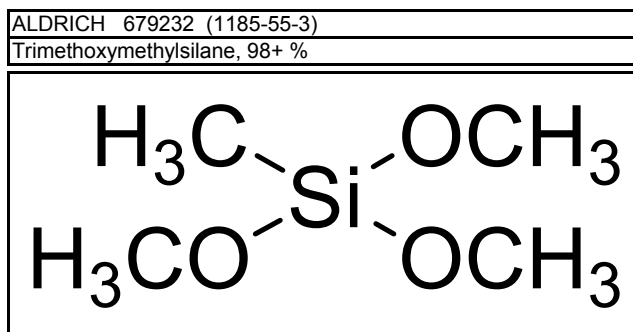
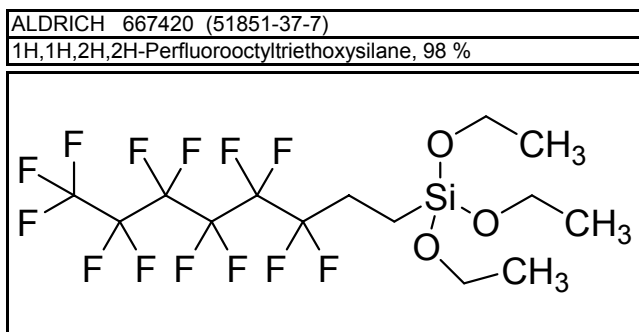
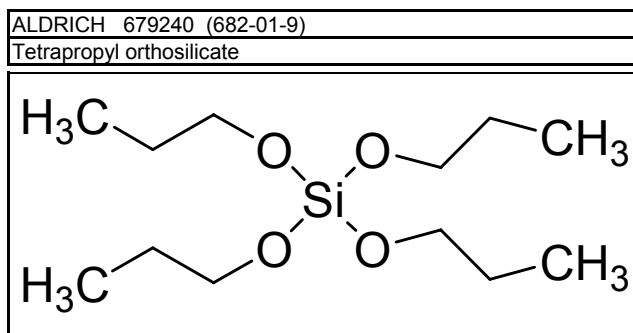
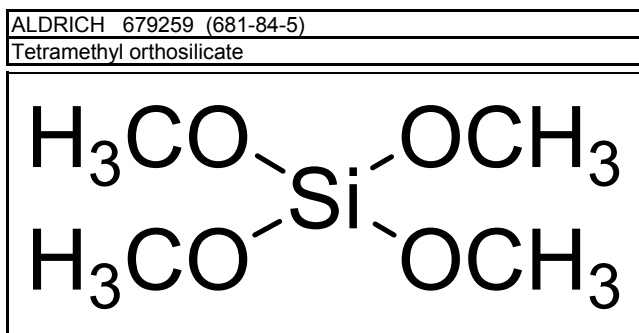
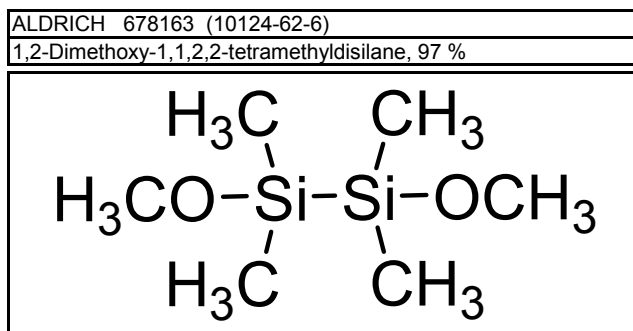
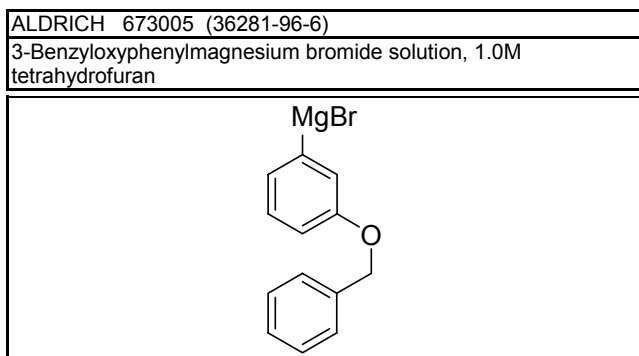
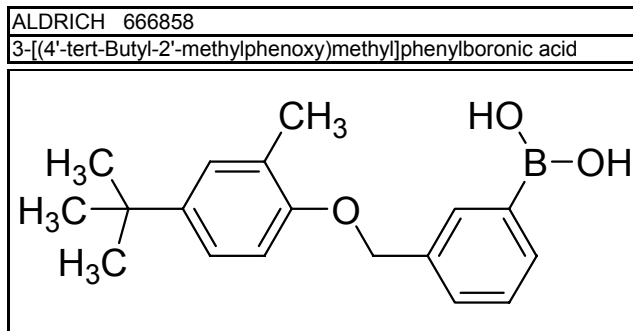
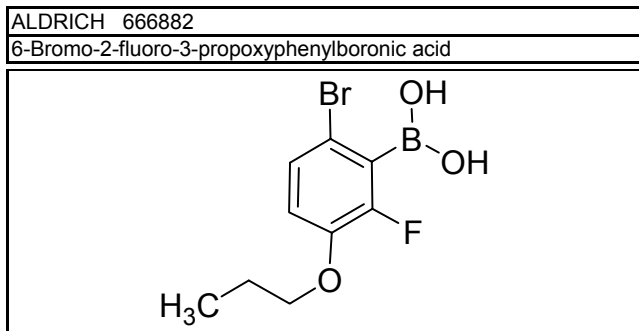
4-(4'-Isopentyloxyphenyl)phenylboronic acid



ALDRICH 652121

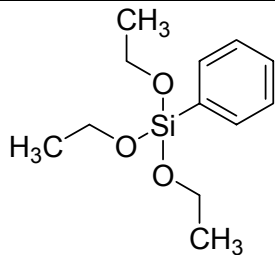
3,5-Dimethyl-4-(3',5'-dimethoxybenzyloxy)phenylboronic acid





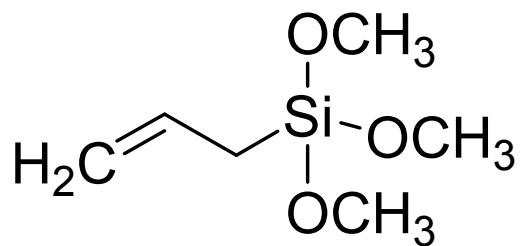
ALDRICH 679291 (780-69-8)

Triethoxyphenylsilane, 98+ %



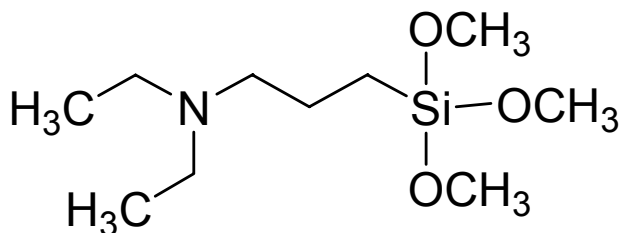
ALDRICH 679267 (2551-83-9)

Allyltrimethoxysilane



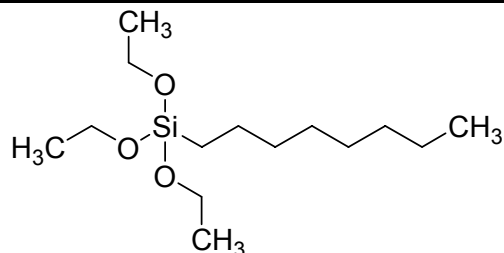
ALDRICH 679356 (41051-80-3)

[3-(Diethylamino)propyl]trimethoxysilane



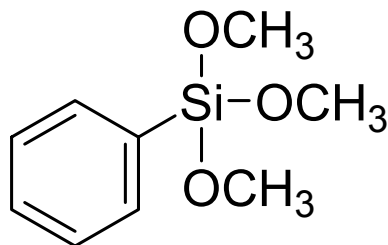
ALDRICH 679305 (2943-75-1)

Triethoxy(octyl)silane



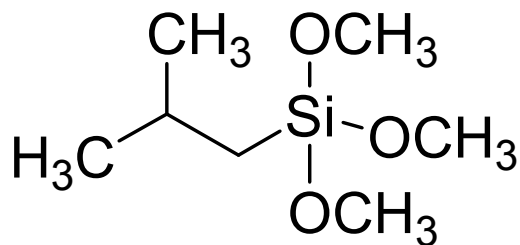
ALDRICH 679313 (2996-92-1)

Trimethoxyphenylsilane



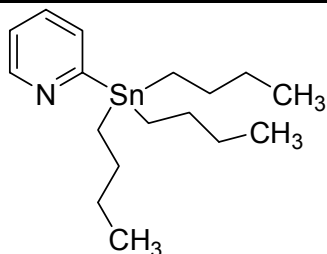
ALDRICH 679364 (18395-30-7)

Isobutyl(trimethoxy)silane



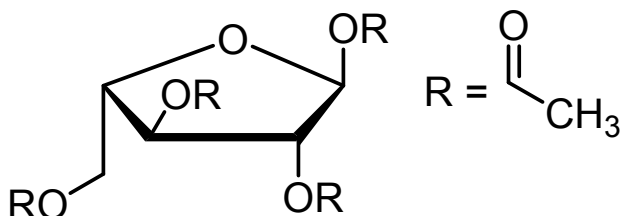
ALDRICH 678333 (17997-47-6)

2-(Tributylstannyl)pyridine, 90 %



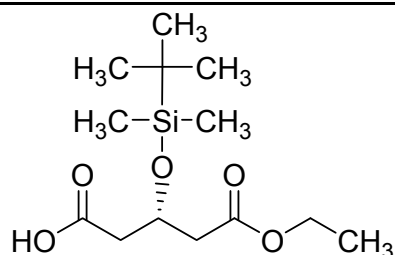
ALDRICH 637440 (79120-81-3)

1,2,3,5-Tetra-O-acetyl-alpha-L-arabinofuranose



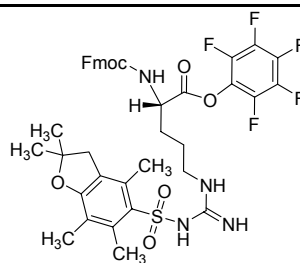
FLUKA 42015

Ethyl (S)-3-(tert-butyl(dimethylsiloxy)glutarate, purum, >= 95.0 % GC

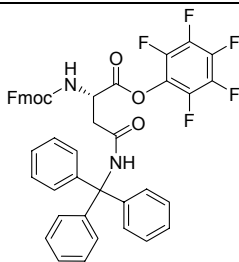


FLUKA 89048

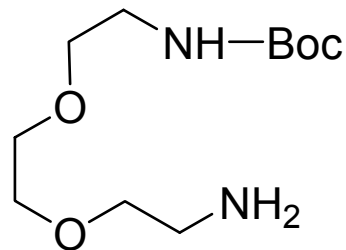
Fmoc-Arg(Pbf)-OPfp, technical, >= 80 % HPLC



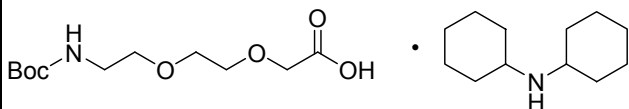
FLUKA 67867 (132388-64-8)
Fmoc-Asn(Trt)-OPfp, purum, >= 98.0 % HPLC



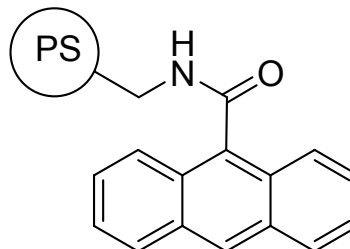
FLUKA 89761 (153086-78-3)
N-Boc-2,2'-(ethylenedioxy)diethylamine, purum, >= 95.0 % NT



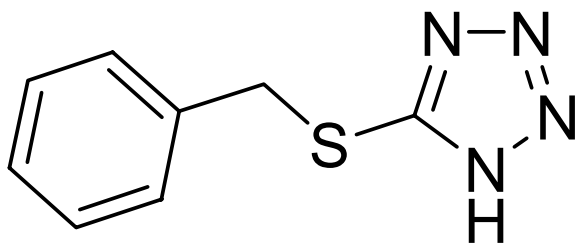
FLUKA 14766
2-[2-(Boc-amino)ethoxy]ethoxyacetic acid dicyclohexylamine salt, purum, >= 98.0 % TLC



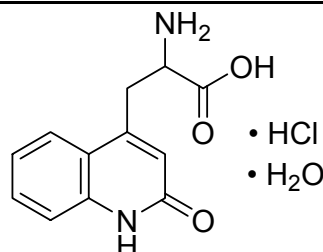
ALDRICH 677868
9-Anthracenecarboxamide, polymer bound



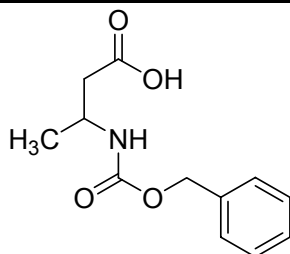
FLUKA 43323 (21871-47-6)
5-(Benzylthio)-1H-tetrazole solution, for DNA synthesis, 0.25M acetonitrile



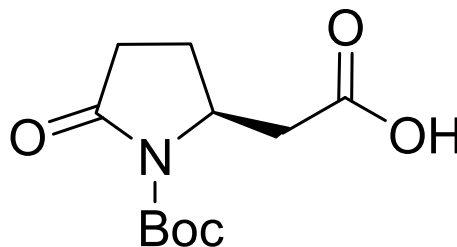
ALDRICH 677337
3-(2-Oxo-1,2-dihydro-4-quinoliny)alanine hydrochloride monohydrate, 96 %



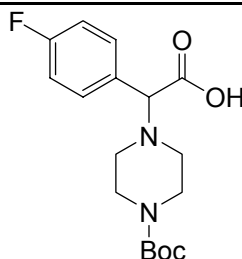
FLUKA 39599
Z-DL-beta-Homoalanine, purum, >= 98.0 % HPLC



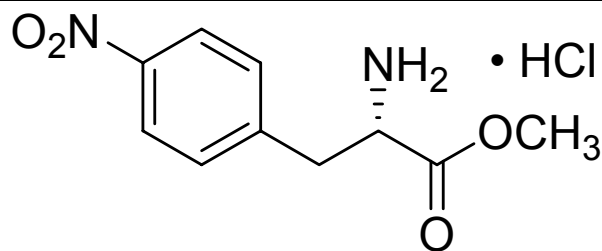
ALDRICH 670618
Boc-beta-Homopyr-OH, purum, >= 99.0 % HPLC



FLUKA 44264
2-(4-Boc-piperazino)-2-(4-fluorophenyl)acetic acid, purum, >= 97.0 % HPLC

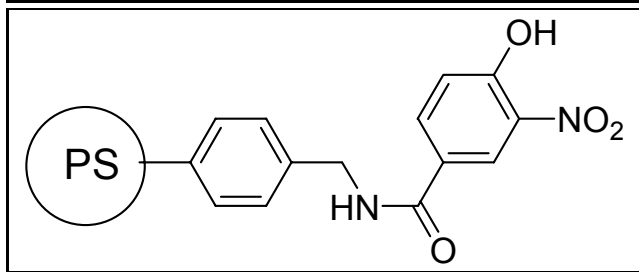


ALDRICH 658421 (17193-40-7)
(S)-(+)-4-Nitrophenylalanine methyl ester hydrochloride, 97 %



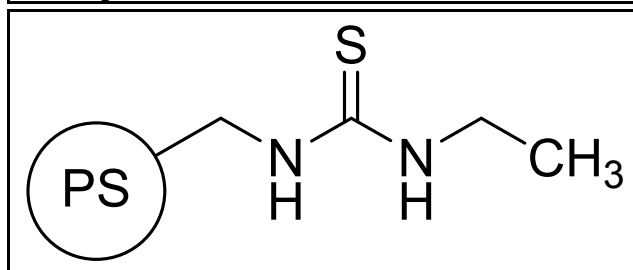
FLUKA 50598

2-Nitrophenol polymer bound, loading ~ 1.3mmol/g



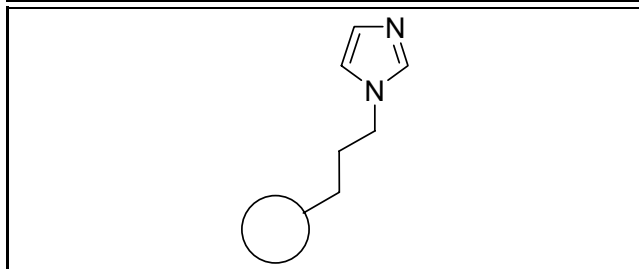
ALDRICH 663360

N-Ethylthiourea, polymer-bound, 70-90mesh, loading 1.5-2.0mmol/g S, 1 %



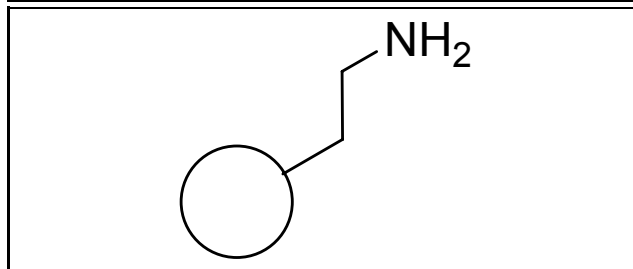
ALDRICH 668605

QuadraPure(R) BDZ, 400-750mum, loading 1.0-1.4mmol/g, macroporous



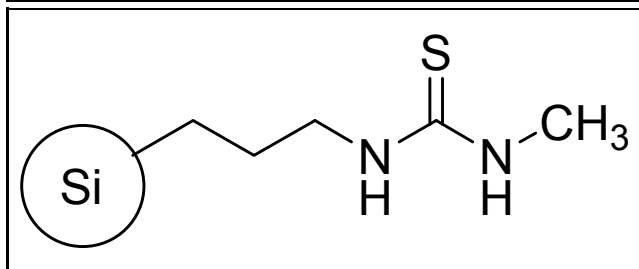
ALDRICH 668583

QuadraPure(R) EDA, 500-800mum, loading 1.5-2.0mmol/g, macroporous



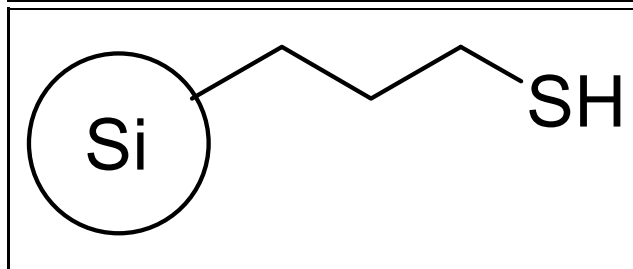
ALDRICH 679518

QuadraSil(R) MTU



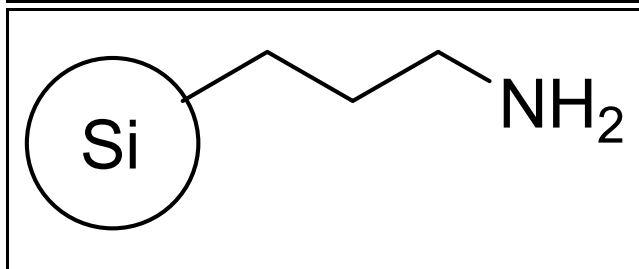
ALDRICH 679526

QuadraSil(R) MP



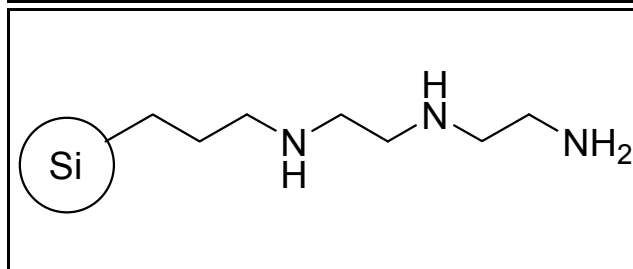
ALDRICH 679534

QuadraSil(R) AP



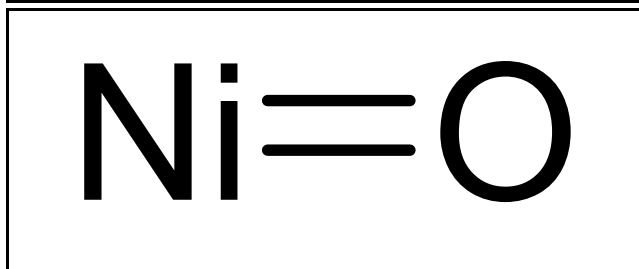
ALDRICH 679496

QuadraSil(R) TA



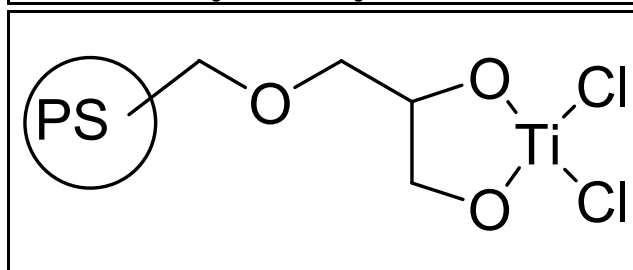
ALDRICH 675172

Nickel oxide supported on silica, 60wt. % NiO

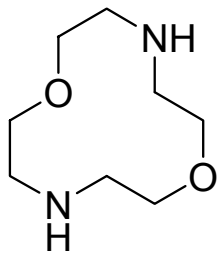


ALDRICH 668214

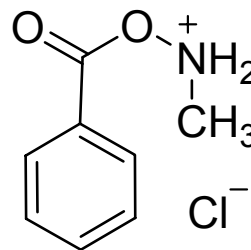
3-Hydroxy-1,2-propanedioxytitanium dichloride, polymer-bound, 100-200mesh, loading 0.5-1.5mmol/g Ti



FLUKA 73753 (294-92-8)
1,7-Diaza-12-crown-4, purum, >= 97.0 % GC



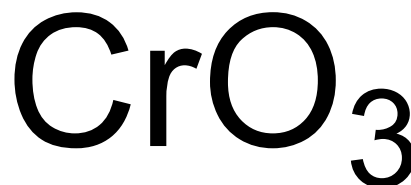
ALDRICH 669393
N-Methyl-O-benzoylhydroxylamine hydrochloride, 97 %



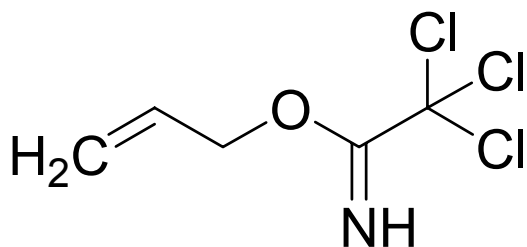
ALDRICH 678732 (10294-34-5)
Boron trichloride solution, 1.0M toluene



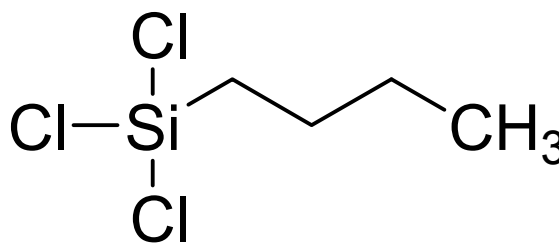
ALDRICH 675644 (1333-82-0)
Chromium(VI) oxide, 99.99 %



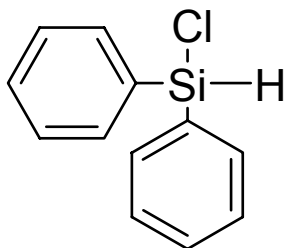
ALDRICH 678414 (51479-73-3)
O-Allyl 2,2,2-trichloroacetimidate, 96 %



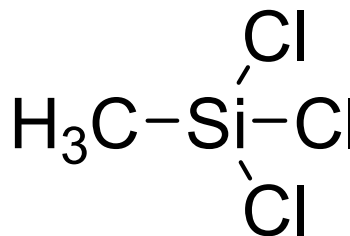
ALDRICH 679224 (7521-80-4)
Butyltrichlorosilane



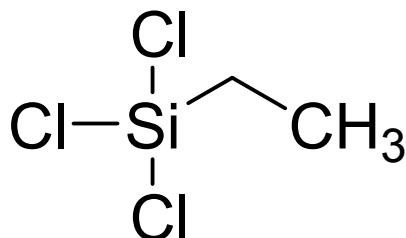
ALDRICH 673935 (1631-83-0)
Chlorodiphenylsilane, technical grade, 90 %



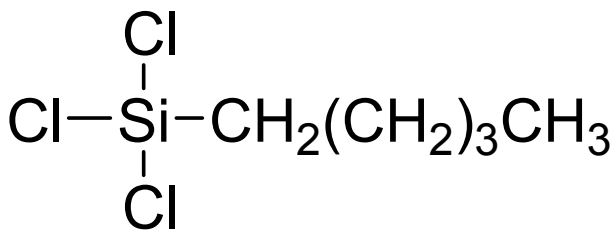
ALDRICH 679208 (75-79-6)
Methyltrichlorosilane, deposition grade, >= 98 % GC, 99.99+ % as metals



ALDRICH 679216 (115-21-9)
Ethyltrichlorosilane, deposition grade, >= 98 % GC, 99.99+ % as metals

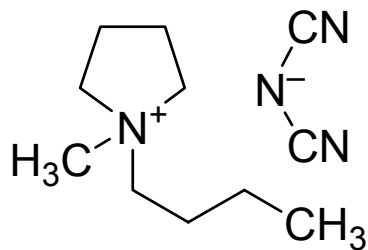


ALDRICH 679194 (107-72-2)
Pentyltrichlorosilane



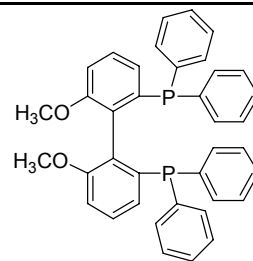
FLUKA 50893 (370865-80-8)

1-Butyl-1-methylpyrrolidinium dicyanamide, $\geq 97.0\%$ T, purum



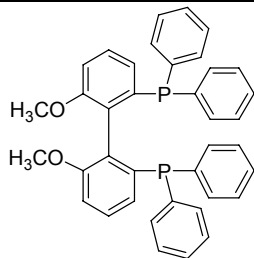
ALDRICH 29510 (133545-16-1)

(R)-(+)-(6,6'-Dimethoxybiphenyl-2,2'-diyl)bis(diphenylphosphine), purum, $\geq 97.0\%$ CHN



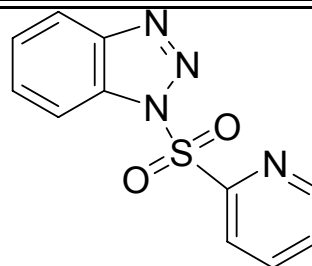
ALDRICH 29511 (133545-17-2)

(S)-(-)-(6,6'-Dimethoxybiphenyl-2,2'-diyl)bis(diphenylphosphine), purum, $\geq 97.0\%$



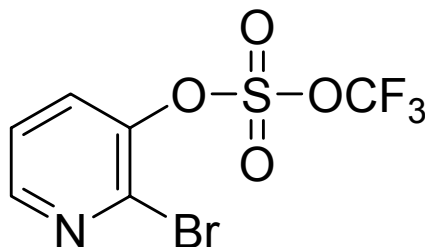
ALDRICH 647349 (678173-42-7)

1-(2-Pyridinylsulfonyl)-1H-benzotriazole, 97 %



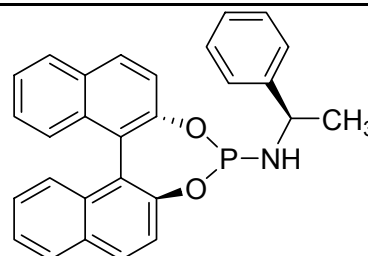
ALDRICH 649945 (157373-97-2)

2-Bromo-3-pyridyl trifluoromethanesulfonate, 96 %



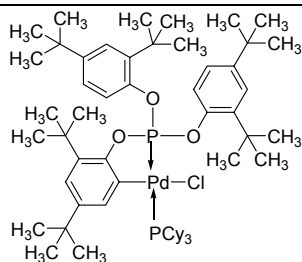
ALDRICH 665347 (422509-53-3)

(S,R)-(+)-(3,5-Dioxa-4-phospha-cyclohepta[2,1-a;4-a']dinaphthalen-4-yl)-(1-phenylethyl)amine, 96 %



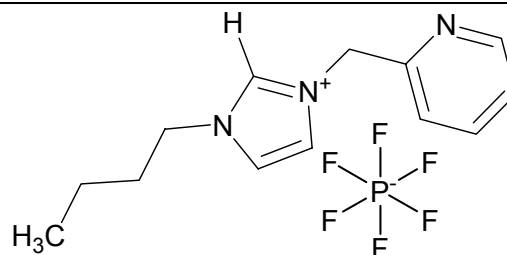
ALDRICH 667315

Chloro(eta2-P,C-tris(2,4-di-tert-butylphenyl)phosphite)(tricyclohexylphosphine)palladium(II)



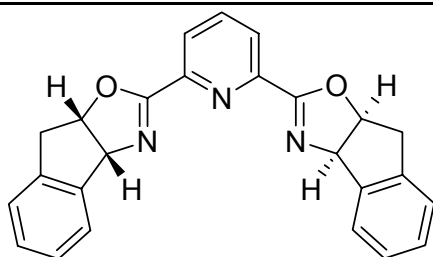
ALDRICH 669121 (873788-09-1)

1-Butyl-3-(2-pyridinylmethyl)-1H-imidazolium hexafluorophosphate, 95 %



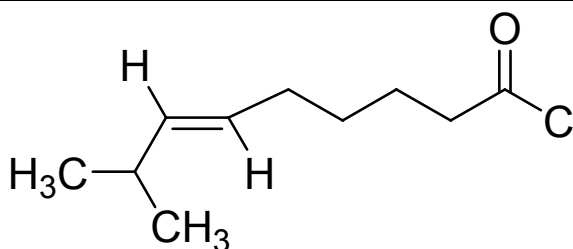
ALDRICH 673986 (357209-32-6)

2,6-Bis[(3aR,8aS)-(+)-8H-indeno[1,2-d]oxazolin-2-yl]pyridine

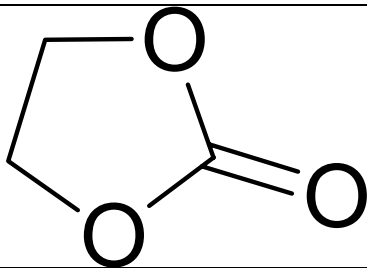


ALDRICH 675083 (95636-02-5)

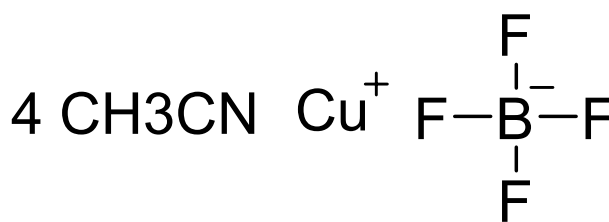
8-Methyl-6-nonenyl chloride, 97 %, predominantly trans



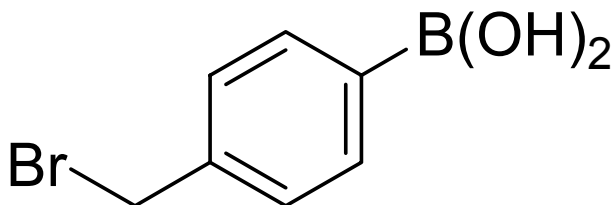
ALDRICH 676802 (96-49-1)
Ethylene carbonate, anhydrous, 99 %



ALDRICH 677892
Tetrakis(acetonitrile)copper(I) tetrafluoroborate, 97 %



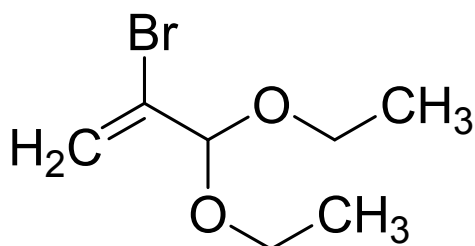
ALDRICH 679437 (68162-47-0)
4-(Bromomethyl)phenylboronic acid



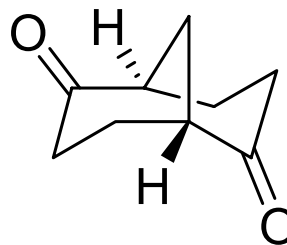
ALDRICH 679488
Rhodium entrapped in aluminum hydroxide matrix preparation, nanoparticles, loading 5wt. %

Rh

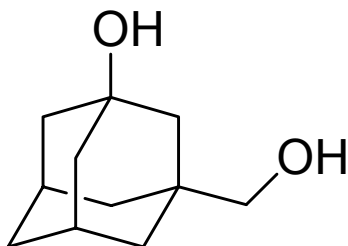
ALDRICH 679593 (17592-40-4)
2-Bromopropenal diethyl acetal, 90 %



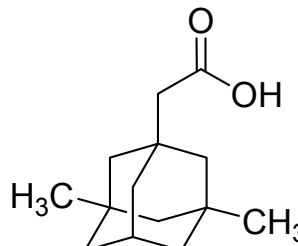
ALDRICH 679933 (16473-11-3)
Bicyclo[3.3.1]nonane-2,6-dione, 97 %



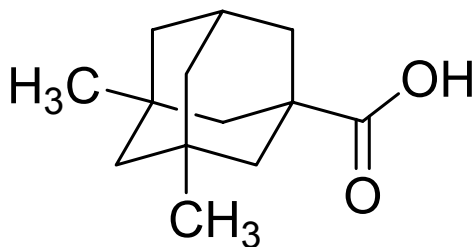
ALDRICH 679968 (38584-37-1)
3-(Hydroxymethyl)-1-adamantol, 97 %



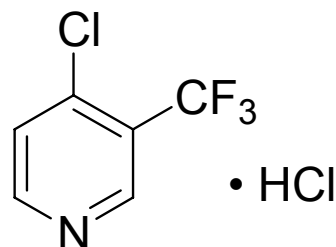
ALDRICH 679976 (14202-14-3)
3,5-Dimethyladamantane-1-acetic acid, 97 %



ALDRICH 679984 (14670-94-1)
3,5-Dimethyladamantane-1-carboxylic acid, 97 %

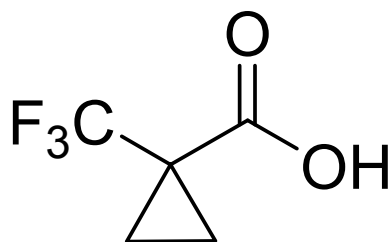


ALDRICH 680036 (732306-24-0)
4-Chloro-3-(trifluoromethyl)pyridine hydrochloride, 97 %



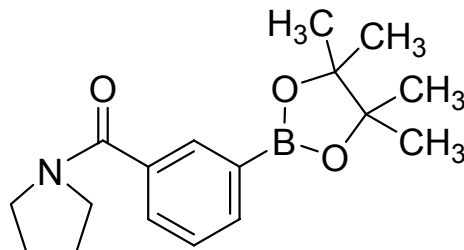
ALDRICH 680141 (277756-46-4)

1-(Trifluoromethyl)cyclopropane-1-carboxylic acid, 97 %



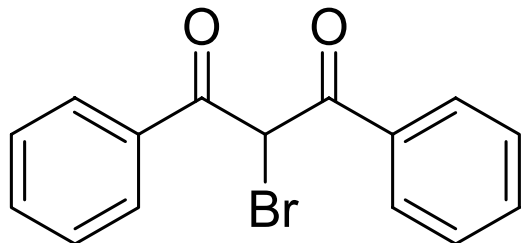
ALDRICH 680222

3-(Pyrrolidine-1-carbonyl)phenylboronic acid pinacol ester, 96 %



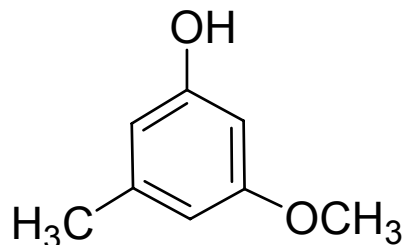
ALDRICH 680443 (728-84-7)

2-Bromo-1,3-diphenylpropane-1,3-dione, 97 %



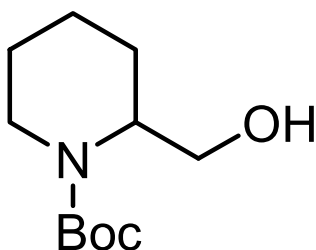
ALDRICH 680893 (3209-13-0)

3-Methoxy-5-methylphenol, 97 %



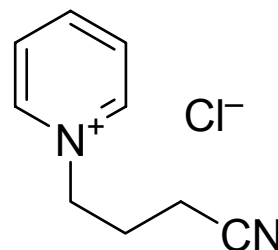
ALDRICH 681296 (157634-00-9)

N-Boc-piperidine-2-methanol, 97 %



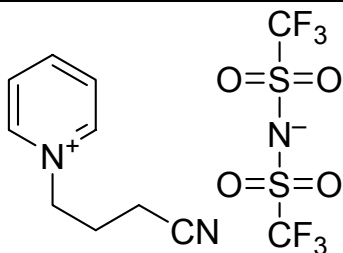
FLUKA 12136

1-(3-Cyanopropyl)pyridinium chloride, purum, >= 98.5 % T



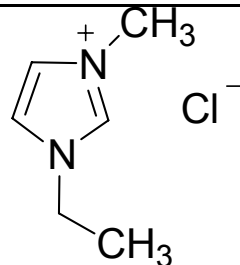
FLUKA 16819

1-(3-Cyanopropyl)pyridinium bis(trifluoromethylsulfonyl)imide, purum, >= 98.5 % T



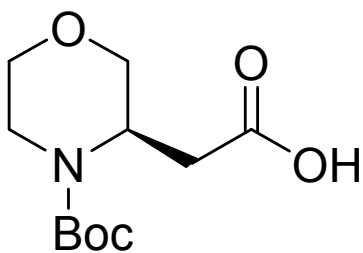
FLUKA 53096 (65039-09-0)

1-Ethyl-3-methylimidazolium chloride, purum, dry, >= 98.0 % T



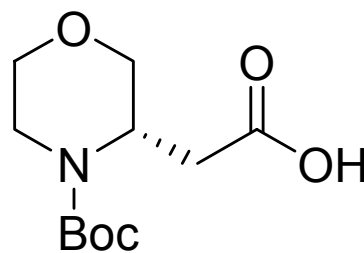
ALDRICH 669830 (761460-03-1)

(R)-4-Boc-3-morpholineacetic acid, 98 %



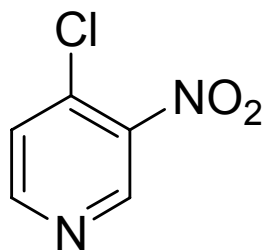
ALDRICH 669954 (839710-38-2)

(S)-4-Boc-3-morpholineacetic acid, 98 %



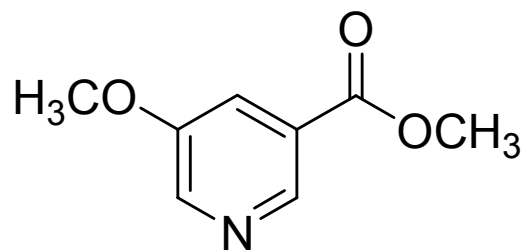
ALDRICH 674079 (13091-23-1)

4-Chloro-3-nitropyridine, 90 %



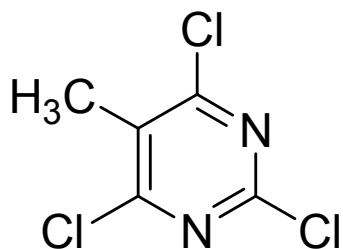
ALDRICH 677884 (29681-46-7)

Methyl 5-methoxy-3-pyridinecarboxylate, 90 %



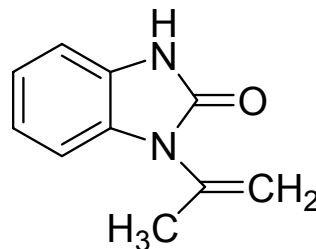
ALDRICH 679054 (1780-36-5)

2,4,6-Trichloro-5-methylpyrimidine, 97 %



ALDRICH 679402 (52099-72-6)

1-Isopropenyl-2-benzimidazolidinone, 97 %



FLUKA 28964 (91599-79-0)

(S)-1-Benzyl-3-hydroxypiperidine, purum, ≥ 97.0 % GC

