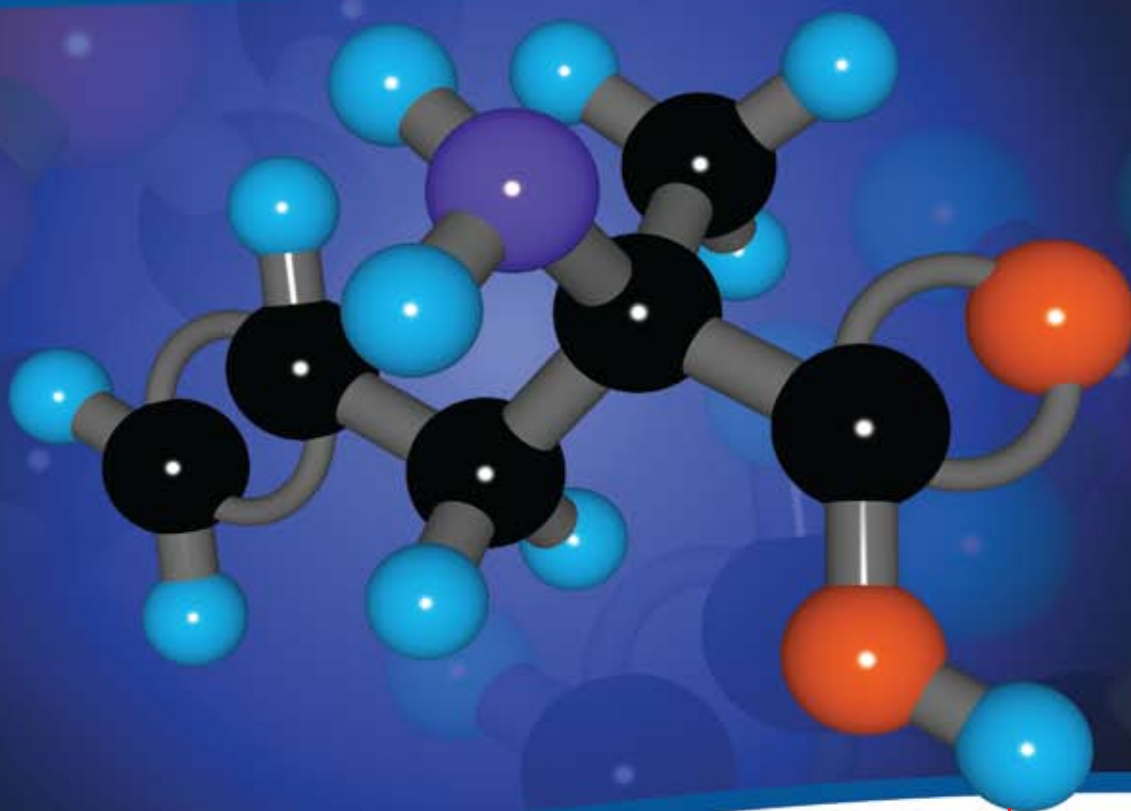


Unnatural Amino Acids



(R)-Allylalanine: a versatile α,α -disubstituted amino acid for peptide synthesis.

Unnatural Amino Acids

Natural Amino Acid
Building Blocks for
Peptide Synthesis

New Tools for PEGylation

Introduction

Natural as well as unnatural amino acids are utilized as chiral building blocks, conformational restraints, molecular scaffolds, and active pharmaceutical ingredients. They represent a nearly infinite array of diverse structural elements for the development of new therapeutic drugs or functional materials. With an increased number of therapeutic targets and novel improved drug delivery methodologies the market for peptide drugs is growing considerably faster than overall pharmaceuticals. Currently hundreds of these therapeutic peptides and peptidomimetics are being tested in clinical phases and a steadily growing number are finally reaching the market.

Sigma-Aldrich® is proud to offer you a choice of more than 2400 amino acid derivatives to construct and tailor your peptides. Furthermore, you will be able to obtain all of the necessary tools for solution or solid phase peptide synthesis from a convenient, single source.

This issue of ChemFiles can only display a small fraction of the broad Sigma-Aldrich portfolio and will focus on recent additions to our unnatural amino acids offering, a convenient listing of most needed, essential amino acid building blocks, and innovative new tools for PEGylation. For a complete list of available building blocks and reagents, please visit sigma-aldrich.com/chemprod. If you are unable to find the specific building block or reagent for your needs, "Please Bother Us" with your suggestions at matthias.junkers@sial.com, or contact your local Sigma-Aldrich office (see back cover).



Matthias Junkers
Product Manager

Table of Contents

Introduction.....	2
Unnatural Amino Acids	3
Glycine & Phenylglycine Derivatives.....	3
Alanine Derivatives	4
Phenylalanine Derivatives.....	4
Miscellaneous Amino Acids	5
Nagase α,α -disubstituted Amino Acids	7
Proline Derivatives.....	8
β -Amino Acids and Homologs.....	11
Natural Amino Acid Building Blocks for Peptide Synthesis	22
BOC-protected Amino Acids.....	22
FMOC-protected Amino Acids	24
New Tools for PEGylation.....	26
New PEG derivatives	26

About Our Cover

The cover structure depicts the α,α -disubstituted amino acid (*R*)-allylalanine (Aldrich **690902**) as a representative example of the innovative new additions to Sigma-Aldrich's portfolio of amino acid building blocks. The olefin moiety in the side chain of this unnatural amino acid can be used to synthesize a plethora of different products. Incorporated into a peptide the olefin moiety offers a unique handle with an orthogonal functionality.

ChemFiles

Vol. 8 No. 7

Aldrich Chemical Co., Inc.
Sigma-Aldrich Corporation
6000 N. Teutonia Ave.
Milwaukee, WI 53209, USA

To Place Orders

Telephone	800-325-3010 (USA)
FAX	800-325-5052 (USA)

Customer & Technical Services

Customer Inquiries	800-325-3010
Technical Service	800-231-8327
SAFC®	800-244-1173
Custom Synthesis	800-244-1173
Flavors & Fragrances	800-227-4563
International	414-438-3850
24-Hour Emergency	414-438-3850
Web Site	sigma-aldrich.com
Email	aldrich@sial.com

Subscriptions

To request your **FREE** subscription to *ChemFiles*, please contact us by:

Phone: 800-325-3010 (USA)

Mail: **Attn: Marketing Communications**
Aldrich Chemical Co., Inc.
Sigma-Aldrich Corporation
P.O. Box 355
Milwaukee, WI 53201-9358

Email: sams-usa@sial.com

International customers, please contact your local Sigma-Aldrich office. For worldwide contact information, please see back cover.

ChemFiles are also available in PDF format on the Internet at sigma-aldrich.com/chemfiles.

Aldrich brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product for its particular use. See reverse side of invoice or packing slip for additional terms and conditions of sale.

All prices are subject to change without notice.

ChemFiles (ISSN 1933-9658) is a publication of Aldrich Chemical Co., Inc. Aldrich is a member of the Sigma-Aldrich Group. © 2008 Sigma-Aldrich Co.

Unnatural Amino Acids

Unnatural amino acids, the non-proteinogenic amino acids that either occur naturally or are chemically synthesized, are becoming more and more important as tools for modern drug discovery research. Due to their unlimited structural diversity and functional versatility, they are widely used as chiral building blocks and molecular scaffolds in constructing combinatorial libraries.¹ They represent a powerful tool in drug discovery when incorporated into therapeutic peptidomimetics and peptide analogs. Used as molecular probes they can help to better understand the function of biological systems.²

Peptidomimetics have found widespread recognition as surrogates for therapeutic peptides derived from proteinaceous amino acids because they may show improved *in vivo* stability, enhanced potency, better oral absorption, improved tissue distribution, and increased selectivity of biological response. In this context, unnatural amino acids can be used to alter the tertiary structure of a peptide. Side chain modifications allow the design of drug candidates with a perfect match to their biological target.

Sigma-Aldrich® is dedicated to becoming a strong partner for your Drug Discovery research by offering a broad and unique range of unnatural amino acid building blocks. In this section you will find recent additions to our amino acid portfolio including derivatives of glycine, phenylglycine, alanine, phenylalanine, and innovative α,α -disubstituted amino acids, as well as a comprehensive listing of proline derivatives and β -amino acids.

References: (1) Ma, J.S. *Chem. Today* **2003**, 65. (2) Dougherty, D.A. *Curr Opin Chem Biol.* **2000**, 6, 645.

Glycine & Phenylglycine Derivatives

Boc-(2-indanyl)-Gly-OH, $\geq 98.0\%$ (HPLC) NEW

Boc-(2-indanyl)-L-glycine

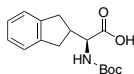
[181227-47-4]

$C_{16}H_{21}NO_4$

FW 291.34

669792-500MG

500 mg



Fmoc-(2-indanyl)-Gly-OH, $\geq 98.0\%$ (HPLC) NEW

Fmoc-(2-indanyl)-L-glycine

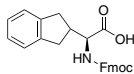
[205526-39-2]

$C_{26}H_{23}NO_4$

FW 413.47

669903-500MG

500 mg



Fmoc-N-(2-Boc-aminoethyl)-Gly-OH, $\geq 97.0\%$ (HPLC)

Fmoc-N-(2-Boc-aminoethyl)glycine

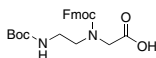
[141743-15-9]

$C_{24}H_{28}N_2O_6$

FW 440.49

670316-500MG

500 mg



Fmoc-N-(4-Boc-aminobutyl)-Gly-OH, $\geq 98.0\%$ (HPLC) NEW

Fmoc-N-(4-Boc-aminobutyl)glycine;

Fmoc-Nlys(Boc)-OH

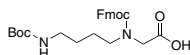
[171856-09-0]

$C_{26}H_{32}N_2O_6$

FW 468.54

670413-500MG

500 mg



Fmoc-N-(2,4-dimethoxybenzyl)-Gly-OH, $\geq 98.0\%$ (HPLC) NEW

Fmoc-N-(2,4-dimethoxybenzyl)glycine;

Fmoc-Dmb-Gly-OH

[166881-42-1]

$C_{26}H_{25}NO_6$

FW 447.48

689459-250MG

250 mg

689459-1G

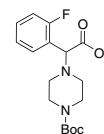
1 g



2-(4-Boc-piperazino)-2-(2-fluorophenyl)acetic acid, $\geq 97.0\%$ (HPLC)

$C_{17}H_{23}FN_2O_4$

FW 338.37



12234-250MG-F

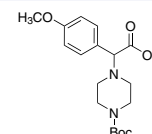
250 mg

12234-1G-F

1 g

2-(4-Boc-piperazino)-2-(4-methoxyphenyl)acetic acid, $\geq 97.0\%$ (HPLC)

[868260-17-7]



16297-250MG-F

250 mg

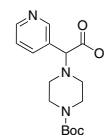
16297-1G-F

1 g

2-(4-Boc-piperazino)-2-(3-pyridyl)acetic acid, $\geq 97.0\%$ (HPLC)

$C_{16}H_{23}N_3O_4$

FW 321.37



16762-250MG-F

250 mg

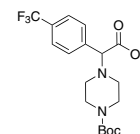
16762-1G-F

1 g

2-(4-Boc-piperazino)-2-[4-(trifluoromethyl)phenyl]acetic acid, $\geq 97.0\%$ (HPLC)

$C_{18}H_{23}F_3N_2O_4$

FW 388.38



28963-250MG-F

250 mg

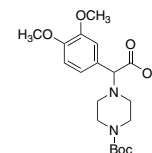
28963-1G-F

1 g

2-(4-Boc-piperazino)-2-(3,4-dimethoxyphenyl)acetic acid, $\geq 97.0\%$ (HPLC)

$C_{19}H_{28}N_2O_6$

FW 380.44

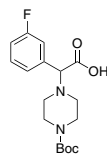
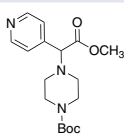
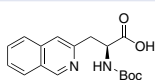


30096-250MG-F

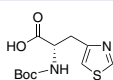
250 mg

30096-1G-F

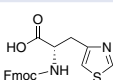
1 g

2-(4-Boc-piperazino)-2-(3-fluorophenyl)acetic acid, ≥97.0% (HPLC)C₁₇H₂₃FN₂O₄
FW 338.3739563-250MG-F 250 mg
39563-1G-F 1 g**Methyl 2-(4-Boc-piperazino)-2-(4-pyridyl)acetate, ≥90% (HPLC)**2-(4-Boc-piperazino)-2-(4-pyridyl)acetic acid methyl ester
C₁₇H₂₅N₃O₄
FW 335.4068892-250MG-F 250 mg
68892-1G-F 1 g**Alanine Derivatives****Boc-β-(2-quinolyl)-Ala-OH, ≥98.0% (HPLC)**Boc-3-(2-quinolyl)-L-alanine;
(S)-2-(Boc-amino)-3-(2-quinolyl)propionic acid
[161453-37-8]
C₁₇H₂₀N₂O₄
FW 316.35

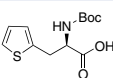
670456-500MG 500 mg

Boc-β-(4-thiazolyl)-Ala-OH, ≥98.0% (HPLC)Boc-3-(4-thiazolyl)-L-alanine;
(S)-2-(Boc-amino)-3-(4-thiazolyl)propionic acid
[119434-75-2]
C₁₁H₁₆N₂O₄S
FW 272.32

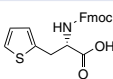
670553-1G 1 g

Fmoc-β-(4-thiazolyl)-Ala-OH, ≥98.0% (HPLC)(S)-2-(Fmoc-amino)-3-(4-thiazolyl)propionic acid;
Fmoc-3-(4-thiazolyl)-L-alanine
[205528-32-1]
C₂₁H₁₈N₂O₄S
FW 394.44

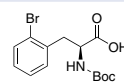
670677-500MG 500 mg

Boc-β-(2-thienyl)-D-Ala-OH, ≥98.0% (HPLC)(R)-2-(Boc-amino)-3-(2-thienyl)propionic acid;
Boc-3-(2-thienyl)-D-alanine
[78452-55-8]
C₁₂H₁₇NO₄S
FW 271.33

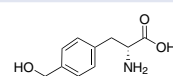
670782-1G 1 g

Fmoc-β-(2-thienyl)-Ala-OH, ≥98.0% (HPLC)(S)-2-(Fmoc-amino)-3-(2-thienyl)propionic acid;
Fmoc-3-(2-thienyl)-L-alanine
[130309-35-2]
C₂₂H₁₉NO₄S
FW 393.46

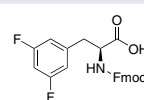
670898-1G 1 g

Phenylalanine Derivatives**Boc-Phe(2-Br)-OH, ≥98.0% (HPLC)**Boc-2-bromo-L-phenylalanine
[261165-02-0]
C₁₄H₁₈BrNO₄
FW 344.20

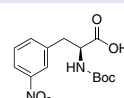
672203-500MG 500 mg

4-(Hydroxymethyl)-D-phenylalanine, ≥97.0% (HPLC)(R)-2-Amino-3-[4-(hydroxymethyl)phenyl]propionic acid
[15720-17-9]
C₁₀H₁₃NO₃
FW 195.22

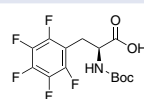
43667-500MG 500 mg

Fmoc-Phe(3,5-F₂)-OH, ≥98.0% (HPLC)Fmoc-3,5-difluoro-L-phenylalanine
[205526-24-5]
C₂₄H₁₉F₂NO₄
FW 423.41

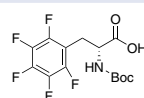
669687-500MG 500 mg

Boc-Phe(3-NO₂)-OH, ≥98.0% (HPLC)Boc-3-nitro-L-phenylalanine
[131980-29-5]
C₁₄H₁₈N₂O₆
FW 310.30

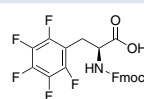
670014-1G 1 g

Boc-pentafluoro-L-phenylalanine, ≥98.0% (HPLC)(S)-2-(Boc-amino)-3-(pentafluorophenyl)propionic acid
[34702-60-8]
C₁₄H₁₄F₅NO₄
FW 355.26

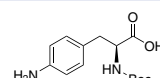
670138-1G 1 g

Boc-pentafluoro-D-phenylalanine, ≥98.0% (HPLC)(R)-2-(Boc-amino)-3-(pentafluorophenyl)propionic acid
[136207-26-6]
C₁₄H₁₄F₅NO₄
FW 355.26

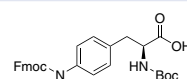
670235-1G 1 g

Fmoc-pentafluoro-L-phenylalanine, ≥98.0% (HPLC)(S)-2-(Fmoc-amino)-3-(pentafluorophenyl)propionic acid
[205526-32-5]
C₂₄H₁₆F₅NO₄
FW 477.38

670340-500MG 500 mg

Boc-Phe(4-NH₂)-OH, ≥98.0% (HPLC)Boc-4-amino-L-phenylalanine
[55533-24-9]
C₁₄H₂₀N₂O₄
FW 280.32

671770-500MG 500 mg

Boc-4-(Fmoc-amino)-L-phenylalanine, ≥98.0% (HPLC)[114346-31-5]
C₂₅H₃₀N₂O₆
FW 502.56

671886-500MG 500 mg

Fmoc-Phe(2-Br)-OH, ≥98.0% (HPLC)

Fmoc-2-bromo-L-phenylalanine

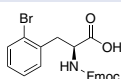
[220497-47-2]

C₂₄H₂₀BrNO₄

FW 466.32

672319-500MG

500 mg

**Boc-Phe(4-Br)-OH, ≥98.0% (HPLC)**

Boc-4-bromo-L-phenylalanine

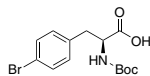
[62129-39-9]

C₁₄H₁₈BrNO₄

FW 344.20

672424-1G

1 g

**Boc-D-Phe(4-Br)-OH, ≥98.0% (HPLC)**

Boc-4-bromo-D-phenylalanine

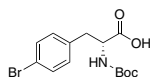
[79561-82-3]

C₁₄H₁₈BrNO₄

FW 344.20

672521-1G

1 g

**Fmoc-Phe(4-Br)-OH, ≥98.0% (HPLC)**

Fmoc-4-bromo-L-phenylalanine

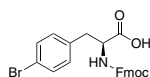
[198561-04-5]

C₂₄H₂₀BrNO₄

FW 466.32

672637-1G

1 g

**Boc-D-Phe(3-Cl)-OH, ≥98.0% (HPLC)**

Boc-3-chloro-D-phenylalanine

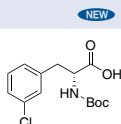
[80102-25-6]

C₁₄H₁₈ClNO₄

FW 299.75

672750-1G

1 g

**Boc-Phe(3,5-F₂)-OH, ≥98.0% (HPLC)**

Boc-3,5-difluoro-L-phenylalanine

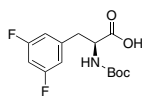
[205445-52-9]

C₁₄H₁₇F₂NO₄

FW 301.29

669563-500MG

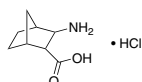
500 mg

**Miscellaneous Amino Acids****(*cis*)-3-Aminobicyclo[2.2.1]heptane-2-carboxylic acid hydrochloride, ≥98.0% (TLC)**(*cis*)-2-Amino-3-carboxybicyclo[2.2.1]heptane hydrochlorideC₈H₁₃NO₂ · HCl

FW 191.66

08356-1G-F

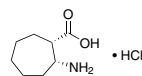
1 g

***cis*-2-Aminocycloheptanecarboxylic acid hydrochloride, ≥98.0% (TLC)**C₈H₁₅NO₂ · HCl

FW 193.67

11252-500MG-F

500 mg

**Boc-4,5-dehydro-Leu-OH (dicyclohexylammonium) salt, ≥96.0% (HPLC)**

Boc-4,5-dehydro-L-leucine

(dicyclohexylammonium) salt;

(S)-2-(Boc-amino)-4-methyl-4-pentenoic acid (dicyclohexylammonium) salt

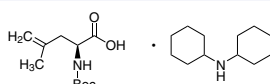
[87720-54-5]

C₁₁H₁₉NO₄ · C₁₂H₂₃N

FW 410.59

11578-50MG

50 mg

***cis*-2-Amino-2-methylcyclohexanecarboxylic acid hydrochloride, ≥98.0% (CHN)**

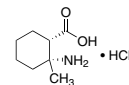
[202921-88-8]

C₈H₁₅NO₂ · HCl

FW 193.67

30254-500MG-F

500 mg

***cis*-2-Amino-2-methylcyclopentanecarboxylic acid hydrochloride, ≥98.0% (CHN)**

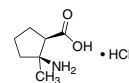
[156292-34-1]

C₇H₁₃NO₂ · HCl

FW 179.64

39927-500MG-F

500 mg

**2-(4-Boc-piperazino)-2-(4-fluorophenyl)acetic acid, ≥97.0% (HPLC)**C₁₇H₂₃FN₂O₄

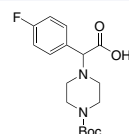
FW 338.37

44264-250MG-F

250 mg

44264-1G-F

1 g

***cis*-6-Amino-3-cyclohexene-1-carboxylic acid hydrochloride, ≥97.0% (AT)***cis*-2-Amino-4-cyclohexene-1-carboxylic acid hydrochloride

[57266-56-5]

C₇H₁₁NO₂ · HCl

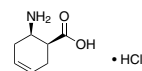
FW 177.63

54105-250MG

250 mg

54105-1G

1 g

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

(R)-1-Boc-morpholin-3-ylacetic acid;

(R)-N-Boc-3-carboxymethyl-morpholine

[761460-03-1]

C₁₁H₁₉NO₅

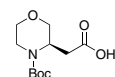
FW 245.27

669830-250MG

250 mg

669830-1G

1 g

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

(S)-1-Boc-morpholin-3-ylacetic acid;

(S)-N-Boc-3-carboxymethyl-morpholine

[839710-38-2]

C₁₁H₁₉NO₅

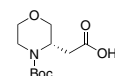
FW 245.27

669954-250MG

250 mg

669954-1G

1 g

**(*S*)-N-Boc-1,2,3,6-tetrahydro-2-pyridinecarboxylic acid, ≥98.0% (HPLC)**

(S)-N-Boc-1,2,3,6-Tetrahydro-2-picolinic acid

[417726-36-4]

C₁₁H₁₇NO₄

FW 227.26

670286-250MG

250 mg

670286-1G

1 g

**Boc-beta-Homopyr-OH, ≥99.0% (HPLC)**

(S)-(1-Boc-5-oxo-pyrrolidin-2-yl)acetic acid

[948594-96-5]

C₁₁H₁₇NO₅

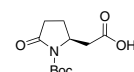
FW 243.26

670618-250MG

250 mg

670618-1G

1 g



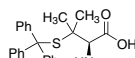
Fmoc-Pen(Trt)-OH, 96%

Fmoc-S-trityl-L-penicillamine;
Fmoc-β,β-dimethyl-Cys(Trt)-OH
[201531-88-6]

C₃₉H₃₅NO₄S
FW 613.76
≥95.5% (HPLC)

670642-500MG

500 mg

**(S)-2-(Boc-amino)octanedioic acid, 97%**

Boc-L-2-aminosuberic acid
[66713-87-9]

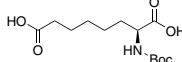
C₁₃H₂₃NO₆
FW 289.32
≥96.5% (GCF)

671983-250MG

250 mg

671983-1G

1 g

**(R)-2-(Boc-amino)octanedioic acid, 99%**

Boc-D-2-aminosuberic acid
[75113-71-2]

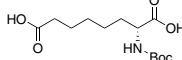
C₁₃H₂₃NO₆
FW 289.32
≥98.5% (GC)

672092-250MG

250 mg

672092-1G

1 g

**exo-cis-3-Aminobicyclo[2.2.1]hept-5-ene-2-carboxylic acid hydrochloride, ≥97.0% (AT)**

[947601-81-2]
C₈H₁₁NO₂ · HCl
FW 189.64

94927-250MG

250 mg

94927-1G

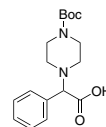
1 g

**2-(4-Boc-piperazino)-2-phenylacetic acid, ≥97.0% (HPLC)**

C₁₇H₂₄N₂O₄
FW 320.38

16298-500MG-F

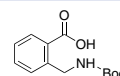
500 mg

**2-(Boc-aminomethyl)benzoic acid, 98%**

Boc-Oamb
[669713-61-5]
C₁₃H₁₇NO₄
FW 251.28
≥97.5% (HPLC)

711659-500MG

500 mg

**Let *Discovery^{CPR}* Put High Throughput Back in Your Operation!****The New Standard in Reagent Management for Medicinal Chemistry and Organic Synthesis****Flexible | Efficient | Convenient**

- Internet-based reagent database and procurement
- Powerful new batch-search and reporting capabilities
- Determine price and availability from your desktop
- From micromoles to grams
- 24–48 hour turnaround time for most compounds
- Reduce waste; eliminate on-site stocking and inventory management
- Specify vial type, atmosphere, labeling/ bar-coding, packaging
- No minimum order required



Custom Packaged Reagents from Sigma-Aldrich puts high throughput back into your chemistry!

When projects demand custom arrays of reagents, *Discovery^{CPR}* can meet the challenge.



To register for an online ordering account or to submit inquiries:

DiscoveryCPR.com

sigma-aldrich.com

SIGMA-ALDRICH®

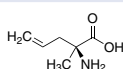
Nagase α,α -disubstituted Amino Acids

Chiral quaternary ammonium catalysts can be particularly useful in the field of asymmetric synthesis. Keji Maruoka and co-workers from Kyoto University have reported potent C_2 -symmetric ammonium salts that catalyze alkylation reactions under remarkably low catalyst loadings.³ Nagase, in close cooperation with Prof. Maruoka has applied his methodology in industrially viable procedures to synthesize α -monosubstituted and novel α,α -disubstituted α -amino acids by alkylation of alanine- (Scheme 1) and glycine-derived (Scheme 2) Schiff bases with alkyl halides. Research quantities of novel Nagase α,α -disubstituted amino acids are now available through Sigma-Aldrich®. Enantiopure allylalanines for example will open pathways to completely new structures. Rutjes and co-workers have already shown that the comparable allylglycine is a remarkable precursor for the synthesis of a broad range of building blocks.⁴

References: (3)(a) Kitamura, M. et al. *Angew. Chem. Int. Ed.* **2005**, *44*, 1549. (b) Ooi, T.; Maruoka, K. *Aldrichimica Acta* **2007**, *40*, 77. (4)(a) Rutjes, F.P.J.T. et al. *J. Org. Biol. Chem.* **2005**, *3*, 3435. (b) Rutjes, F.P.J.T. et al. *J. Chem. Soc., Perkin Trans. 1* **2000**, 4197.

(S)-(-)- α -Allylalanine

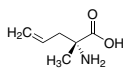
(S)-2-Amino-2-methyl-4-pentenoic acid
[96886-55-4]
 $C_6H_{11}NO_2$
FW 129.16



691011-100MG 100 mg
691011-500MG 500 mg

(R)-(+)- α -Allylalanine

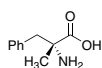
(R)-2-Amino-2-methyl-4-pentenoic acid
[96886-56-5]
 $C_6H_{11}NO_2$
FW 129.16



690902-100MG 100 mg
690902-500MG 500 mg

α -Methyl-D-phenylalanine, $\geq 98.0\%$ (HPLC)

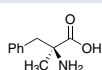
(R)-(+)-2-Amino-2-methyl-3-phenylpropionic acid;
 α -Me-D-Phe-OH
[17350-84-4]
 $C_{10}H_{13}NO_2$
FW 179.22



690791-250MG 250 mg

α -Methyl-L-phenylalanine, $\geq 98.0\%$ (HPLC)

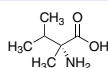
(S)-(-)-2-Amino-2-methyl-3-phenylpropionic acid;
 α -Me-L-Phe-OH
[23239-35-2]
 $C_{10}H_{13}NO_2$
FW 179.22



690694-100MG 100 mg
690694-500MG 500 mg

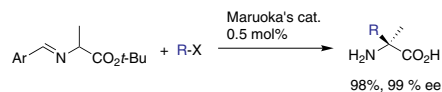
(R)-(+)- α -Methylvaline, 99%

α -Me-D-Val-OH; α -Methyl-D-valine
[53940-82-2]
 $C_6H_{13}NO_2$
FW 131.17

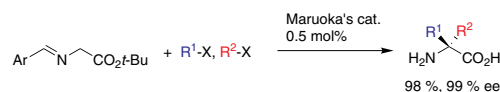


► $\geq 98.5\%$ (TLC)

690589-100MG 100 mg
690589-500MG 500 mg



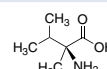
Scheme 1



Scheme 2

(S)-(-)- α -Methylvaline, 99%

α -Methyl-L-valine; α -Me-Val-OH
[53940-83-3]
 $C_6H_{13}NO_2$
FW 131.17

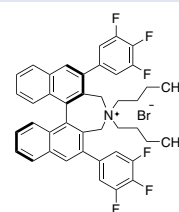


► $\geq 98.5\%$ (TLC)

690465-500MG 500 mg

(11bR)-(-)-4,4-Dibutyl-4,5-dihydro-2,6-bis(3,4,5-trifluorophenyl)-3H-dinaphth[2,1-c:1',2'-e]azepinium bromide

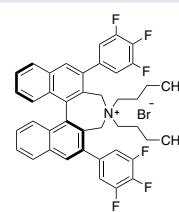
[887938-70-7]
 $C_{42}H_{36}BrF_6N$
FW 748.64



687596-50MG 50 mg

(11bS)-(+)-4,4-Dibutyl-4,5-dihydro-2,6-bis(3,4,5-trifluorophenyl)-3H-dinaphth[2,1-c:1',2'-e]azepinium bromide

[851942-89-7]
 $C_{42}H_{36}BrF_6N$
FW 748.64



677086-50MG 50 mg

Proline Derivatives

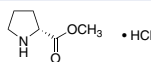
Proline is a non-polar, natural amino acid that forms a tertiary amide when incorporated into peptides. Thus, it is the only proteinogenic amino acid that does not act as a hydrogen bond donor in a peptide chain. Proline is known as a classical breaker of both the α -helical and β -sheet secondary structures in proteins and peptides, and it plays a crucial role in protein folding. Synthetic proline derivatives, mimetics and analogs offer further options to tune the biological, pharmaceutical, or physicochemical properties of peptides and proteins.

In recent years proline derivatives and analogs have also found increasing popularity as organocatalysts in asymmetric synthesis.⁵

References: (5)(a) Vignola, N.; List, B. *J. Am. Chem. Soc.* **2003**, *125*, 450. (b) Dalko, P.L.; Moisan, L.; *Angew. Chem. Int. Ed.* **2004**, *43*, 5138.

D-Proline methyl ester hydrochloride, $\geq 98.0\%$ (AT)

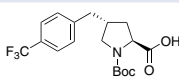
$C_6H_{11}NO_2 \cdot HCl$
FW 165.62



689491-1G 1 g
689491-5G 5 g

Boc-(R)-4-[4-(trifluoromethyl)benzyl]-L-proline, $\geq 98.0\%$ (HPLC)

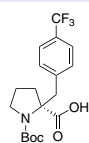
[957311-17-0]
 $C_{18}H_{22}F_3NO_4$
FW 373.37



01336-500MG-F 500 mg

Boc-(S)- α -(4-trifluoromethylbenzyl)-Pro-OH, $\geq 98.0\%$ (HPLC)

Boc- α -(4-trifluoromethylbenzyl)-D-proline; (S)-1-Boc-2-(4-trifluoromethylbenzyl)-2-pyrrolidinecarboxylic acid
 $C_{18}H_{22}F_3NO_4$
FW 373.37



05199-500MG-F 500 mg

Boc-(S)- α -allyl-Pro-OH, $\geq 98.0\%$ (HPLC)

Boc- α -allyl-D-proline;
(S)-2-Allyl-1-Boc-2-pyrrolidinecarboxylic acid
[706806-59-9]
 $C_{13}H_{21}NO_4$
FW 255.31



06486-500MG-F 500 mg

Boc-(R)- α -allyl-Pro-OH, $\geq 98.0\%$ (HPLC)

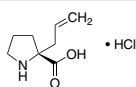
Boc- α -allyl-L-proline;
(R)-2-Allyl-1-Boc-2-pyrrolidinecarboxylic acid
[144085-23-4]
 $C_{13}H_{21}NO_4$
FW 255.31



06538-500MG-F 500 mg

(R)- α -Allyl-proline hydrochloride, $\geq 98.0\%$ (TLC)

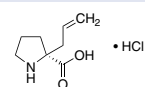
α -Allyl-L-proline hydrochloride;
(R)-2-Allyl-2-pyrrolidinecarboxylic acid hydrochloride
[177206-69-8]
 $C_8H_{13}NO_2 \cdot HCl$
FW 191.66



06541-500MG-F 500 mg

(S)- α -Allyl-proline hydrochloride, $\geq 98.0\%$ (TLC)

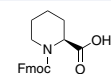
α -Allyl-D-proline hydrochloride;
(S)-2-Allyl-2-pyrrolidinecarboxylic acid hydrochloride
[129704-91-2]
 $C_8H_{13}NO_2 \cdot HCl$
FW 191.66



06594-500MG-F 500 mg

Fmoc-Pip-OH, $\geq 98.0\%$ (HPLC)

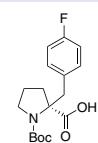
(S)-1-Fmoc-piperidine-2-carboxylic acid;
N-Fmoc-L-pipecolic acid
[86069-86-5]
 $C_{21}H_{21}NO_4$
FW 351.40



09777-1G 1 g

Boc-(S)- α -(4-fluorobenzyl)-Pro-OH, $\geq 98.0\%$ (HPLC)

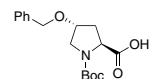
Boc- α -(4-fluorobenzyl)-D-proline;
(S)-1-Boc-2-(4-fluorobenzyl)-2-pyrrolidinecarboxylic acid
[706806-65-7]
 $C_{17}H_{22}FNO_4$
FW 323.36



14931-500MG-F 500 mg

Boc-Hyp(Bzl)-OH, $\geq 98.0\%$ (HPLC)

Boc-O-benzyl-L-hydroxyproline
[54631-81-1]
 $C_{17}H_{23}NO_5$
FW 321.37



15535-5G 5 g

α -Methyl-L-proline, $\geq 98.0\%$ (TLC)

[42856-71-3]
 $C_6H_{11}NO_2$
FW 129.16



17249-250MG 250 mg

17249-1G 1 g

L-Pipecolic acid, 99% (titration)

L-Homoproline; (S)-(-)-2-Piperidinecarboxylic acid
[3105-95-1]
 $C_6H_{11}NO_2$
FW 129.16



P2519-100MG 100 mg

P2519-250MG 250 mg

P2519-1G 1 g

D-Pipecolic acid, 99%

(R)-(+)-2-Piperidinecarboxylic acid;
(R)-Piperidine-2-carboxylic acid; D-Homoproline
[1723-00-8]
 $C_6H_{11}NO_2$
FW 129.16

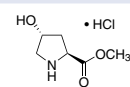


268062-25MG 25 mg

268062-100MG 100 mg

L-4-Hydroxyproline methyl ester hydrochloride, $\geq 98.0\%$ (TLC)

[40216-83-9]
 $C_6H_{11}NO_3 \cdot HCl$
FW 181.62



30681-5G 5 g

Boc- α -(diphenylmethyl)-DL-Pro-OH, $\geq 90\%$ (HPLC)

Boc- α -(diphenylmethyl)-DL-proline; Boc- α -(benzhydryl)-DL-proline; 1-Boc-2-(diphenylmethyl)-2-pyrrolidinecarboxylic acid [351002-64-7]

$C_{23}H_{27}NO_4$
FW 381.46



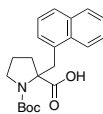
30763-500MG

500 mg

Boc- α -(1-naphthylmethyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -(1-naphthylmethyl)-DL-proline; 1-Boc-2-(1-naphthylmethyl)-2-pyrrolidinecarboxylic acid [351002-65-8]

$C_{23}H_{25}NO_4$
FW 355.43

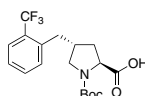


36748-500MG

500 mg

Boc-(R)-4-[2-(trifluoromethyl)benzyl]-L-proline, $\geq 98.0\%$ (HPLC)

$C_{18}H_{22}F_3NO_4$
FW 373.37



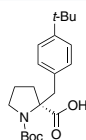
38455-500MG-F

500 mg

Boc-(S)- α -(4-tert-butylbenzyl)-Pro-OH, $\geq 98.0\%$ (HPLC)

Boc- α -(4-tert-butylbenzyl)-D-proline; (S)-1-Boc-2-(4-tert-butylbenzyl)-2-pyrrolidinecarboxylic acid

$C_{21}H_{31}NO_4$
FW 361.48



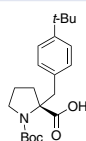
39166-500MG-F

500 mg

Boc-(R)- α -(4-tert-butylbenzyl)-Pro-OH, $\geq 97.0\%$ (HPLC)

(R)-1-Boc-2-(4-tert-butylbenzyl)-2-pyrrolidinecarboxylic acid; Boc- α -(4-tert-butylbenzyl)-L-proline

$C_{21}H_{31}NO_4$
FW 361.48

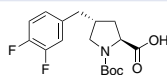


39793-500MG-F

500 mg

Boc-(R)-4-(3,4-difluorobenzyl)-L-proline, $\geq 97.0\%$ (HPLC)

(2S,4R)-1-Boc-4-(3,4-difluorobenzyl)-2-pyrrolidinecarboxylic acid
 $C_{17}H_{21}F_2NO_4$
FW 341.35

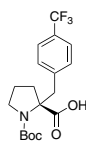


40372-500MG-F

500 mg

Boc-(R)- α -(4-trifluoromethylbenzyl)-Pro-OH, $\geq 98.0\%$ (HPLC)

Boc- α -(4-trifluoromethylbenzyl)-L-proline; (R)-1-Boc-2-(4-trifluoromethylbenzyl)-2-pyrrolidinecarboxylic acid
 $C_{18}H_{22}F_3NO_4$
FW 373.37



42004-500MG-F

500 mg

cis-3-Hydroxy-DL-proline, 95%

(±)-cis-3-Hydroxypyrrolidine-2-carboxylic acid [4298-05-9]

$C_5H_9NO_3$
FW 131.13

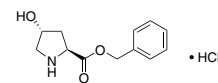


420174-10MG

10 mg

L-4-Hydroxy-proline benzyl ester hydrochloride, $\geq 98.0\%$ (HPLC)

[62147-27-7]
 $C_{12}H_{15}NO_3 \cdot HCl$
FW 257.71

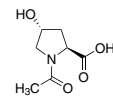


43746-5G

5 g

trans-1-Acetyl-4-hydroxy-L-proline, 99%

N-Acetyl-L-hydroxyproline
[33996-33-7]
 $C_7H_9NO_4$
FW 173.17



441562-500MG

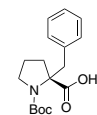
500 mg

441562-10G

10 g

Boc-(R)- α -benzyl-Pro-OH, $\geq 99.0\%$ (HPLC)

Boc- α -benzyl L-proline; (R)-2-Benzyl-1-Boc-2-pyrrolidinecarboxylic acid [706806-60-2]
 $C_{17}H_{23}NO_4$
FW 305.37

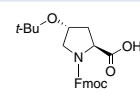


47079-500MG-F

500 mg

Fmoc-Hyp(tBu)-OH, $\geq 98.0\%$ (HPLC)

Fmoc-O-tert-butyl-L-hydroxyproline
[122996-47-8]
 $C_{24}H_{27}NO_5$
FW 409.47

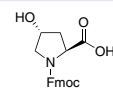


47517-5G-F

5 g

Fmoc-Hyp-OH, $\geq 98.0\%$ (HPLC, sum of enantiomers)

Fmoc-L-4-hydroxyproline
[88050-17-3]
 $C_{20}H_{19}NO_5$
FW 353.37

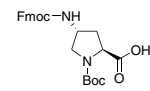


47686-5G

5 g

N-Boc-trans-4-N-Fmoc-amino-L-proline, 97%

(2S,4R)-N-(tert-Butoxycarbonyl)-4-N-(9-fluorenylmethoxycarbonyl)aminopyrrolidine-2-carboxylic acid [176486-63-8]
 $C_{25}H_{28}N_2O_6$
FW 452.50

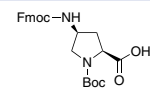


534390-1G

1 g

N-Boc-cis-4-N-Fmoc-amino-L-proline, 97%

(2S,4S)-N-tert-Butoxycarbonyl-4-N-(9-fluorenylmethoxycarbonyl)aminopyrrolidine-2-carboxylic acid [174148-03-9]
 $C_{25}H_{28}N_2O_6$
FW 452.50



534404-1G

1 g

Boc- α -allyl-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -allyl-DL-proline; 2-Allyl-1-Boc-2-pyrrolidinecarboxylic acid [315234-49-2]
 $C_{13}H_{21}NO_4$
FW 255.31

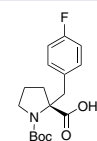


58147-500MG

500 mg

Boc-(R)- α -(4-fluorobenzyl)-Pro-OH, $\geq 98.0\%$ (HPLC)

Boc- α -(4-fluorobenzyl)-L-proline; (R)-1-Boc-2-(4-fluorobenzyl)-2-pyrrolidinecarboxylic acid [706806-64-6]
 $C_{17}H_{22}FNO_4$
FW 323.36



67420-500MG-F

500 mg

Boc- α -Me-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -methyl-DL-proline;
1-Boc-2-methyl-2-pyrrolidinecarboxylic acid
[203869-80-1]
 $C_{11}H_{19}NO_4$
FW 229.27

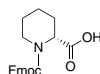


68691-500MG

500 mg

Fmoc-D-Pip-OH, $\geq 98.0\%$ (HPLC)

(*R*)-*N*-Fmoc-piperidine-2-carboxylic acid;
N-Fmoc-D-pipecolic acid
[101555-63-9]
 $C_{21}H_{21}NO_4$
FW 351.40

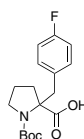


73418-1G

1 g

Boc- α -(4-fluorobenzyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -(4-fluorobenzyl)-DL-proline
[351002-78-3]
 $C_{17}H_{23}FNO_4$
FW 323.36

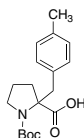


74082-500MG

500 mg

Boc- α -(4-methylbenzyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -(4-methylbenzyl)-DL-proline;
1-Boc-2-(4-methylbenzyl)-2-pyrrolidinecarboxylic acid
[351002-82-9]
 $C_{18}H_{25}NO_4$
FW 319.40



76501-500MG

500 mg

Boc-(*S*)- α -benzyl-Pro-OH, $\geq 97.0\%$ (HPLC)

Boc- α -benzyl-D-proline;
(*S*)-2-Benzyl-1-Boc-2-pyrrolidinecarboxylic acid
[706806-61-3]
 $C_{17}H_{23}NO_4$
FW 305.37

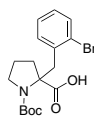


76896-500MG-F

500 mg

Boc- α -(2-bromobenzyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -(2-bromobenzyl)-DL-proline;
1-Boc-2-(2-bromobenzyl)-2-pyrrolidinecarboxylic acid
[351002-85-2]
 $C_{17}H_{22}BrNO_4$
FW 384.26

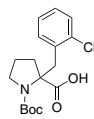


90682-500MG

500 mg

Boc- α -(2-chlorobenzyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -(2-chlorobenzyl)-DL-proline;
1-Boc-2-(2-chlorobenzyl)-2-pyrrolidinecarboxylic acid
[351002-86-3]
 $C_{17}H_{22}ClNO_4$
FW 339.81

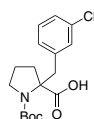


90683-500MG

500 mg

Boc- α -(3-chlorobenzyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

1-Boc-2-(3-chlorobenzyl)-2-pyrrolidinecarboxylic acid;
Boc- α -(3-chlorobenzyl)-DL-proline
[351002-87-4]
 $C_{17}H_{22}ClNO_4$
FW 339.81

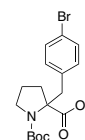


90684-500MG

500 mg

Boc- α -(4-bromobenzyl)-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -(4-bromobenzyl)-DL-proline;
1-Boc-2-(4-bromobenzyl)-2-pyrrolidinecarboxylic acid
[336817-91-5]
 $C_{17}H_{22}BrNO_4$
FW 384.26



94866-500MG

500 mg

Boc- α -propyl-DL-Pro-OH, $\geq 90\%$ (HPLC)

Boc- α -propyl-DL-proline
[351002-88-5]
 $C_{13}H_{23}NO_4$
FW 257.33



95566-500MG

500 mg

Boc- α -benzyl-DL-Pro-OH, $\geq 96.0\%$ (HPLC)

Boc- α -benzyl-DL-proline;
1-Boc-2-benzyl-2-pyrrolidinecarboxylic acid
[351002-72-7]
 $C_{17}H_{23}NO_4$
FW 305.37

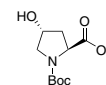


52969-500MG

500 mg

Boc-Hyp-OH, $\geq 98.0\%$ (TLC)

Boc-L-hydroxyproline;
trans-*N*-(*tert*-Butoxycarbonyl)-4-hydroxy-L-proline
[13726-69-7]
 $C_{10}H_{17}NO_5$
FW 231.25



15544-5G

5 g

15544-25G

25 g

***N*-Boc-2-piperidinecarboxylic acid, 98%**

Boc-DL-Pip-OH; ($\pm$)-1-Boc-piperidine-2-carboxylic acid;
N-Boc-DL-pipecolic acid
[98303-20-9]
 $C_{11}H_{19}NO_4$
FW 229.27



495875-10G

10 g

495875-50G

50 g

(*R*)-(+)-*N*-Boc-2-piperidinecarboxylic acid, 98%

Boc-D-Pip-OH; *N*-Boc-D-pipecolic acid; (+)-*N*-Boc-(*R*)-pipecolic acid; (*R*)-*N*-Boc-piperidine-2-carboxylic acid;
(*R*)-(+)-1-(*tert*-Butoxycarbonyl)-2-piperidinecarboxylic acid
[28697-17-8]
 $C_{11}H_{19}NO_4$
FW 229.27

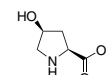


516341-5G

5 g

***cis*-4-Hydroxy-L-proline**

(2*S*,4*S*)-(-)-4-Hydroxy-2-pyrrolidinecarboxylic acid
[618-27-9]
 $C_5H_9NO_3$
FW 131.13



H1637-50MG

50 mg

H1637-250MG

250 mg

H1637-1G

1 g

Boc-Pip-OH, 98%

(*S*)-1-Boc-piperidine-2-carboxylic acid; *N*-Boc-L-pipecolic acid;
(*S*)-(-)-1-(*tert*-Butoxycarbonyl)-2-piperidinecarboxylic acid
[26250-84-0]
 $C_{11}H_{19}NO_4$
FW 229.27



516368-1G

1 g

516368-5G

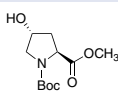
5 g

516368-25G

25 g

***N*-Boc-*trans*-4-hydroxy-L-proline methyl ester, 97%**

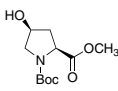
[74844-91-0]
 $C_{11}H_{19}NO_5$
 FW 245.27



639737-1G 1 g
 639737-5G 5 g

***N*-Boc-*cis*-4-hydroxy-L-proline methyl ester, 97%**

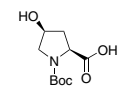
[102195-79-9]
 $C_{11}H_{19}NO_5$
 FW 245.27



654000-1G 1 g
 654000-5G 5 g

***N*-Boc-*cis*-4-Hydroxy-L-proline, 97%**

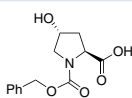
N-Boc-*cis*-4-hydroxypyrrolidine-2-carboxylic acid
 [87691-27-8]
 $C_{10}H_{17}NO_5$
 FW 231.25



654019-1G 1 g
 654019-5G 5 g

Z-Hyp-OH, ≥99.0% (T)

Z-L-4-Hydroxyproline; *N*-Cbz-Hydroxy-L-proline
 [13504-85-3]
 $C_{13}H_{15}NO_5$
 FW 265.26



96310-5G 5 g
 96310-25G 25 g

Pipecolic acid, 98%

Homoproline; 2-Piperidinecarboxylic acid;
 Pipecolic acid free base; (±)-Piperidine-2-carboxylic acid
 [535-75-1]
 $C_6H_{11}NO_2$
 FW 129.16



P45850-25G 25 g
 P45850-100G 100 g

3,4-Dehydro-DL-proline, 98%

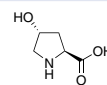
(±)-3-Pyrroline-2-carboxylic acid;
 (±)-2,5-Dihydro-1*H*-pyrrole-2-carboxylic acid
 [3395-35-5]
 $C_5H_7NO_2$
 FW 113.11



862126-100MG 100 mg

***trans*-4-Hydroxy-L-proline, ≥99%**

Hyp; (2*S*,4*R*)-4-Hydroxypyrrolidine-2-carboxylic acid
 [51-35-4]
 $C_5H_9NO_3$
 FW 131.13



H54409-2.5G 2.5 g
 H54409-10G 10 g
 H54409-25G 25 g
 H54409-100G 100 g

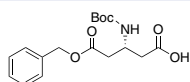
β-Amino Acids and Homologs

Although they are less abundant than their α-analogues, β-amino acids occur in nature both in free form and bound to peptides. Oligomers composed exclusively of β-amino acids (so-called β-peptides) might be the most thoroughly investigated peptidomimetics. Besides being remarkably stable to metabolism, exhibiting slow microbial degradation, and inherently stable to proteases and peptidases, they fold into well-ordered secondary structures consisting of helices, turns, and sheets. In this respect, the most intriguing effects have been observed when β²-amino acids are present in the β-peptide backbone.⁶ A whole new "world" has emerged from the design of fascinating new peptidic macromolecules from β- and γ-homologated proteinogenic amino acids and other components.⁷ Sigma-Aldrich® is proud to be a leading supplier of β-amino acids. This ChemFiles presents a comprehensive list of available products.

References: (6) Lelais, G.; Seebach, D. *Biopolymers* **2004**, 76, 206. (7) Seebach, D.; Beck, A.K.; Bierbaum, D.J. *Chem. & Biodiv.* **2004**, 1, 1111.

Boc-β-Glu(OBzl)-OH, ≥98.0% (HPLC)

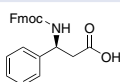
'Boc-L-β-homoaspartic acid 5-benzyl ester';
 Boc-L-β-glutamic acid 5-benzyl ester
 [254101-10-5]
 $C_{17}H_{23}NO_6$
 FW 337.37



03691-1G 1 g

Fmoc-β-D-Phe-OH, ≥98.0% (HPLC)

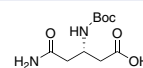
(*S*)-3-(Fmoc-amino)-3-phenylpropionic acid;
 Fmoc-D-β-phenylalanine
 [209252-15-3]
 $C_{24}H_{21}NO_4$
 FW 387.43



00396-100MG 100 mg

Boc-β-Gln-OH, ≥98.0% (TLC)

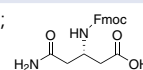
'*N*_ε-Boc-L-β-homoasparagine'; *N*_ε-Boc-L-β-glutamine;
 (*S*)-3-(Boc-amino)pentanoic acid 5-amide
 $C_{10}H_{18}N_2O_5$
 FW 246.26



03651-1G 1 g

Fmoc-β-Gln-OH, ≥97.0% (HPLC)

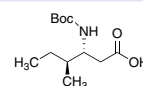
*N*_ε-Fmoc-L-β-homoasparagine; *N*_ε-Fmoc-L-β-glutamine;
 (*S*)-3-(Fmoc-amino)pentanoic acid 5-amide
 [283160-18-9]
 $C_{20}H_{20}N_2O_5$
 FW 368.38



03652-1G 1 g

Boc-β-Homoile-OH, ≥98.0% (TLC)

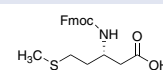
(3*R*,4*S*)-3-(Boc-amino)-4-methylhexanoic acid;
 Boc-L-β-homoisoleucine
 [218608-82-3]
 $C_{12}H_{23}NO_4$
 FW 245.32



03654-250MG 250 mg

Fmoc-β-Homomet-OH, ≥97.0% (HPLC)

(*R*)-3-(Fmoc-amino)-5-(methylthio)pentanoic acid;
 Fmoc-L-β-homomethionine
 [266359-48-2]
 $C_{21}H_{23}NO_4S$
 FW 385.48



03658-1G 1 g

(*S*)-(+)-Pyrrolidine-3-carboxylic acid, ≥98.0% (NT)

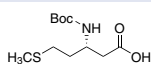
(*S*)-β-Proline
 [72580-53-1]
 $C_5H_9NO_2$
 FW 115.13



670871-500MG 500 mg

Boc-β-Homomet-OH, ≥98.0% (TLC)

(*R*)-3-(Boc-amino)-5-(methylthio)pentanoic acid;
Boc-L-β-homomethionine
[244251-20-5]
C₁₁H₂₁NO₄S
FW 263.35

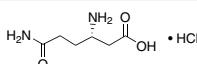


03661-1G

1 g

L-β-Homoglutamine hydrochloride, ≥98.0% (TLC)

(*S*)-3-Aminoadipic acid 6-amide hydrochloride
C₆H₁₂N₂O₃ · HCl
FW 196.63

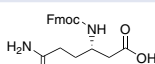


03663-1G

1 g

Fmoc-β-Homogln-OH, ≥98.0% (HPLC)

(*S*)-3-(Fmoc-amino)adipic acid 6-amide;
N_β-Fmoc-L-β-homoglutamine
[283160-17-8]
C₂₁H₂₂N₂O₅
FW 382.41

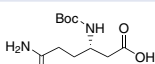


03666-1G

1 g

Boc-β-Homogln-OH, ≥98.0% (TLC)

(*S*)-3-(Boc-amino)adipic acid 6-amide;
N_β-Boc-L-β-homoglutamine
C₁₁H₂₀N₂O₅
FW 260.29

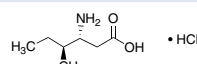


03667-1G

1 g

L-β-Homoisoleucine hydrochloride, ≥98.0% (TLC)

(3*R*,4*S*)-3-Amino-4-methylhexanoic acid hydrochloride
[219310-10-8]
C₇H₁₅NO₂ · HCl
FW 181.66

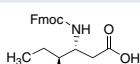


03669-250MG

250 mg

Fmoc-β-Homoile-OH, ≥98.0% (HPLC)

(3*R*,4*S*)-3-(Fmoc-amino)-4-methylhexanoic acid;
Fmoc-L-β-homoisoleucine
[193954-27-7]
C₂₂H₂₅NO₄
FW 367.44

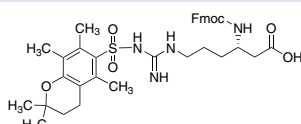


03671-1G

1 g

Fmoc-β-Homoarg(Pmc)-OH, ≥95.0% (HPLC)

N_β-Fmoc-N_ω-(2,2,5,7,8-pentamethylchromane-6-sulfonyl)-L-β-homoarginine
C₃₆H₄₄N₄O₇S
FW 676.82



03673-100MG

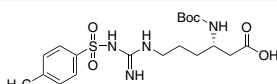
100 mg

03673-500MG

500 mg

Boc-β-Homoarg(Tos)-OH, ≥98.0% (TLC)

N_β-Boc-N_ω-tosyl-L-β-homoarginine
[136271-81-3]
C₁₉H₃₀N₄O₆S
FW 442.53

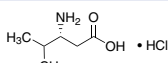


03674-1G

1 g

L-β-Leucine hydrochloride, ≥98.0% (TLC)

'L-β-Homovaline' hydrochloride; (*R*)-3-Amino-4-methylpentanoic acid hydrochloride
[219310-09-5]
C₆H₁₃NO₂ · HCl
FW 167.63



03675-250MG

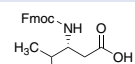
250 mg

03675-1G

1 g

Fmoc-β-Leu-OH, ≥97.0% (HPLC)

Fmoc-L-β-homovaline; Fmoc-L-β-leucine;
(*R*)-3-(Fmoc-amino)-4-methylpentanoic acid
[172695-33-9]
C₂₁H₂₃NO₄
FW 353.41

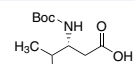


03676-1G

1 g

Boc-β-Leu-OH, ≥98.0% (TLC)

Boc-L-β-homovaline; Boc-L-β-leucine;
(*R*)-3-(Boc-amino)-4-methylpentanoic acid
[183990-64-9]
C₁₁H₂₁NO₄
FW 231.29

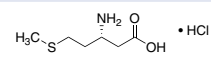


03678-1G

1 g

L-β-Homomethionine hydrochloride, ≥98.0% (TLC)

R-3-Amino-5-(methylthio)pentanoic acid hydrochloride
C₆H₁₃NO₂S · HCl
FW 199.70

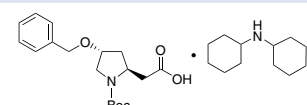


03681-1G

1 g

Boc-β-Homohyp(Bzl)-OH (dicyclohexylammonium) salt, ≥85% (HPLC)

Boc-*O*-benzyl-L-β-homohydroxyproline (dicyclohexylammonium) salt;
(2*S*,4*R*)-1-Boc-4-benzyloxy-pyrrolidine-2-acetic acid (dicyclohexylammonium) salt
C₁₈H₂₅NO₅ · C₁₂H₂₃N
FW 516.71

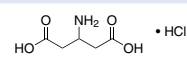


03683-1G

1 g

β-Glutamic acid hydrochloride, ≥98.0% (TLC)

3-Aminoglutaric acid hydrochloride;
'β-Homoaspartic acid' hydrochloride
C₅H₉NO₄ · HCl
FW 183.59

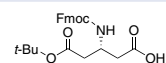


03688-500MG

500 mg

Fmoc-β-Glu(OtBu)-OH, ≥98.0% (HPLC)

Fmoc-L-β-glutamic acid 5-*tert*-butyl ester;
Fmoc-L-β-homoaspartic acid 5-*tert*-butyl ester
[209252-17-5]
C₂₄H₂₇NO₆
FW 425.47

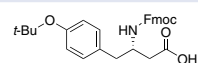


03689-1G

1 g

Fmoc-β-Homotyr(tBu)-OH, ≥98.0% (HPLC)

Fmoc-*O*-*tert*-butyl-L-β-homotyrosine
[219967-69-8]
C₂₉H₃₁NO₅
FW 473.56

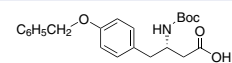


03692-1G

1 g

Boc-β-Homotyr(Bzl)-OH, ≥97.0% (HPLC)

Boc-*O*-benzyl-L-β-homotyrosine
[126825-16-9]
C₂₂H₂₇NO₅
FW 385.45

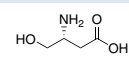


03693-1G

1 g

L-β-Homoserine, ≥98.0% (TLC)

(*R*)-3-Amino-4-hydroxybutyric acid
[16504-56-6]
C₄H₉NO₃
FW 119.12



03694-100MG

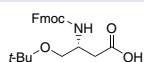
100 mg

03694-500MG

500 mg

Fmoc-β-Homoser(tBu)-OH, ≥97.5% (HPLC)

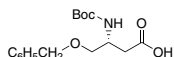
Fmoc-*O*-*tert*-butyl-L-β-homoserine
[203854-51-7]
C₂₃H₂₇NO₅
FW 397.46



03696-1G 1 g

Boc-β-Homoser(Bzl)-OH, ≥98.0% (HPLC)

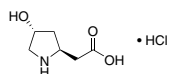
Boc-*O*-benzyl-L-β-homoserine
[218943-31-8]
C₁₆H₂₃NO₅
FW 309.36



03697-1G 1 g

L-β-Homohydroxyproline hydrochloride, ≥98.0% (TLC)

(2*S*,4*R*)-4-Hydroxy-2-pyrrolidinylacetic acid hydrochloride
C₆H₁₁NO₃ · HCl
FW 181.62

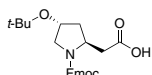


03698-100MG 100 mg

03698-500MG 500 mg

Fmoc-β-Homohyp(tBu)-OH, ≥98.0% (HPLC)

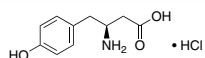
Fmoc-*O*-*tert*-butyl-L-β-homohydroxyproline
[957509-29-4]
C₂₅H₂₉NO₅
FW 423.50



03751-500MG 500 mg

L-β-Homotyrosine hydrochloride, ≥97.0% (HPLC)

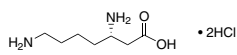
(*S*)-3-Amino-4-(4-hydroxyphenyl)butyric acid hydrochloride
C₁₀H₁₃NO₃ · HCl
FW 231.68



03758-100MG 100 mg

L-β-Homolysine dihydrochloride, ≥98.0% (TLC)

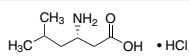
(*S*)-3,7-Diaminoheptanoic acid dihydrochloride
C₇H₁₆N₂O₂ · 2HCl
FW 233.14



03759-100MG 100 mg

L-β-Homoleucine hydrochloride, ≥98.0% (TLC)

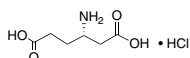
(*S*)-3-Amino-5-methylhexanoic acid hydrochloride
[96386-92-4]
C₇H₁₅NO₂ · HCl
FW 181.66



03764-1G 1 g

L-β-Homoglutamic acid hydrochloride, ≥98.0% (TLC)

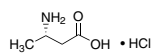
(*S*)-3-Amino adipic acid hydrochloride;
(*S*)-3-Aminohexanedioic acid hydrochloride
[61884-74-0]
C₆H₁₁NO₄ · HCl
FW 197.62



03765-100MG 100 mg

L-β-Homoalanine hydrochloride, ≥98.0% (TLC)

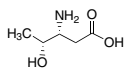
(*S*)-3-Aminobutyric acid hydrochloride
[58610-41-6]
C₄H₉NO₂ · HCl
FW 139.58



03766-250MG 250 mg

L-β-Homothreonine, ≥98.0% (TLC)

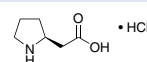
(3*R*,4*R*)-3-Amino-4-hydroxypentanoic acid
C₅H₁₁NO₃
FW 133.15



03767-100MG 100 mg

'L-β-Homoproline' hydrochloride, ≥98.0% (TLC)

(*S*)-2-(2-Pyrrolidinyl)acetic acid hydrochloride
[53912-85-9]
C₆H₁₁NO₂ · HCl
FW 165.62

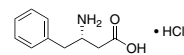


03768-250MG 250 mg

03768-1G 1 g

L-β-Homophenylalanine hydrochloride, ≥98.0% (TLC)

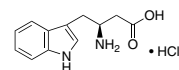
(*S*)-3-Amino-4-phenylbutyric acid hydrochloride
[138165-77-2]
C₁₀H₁₃NO₂ · HCl
FW 215.68



03769-250MG 250 mg

L-β-Homotryptophan hydrochloride, ≥98.0% (TLC)

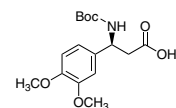
(*S*)-3-Amino-4-(3-indolyl)butyric acid hydrochloride
C₁₂H₁₄N₂O₂ · HCl
FW 254.71



03790-100MG 100 mg

(*S*)-Boc-3,4-dimethoxy-β-Phe-OH, ≥98.0% (HPLC)

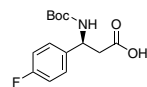
(*S*)-3-(Boc-amino)-3-(3,4-methoxyphenyl)propionic acid;
Boc-3,4-dimethoxy-D-β-phenylalanine
[499995-84-5]
C₁₆H₂₃NO₆
FW 325.36



04542-500MG-F 500 mg

(*S*)-Boc-4-fluoro-β-Phe-OH, ≥98.0% (HPLC)

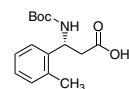
(*S*)-3-(Boc-amino)-3-(4-fluorophenyl)propionic acid;
Boc-4-fluoro-D-β-phenylalanine
[479064-88-5]
C₁₄H₁₈FO₄
FW 283.30



05102-500MG-F 500 mg

(*R*)-Boc-2-methyl-β-Phe-OH, ≥98.0% (HPLC)

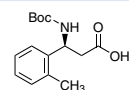
(*R*)-3-(Boc-amino)-3-(2-methylphenyl)propionic acid;
Boc-2-methyl-L-β-phenylalanine
[500770-86-5]
C₁₅H₂₁NO₄
FW 279.33



06731-500MG-F 500 mg

(*S*)-Boc-2-methyl-β-Phe-OH, ≥98.0% (HPLC)

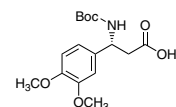
(*S*)-3-(Boc-amino)-3-(2-methylphenyl)propionic acid;
Boc-2-methyl-D-β-phenylalanine
[499995-74-3]
C₁₅H₂₁NO₄
FW 279.33



06939-500MG-F 500 mg

(*R*)-Boc-3,4-dimethoxy-β-Phe-OH, ≥98.0% (HPLC)

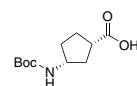
(*R*)-3-(Boc-amino)-3-(3,4-methoxyphenyl)propionic acid;
Boc-3,4-dimethoxy-L-β-phenylalanine
[500788-93-2]
C₁₆H₂₃NO₆
FW 325.36



08009-500MG-F 500 mg

(1*S*,3*R*)-(+)-3-(Boc-amino)cyclopentanecarboxylic acid, ≥96.0% (TLC)

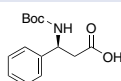
[261165-05-3]
C₁₁H₁₉NO₄
FW 229.27



09784-500MG 500 mg

Boc-D-β-Phe-OH, ≥98.0% (HPLC)

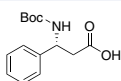
(S)-3-(Boc-amino)-3-phenylpropionic acid;
Boc-D-β-phenylalanine
[103365-47-5]
C₁₄H₁₉NO₄
FW 265.30



09793-1G 1 g

Boc-β-Phe-OH, ≥96.0% (HPLC)

Boc-L-β-phenylalanine;
(R)-3-(Boc-amino)-3-phenylpropionic acid
[161024-80-2]
C₁₄H₁₉NO₄
FW 265.30

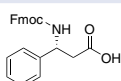


09794-250MG 250 mg

09794-1G 1 g

Fmoc-β-Phe-OH, ≥98.0% (HPLC)

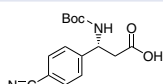
(R)-3-(Fmoc-amino)-3-phenylpropionic acid;
Fmoc-L-β-phenylalanine
[220498-02-2]
C₂₄H₂₁NO₄
FW 387.43



09795-500MG 500 mg

(R)-Boc-4-cyano-β-Phe-OH, ≥98.0% (HPLC)

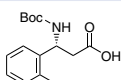
(R)-3-(Boc-amino)-3-(4-cyanophenyl)propionic acid;
Boc-4-cyano-L-β-phenylalanine
[501015-22-1]
C₁₅H₁₈N₂O₄
FW 290.31



12339-500MG-F 500 mg

(R)-Boc-2-(trifluoromethyl)-β-Phe-OH, ≥98.0% (HPLC)

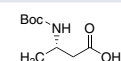
(R)-3-(Boc-amino)-3-[2-(trifluoromethyl)phenyl]propionic acid;
Boc-2-(trifluoromethyl)-L-β-phenylalanine
[501015-17-4]
C₁₅H₁₈F₃NO₄
FW 333.30



14670-500MG-F 500 mg

Boc-β-Homoala-OH, ≥98.0% (TLC)

(S)-3-(Boc-amino)butyric acid; Boc-L-β-homoalanine
[158851-30-0]
C₉H₁₇NO₄
FW 203.24

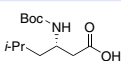


14974-250MG 250 mg

14974-1G 1 g

Boc-β-Homoleu-OH, ≥98.0% (TLC)

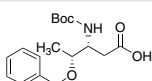
(S)-3-(Boc-amino)-5-methylhexanoic acid;
Boc-L-β-homoleucine
[132549-43-0]
C₁₂H₂₃NO₄
FW 245.32



14975-1G 1 g

Boc-β-Homothr(Bzl)-OH, ≥98.0% (TLC)

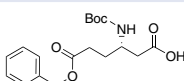
(3R,4R)-4-Benzoyloxy-3-(Boc-amino)pentanoic acid;
Boc-O-benzyl-L-β-homothreonine
C₁₇H₂₅NO₅
FW 323.38



14976-1G 1 g

Boc-β-Homoglu(OBzl)-OH, ≥98.0% (TLC)

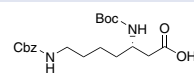
(S)-3-(Boc-amino)adipic acid 6-benzyl ester;
Boc-L-β-homoglutamic acid 6-benzyl ester
[218943-30-7]
C₁₈H₂₅NO₆
FW 351.39



14977-1G 1 g

Boc-β-Homolys(Z)-OH, ≥98.0% (TLC)

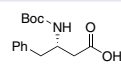
(S)-3-(Boc-amino)-7-(Z-amino)heptanoic acid;
N_β-Boc-N_ω-Z-L-β-homolysine
C₂₀H₃₀N₂O₆
FW 394.46



14978-1G 1 g

Boc-β-Homophe-OH, ≥98.0% (TLC)

(S)-3-(Boc-amino)-4-phenylbutyric acid;
Boc-L-β-homophenylalanine
[51871-62-6]
C₁₅H₂₁NO₄
FW 279.33

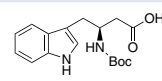


14979-250MG 250 mg

14979-1G 1 g

Boc-β-Homotrp-OH, ≥98.0% (TLC)

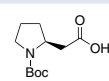
(S)-3-(Boc-amino)-4-(3-indolyl)butyric acid;
N_β-Boc-L-β-homotryptophan
[229639-48-9]
C₁₇H₂₂N₂O₄
FW 318.37



14981-250MG 250 mg

Boc-β³-Homopro-OH, ≥98.0% (TLC)

(S)-2-(1-Boc-2-pyrrolidinyl)acetic acid; Boc-L-β³-homoproline
[56502-01-3]
C₁₁H₁₉NO₄
FW 229.27

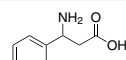


14982-250MG 250 mg

14982-1G 1 g

DL-β-Phenylalanine, 98%

DL-β-Homophenylglycine; ±-3-Amino-3-phenylpropionic acid;
β-Aminohydrocinnamic acid; DL-3-Amino-3-phenylpropionic acid
[614-19-7]
C₉H₁₁NO₂
FW 165.19



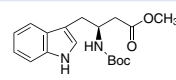
159492-1G 1 g

159492-5G 5 g

159492-25G 25 g

Boc-β-Homotrp-OMe, ≥98.0% (TLC)

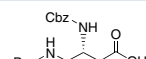
N_β-Boc-L-β-homotryptophan methyl ester
C₁₈H₂₄N₂O₄
FW 332.39



17008-250MG 250 mg

Z-β-Dab(Boc)-OH, ≥98.0% (TLC)

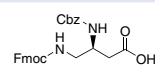
Z-β-DBu(Boc)-OH; N_β-Z-N_γ-Boc-L-3,4-diaminobutyric acid;
(R)-4-(Boc-amino)-3-(Z-amino)butyric acid;
N_γ-Boc-N_β-Z-L-3,4-diaminobutyric acid
[108919-51-3]
C₁₇H₂₄N₂O₆
FW 352.38



17974-1G 1 g

Z-D-β-Dab(Fmoc)-OH, ≥98.0% (HPLC)

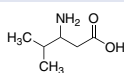
β-Z-γ-Fmoc-D-3,4-diaminobutyric acid; (S)-4-(Fmoc-amino)-3-(Z-amino)butyric acid; γ-Fmoc-β-Z-D-3,4-diaminobutyric acid; Z-D-β-DBu(Fmoc)-OH
C₂₇H₂₆N₂O₆
FW 474.51



17984-1G 1 g

DL-β-Leucine, ≥98.0% (NT)

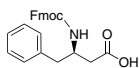
(±)-3-Amino-4-methylpentanoic acid;
(±)-3-Amino-4-methylvaleric acid
[5699-54-7]
C₆H₁₃NO₂
FW 131.17



17988-1G-F 1 g

Fmoc-D-β-Homophe-OH, ≥97.0% (HPLC)

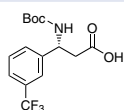
Fmoc-D-β-homophenylalanine;
(R)-3-(Fmoc-amino)-4-phenylbutyric acid
[209252-16-4]
C₂₅H₂₃NO₄
FW 401.45



18074-500MG-F 500 mg

(R)-Boc-3-(trifluoromethyl)-β-Phe-OH, ≥95.0% (HPLC)

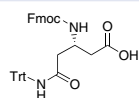
(R)-3-(Boc-amino)-3-[3-(trifluoromethyl)phenyl]propionic acid;
Boc-3-(trifluoromethyl)-L-β-phenylalanine
[501015-18-5]
C₁₅H₁₈F₃NO₄
FW 333.30



18328-500MG-F 500 mg

Fmoc-β-Gln(Trt)-OH, ≥96.0% (HPLC)

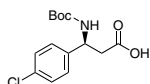
Fmoc-β-Homoasn(Trt)-OH; N_β-Fmoc-N₆-trityl-L-homoasparagine; (S)-3-(Fmoc-amino)-N-trityl-valeric acid 5-amide
[283160-20-3]
C₃₉H₃₄N₂O₅
FW 610.70



18505-1G-F 1 g

(S)-Boc-4-chloro-β-Phe-OH, ≥98.0% (HPLC)

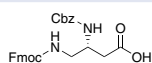
(S)-3-(Boc-amino)-3-(4-chlorophenyl)propionic acid;
Boc-4-chloro-D-β-phenylalanine
[479064-90-9]
C₁₄H₁₈ClNO₄
FW 299.75



19343-500MG-F 500 mg

Z-β-Dab(Fmoc)-OH, ≥98.0% (HPLC)

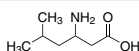
Z-β-Dab(Fmoc)-OH; (R)-4-(Fmoc-amino)-3-(Z-amino)butyric acid; γ-Fmoc-β-Z-L-3,4-diaminobutyric acid;
β-Z-γ-Fmoc-L-3,4-diaminobutyric acid
[349547-09-7]
C₂₇H₂₆N₂O₆
FW 474.51



21622-1G 1 g

DL-β-Homoleucine, ≥99.0% (NT)

(±)-3-Amino-5-methylcaproic acid;
(±)-3-Amino-5-methylhexanoic acid
[3653-34-7]
C₇H₁₅NO₂
FW 145.20

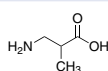


21625-1G 1 g

21625-5G 5 g

DL-3-Aminoisobutyric acid, 98%

α-Methyl-β-alanine
[144-90-1]
C₄H₉NO₂
FW 103.12

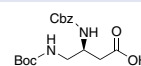


217794-1G 1 g

217794-5G 5 g

Z-D-β-Dab(Boc)-OH, ≥98.0% (TLC)

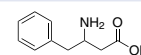
Z-D-β-Dab(Boc)-OH; (S)-4-(Boc-amino)-3-(Z-amino)butyric acid; N_β-Boc-N_γ-Z-D-3,4-diaminobutyric acid;
N_β-Z-N_γ-Boc-D-3,4-diaminobutyric acid
[96186-30-0]
C₁₇H₂₄N₂O₆
FW 352.38



28206-250MG 250 mg

DL-β-Homophenylalanine, ≥98.0% (HPLC)

(±)-3-Amino-4-phenylbutyric acid
[15099-85-1]
C₁₀H₁₃NO₂
FW 179.22

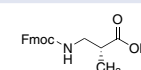


28217-1G-F 1 g

28217-5G-F 5 g

(R)-3-(Fmoc-amino)-2-methylpropionic acid, ≥95.0% (HPLC)

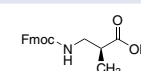
(R)-Fmoc-β²-Homoala-OH
[211682-15-4]
C₁₉H₁₉NO₄
FW 325.36



30975-500MG-F 500 mg

(S)-3-(Fmoc-amino)-2-methylpropionic acid, ≥95.0% (HPLC)

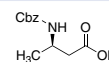
(S)-Fmoc-β²-Homoala-OH
[203854-58-4]
C₁₉H₁₉NO₄
FW 325.36



38811-500MG-F 500 mg

Z-β-D-Homoala-OH, ≥98.0% (HPLC)

Z-β-D-homoalanine; (R)-3-(Z-amino)butyric acid
[67843-72-5]
C₁₂H₁₅NO₄
FW 237.25

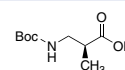


39748-250MG-F 250 mg

39748-1G-F 1 g

(S)-3-(Boc-amino)-2-methylpropionic acid, ≥98.0% (TLC)

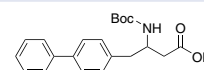
(S)-Boc-β²-Homoala-OH
[190897-47-3]
C₉H₁₇NO₄
FW 203.24



40195-500MG-F 500 mg

(±)-3-(Boc-amino)-4-(4-biphenyl)butyric acid, ≥98.0% (HPLC)

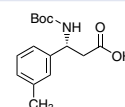
Boc-4'-phenyl-DL-β-homophenylalanine
[683219-74-1]
C₂₁H₂₅NO₄
FW 355.43



41844-500MG-F 500 mg

(R)-Boc-3-methyl-β-Phe-OH, ≥98.0% (HPLC)

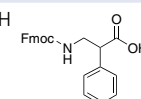
(R)-3-(Boc-amino)-3-(3-methylphenyl)propionic acid;
Boc-3-methyl-L-β-phenylalanine
[464930-76-5]
C₁₅H₂₁NO₄
FW 279.33



42722-500MG-F 500 mg

(±)-3-(Fmoc-amino)-2-phenylpropionic acid, ≥95.0% (HPLC)

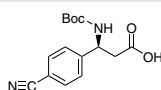
(±)-Fmoc-2-phenyl-β-alanine; Fmoc-β²-DL-Homophe-OH
[683217-60-9]
C₂₄H₂₁NO₄
FW 387.43



43026-500MG-F 500 mg

(S)-Boc-4-cyano-β-Phe-OH, ≥98.0% (HPLC)

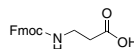
(S)-3-(Boc-amino)-3-(4-cyanophenyl)propionic acid;
Boc-4-cyano-D-β-phenylalanine
[500770-82-1]
C₁₅H₁₈N₂O₄
FW 290.31



44376-500MG-F 500 mg

Fmoc-β-Ala-OH, ≥99.0% (HPLC)

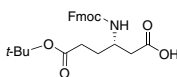
Fmoc-β-alanine
[35737-10-1]
C₁₈H₁₇NO₄
FW 311.33



47587-1G-F 1 g
47587-5G-F 5 g
47587-25G-F 25 g

Fmoc-β-Homoglu(OtBu)-OH, ≥98.0% (HPLC)

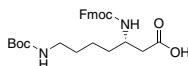
(S)-3-(Fmoc-amino)adipic acid 6-*tert*-butyl ester;
Fmoc-L-β-homoglutamic acid 6-*tert*-butyl ester
[203854-49-3]
C₂₅H₂₉NO₆
FW 439.50



47837-1G 1 g

Fmoc-β-Homolys(Boc)-OH, ≥95.0% (TLC)

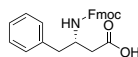
(S)-7-(Boc-amino)-3-(Fmoc-amino)heptanoic acid;
N_β-Fmoc-N_ω-Boc-L-β-homolysine
[203854-47-1]
C₂₇H₃₄N₂O₆
FW 482.57



47874-1G 1 g

Fmoc-β-Homophe-OH, ≥95.0% (HPLC)

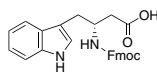
(S)-3-(Fmoc-amino)-4-phenylbutyric acid;
Fmoc-L-β-homophenylalanine
[193954-28-8]
C₂₅H₂₃NO₄
FW 401.45



47878-1G-F 1 g

Fmoc-β-Homotrp-OH, ≥96.0% (HPLC)

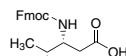
(S)-3-(Fmoc-amino)-4-(3-indolyl)butyric acid;
N_β-Fmoc-L-β-homotryptophan
C₂₇H₂₄N₂O₄
FW 440.49



47901-1G 1 g

Fmoc-β-Homotr(tBu)-OH, ≥96.0% (HPLC)

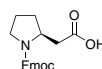
(3*R*,4*R*)-4-*tert*-Butoxy-3-(Fmoc-amino)pentanoic acid;
Fmoc-*O*-*tert*-butyl-L-β-homothreonine
C₂₄H₂₉NO₅
FW 411.49



47911-1G 1 g

Fmoc-L-β³-homoproline, ≥98.0% (HPLC)

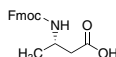
(S)-2-(1-Fmoc-2-pyrrolidinyl)acetic acid
[193693-60-6]
C₂₁H₂₁NO₄
FW 351.40



47912-1G 1 g

Fmoc-β-Homoala-OH, ≥98.0% (HPLC)

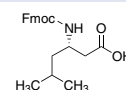
(S)-3-(Fmoc-amino)butyric acid; Fmoc-L-β-homoalanine
[193954-26-6]
C₁₉H₁₉NO₄
FW 325.36



47935-1G 1 g

Fmoc-β-Homoleu-OH, ≥96.0% (HPLC)

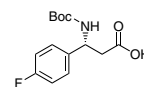
Fmoc-L-β-homoleucine; (S)-3-(Fmoc-amino)-5-methylhexanoic acid
[193887-44-4]
C₂₂H₂₅NO₄
FW 367.44



47946-250MG 250 mg
47946-1G 1 g

(R)-Boc-4-fluoro-β-Phe-OH, ≥98.0% (HPLC)

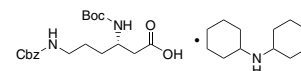
(R)-3-(Boc-amino)-3-(4-fluorophenyl)propionic acid;
Boc-4-fluoro-L-β-phenylalanine
[479064-94-3]
C₁₄H₁₈FO₄
FW 283.30



51006-500MG-F 500 mg

Boc-β-Lys(Z)-OH (dicyclohexylammonium) salt, ≥98.0% (TLC)

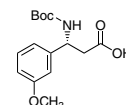
Boc-L-β-Homoorn(Z)-OH
(dicyclohexylammonium) salt;
N_β-Boc-N_ω-Z-L-β-lysine
(dicyclohexylammonium) salt
C₁₉H₂₈N₂O₆ · C₁₂H₂₃N
FW 561.75



52136-1G 1 g

(R)-Boc-3-methoxy-β-Phe-OH, ≥98.0% (HPLC)

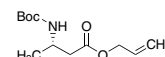
(R)-3-(Boc-amino)-3-(3-methoxyphenyl)propionic acid;
Boc-3-methoxy-L-β-phenylalanine
[500788-86-3]
C₁₅H₂₁NO₅
FW 295.33



52169-500MG-F 500 mg

Boc-β-Homoala-OAlI, ≥98.0% (TLC)

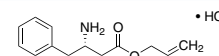
(S)-3-(Boc-amino)butyric acid allyl ester;
Boc-L-β-homoalanine allyl ester
C₁₂H₂₁NO₄
FW 243.30



53458-1G 1 g

L-β-Homophenylalanine allyl ester hydrochloride, ≥98.0% (TLC)

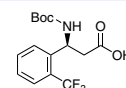
Allyl (S)-3-amino-4-phenylbutyrate
hydrochloride
C₁₃H₁₇NO₂ · HCl
FW 255.74



53575-1G 1 g

(S)-Boc-2-(trifluoromethyl)-β-Phe-OH, ≥98.0% (HPLC)

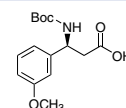
(S)-3-(Boc-amino)-3-[2-(trifluoromethyl)phenyl]propionic acid;
Boc-2-(trifluoromethyl)-D-β-phenylalanine
[500770-77-4]
C₁₅H₁₈F₃NO₄
FW 333.30



55685-500MG-F 500 mg

(S)-Boc-3-methoxy-β-Phe-OH, ≥98.0% (HPLC)

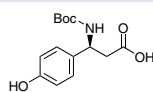
(S)-3-(Boc-amino)-3-(3-methoxyphenyl)propionic acid;
Boc-3-methoxy-D-β-phenylalanine
[499995-77-6]
C₁₅H₂₁NO₅
FW 295.33



55894-500MG-F 500 mg

(S)-Boc-β-Tyr-OH, ≥98.0% (HPLC)

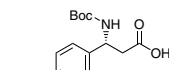
(S)-3-(Boc-amino)-3-(4-hydroxyphenyl)propionic acid;
Boc-D-β-tyrosine
[499995-80-1]
C₁₄H₁₉NO₅
FW 281.30



56996-500MG-F 500 mg

(R)-Boc-4-methoxy-β-Phe-OH, ≥98.0% (HPLC)

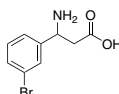
(R)-3-(Boc-amino)-3-(4-methoxyphenyl)propionic acid;
Boc-β-Tyr(Me)-OH; Boc-4-methoxy-L-β-phenylalanine;
Boc-O-methyl-L-β-tyrosine
[500788-87-4]
C₁₅H₂₁NO₅
FW 295.33



56999-500MG-F 500 mg

3-Amino-3-(3-bromophenyl)propionic acid, 97%

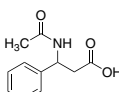
[117391-50-1]
C₉H₁₀BrNO₂
FW 244.09



573450-1G 1 g

N-Acetyl-DL-β-phenylalanine, ≥97.0% (HPLC)

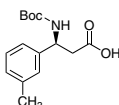
(±)-3-(Acetylamino)-3-phenylpropionic acid
[40638-98-0]
C₁₁H₁₃NO₃
FW 207.23



59891-250MG-F 250 mg

(S)-Boc-3-methyl-β-Phe-OH, ≥98.0% (HPLC)

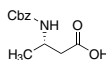
(S)-3-(Boc-amino)-3-(3-methylphenyl)propionic acid;
Boc-3-methyl-D-β-phenylalanine
[499995-75-4]
C₁₅H₂₁NO₄
FW 279.33



59967-500MG-F 500 mg

Z-β-Homoala-OH, ≥98.0% (HPLC)

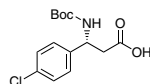
Z-L-β-Homoalanine; (S)-3-(Z-amino)butyric acid
[83509-88-0]
C₁₂H₁₅NO₄
FW 237.25



61669-500MG-F 500 mg

(R)-Boc-4-chloro-β-Phe-OH, ≥98.0% (HPLC)

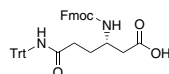
(R)-3-(Boc-amino)-3-(4-chlorophenyl)propionic acid;
Boc-4-chloro-L-β-phenylalanine
[479064-93-2]
C₁₄H₁₈ClNO₄
FW 299.75



61902-500MG-F 500 mg

Fmoc-β-Homogln(Trt)-OH, ≥95.0% (HPLC)

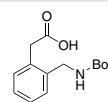
N_β-Fmoc-N_ε-trityl-L-homoglutamine;
(S)-3-(Fmoc-amino)-N-trityl-adipic acid 6-amide
[401915-55-7]
C₄₀H₃₆N₂O₅
FW 624.72



64179-250MG-F 250 mg

2-(Boc-aminomethyl)phenylacetic acid, ≥99.0%

[40851-66-9]
C₁₄H₁₉NO₄
FW 265.30



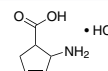
≥99.0% (HPLC)
99.0-101.0% (NaOH 0.1 M, T)

670758-100MG 100 mg

670758-500MG 500 mg

cis-2-Amino-3-cyclopentene-1-carboxylic acid hydrochloride, ≥97.0% (AT)

[122022-92-8]
C₆H₉NO₂ · HCl
FW 163.60

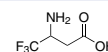


68602-250MG 250 mg

68602-1G 1 g

3-Amino-4,4,4-trifluorobutyric acid, ≥98.0% (TLC)

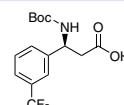
4,4,4-Trifluoro-β-homoalanine
[584-20-3]
C₄H₆F₃NO₂
FW 157.09



69289-500MG-F 500 mg

(S)-Boc-3-(trifluoromethyl)-β-Phe-OH, ≥98.0% (HPLC)

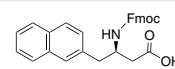
(S)-3-(Boc-amino)-3-[3-(trifluoromethyl)phenyl]propionic acid;
Boc-3-(trifluoromethyl)-D-β-phenylalanine
[500770-78-5]
C₁₅H₁₈F₃NO₄
FW 333.30



69418-500MG-F 500 mg

(R)-3-(Fmoc-amino)-4-(2-naphthyl)butyric acid, ≥95.0% (HPLC)

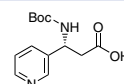
Fmoc-D-β-2-Homonal-OH;
Fmoc-4-(2-naphthyl)-D-β-homoalanine
[269398-91-6]
C₂₉H₂₅NO₄
FW 451.51



70796-500MG-F 500 mg

(R)-Boc-3-(3-pyridyl)-β-Ala-OH, ≥96.0% (HPLC)

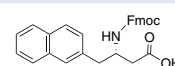
L-Boc-3-(3-pyridyl)-β-alanine;
(R)-3-(Boc-amino)-3-(3-pyridyl)propionic acid
[500788-96-5]
C₁₃H₁₈N₂O₄
FW 266.29



72722-500MG-F 500 mg

(S)-3-(Fmoc-amino)-4-(2-naphthyl)butyric acid, ≥97.0% (HPLC)

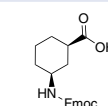
Fmoc-β-2-Homonal-OH;
Fmoc-4-(2-naphthyl)-L-β-homoalanine
[270063-40-6]
C₂₉H₂₅NO₄
FW 451.51



72829-500MG-F 500 mg

cis-3-(Fmoc-amino)cyclohexanecarboxylic acid, ≥98.0% (HPLC)

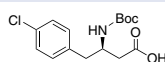
C₂₂H₂₃NO₄
FW 365.42



73024-500MG-F 500 mg

(R)-Boc-4-chloro-β-Homophe-OH, ≥98.0% (HPLC)

(R)-3-(Boc-amino)-4-(4-chlorophenyl)butyric acid;
Boc-4-chloro-D-β-homophenylalanine
[218608-96-9]
C₁₅H₂₀ClNO₄
FW 313.78

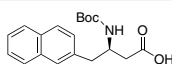


73086-500MG-F

500 mg

(R)-3-(Boc-amino)-4-(2-naphthyl)butyric acid, ≥98.0% (HPLC)

Boc-D-β-2-Homonal-OH;
Boc-4-(2-naphthyl)-D-β-homoalanine
[219297-10-6]
C₁₉H₂₃NO₄
FW 329.39

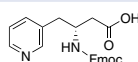


73300-500MG-F

500 mg

(R)-Fmoc-4-(3-pyridyl)-β-Homoala-OH, ≥98.0% (HPLC)

(R)-3-(Fmoc-amino)-4-(3-pyridyl)butyric acid;
Fmoc-4-(3-pyridyl)-D-β-homoalanine
[269396-66-9]
C₂₄H₂₂N₂O₄
FW 402.44

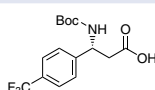


73309-500MG-F

500 mg

(R)-Boc-4-(trifluoromethyl)-β-Phe-OH, ≥98.0% (HPLC)

(R)-3-(Boc-amino)-3-[4-(trifluoromethyl)phenyl]propionic acid;
Boc-4-(trifluoromethyl)-L-β-phenylalanine
[501015-19-6]
C₁₅H₁₈F₃NO₄
FW 333.30

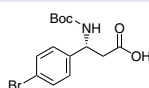


73999-500MG-F

500 mg

(R)-Boc-4-bromo-β-Phe-OH, ≥98.0% (HPLC)

(R)-3-(Boc-amino)-3-(4-bromophenyl)propionic acid;
Boc-4-bromo-L-β-phenylalanine
[261380-20-5]
C₁₄H₁₆BrNO₄
FW 344.20

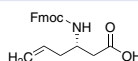


74037-500MG-F

500 mg

(S)-3-(Fmoc-amino)-5-hexenoic acid, ≥95.0% (HPLC)

Fmoc-4-vinyl-L-β-homoalanine
[270263-04-2]
C₂₁H₂₁NO₄
FW 351.40

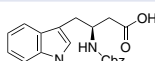


75561-500MG-F

500 mg

Z-β-Homotrp(Boc)-OH, ≥97.0% (HPLC)

N_B-Z-N_{in}-Boc-L-β-homotryptophan;
N_{in}-Boc-N_B-Z-L-β-homotryptophan
C₂₅H₂₈N₂O₆
FW 452.50

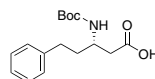


76024-500MG-F

500 mg

(S)-3-(Boc-amino)-5-phenylpentanoic acid, ≥97.0% (C)

Boc-5-phenyl-L-β-norvaline
[218608-84-5]
C₁₆H₂₃NO₄
FW 293.36

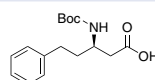


76144-500MG-F

500 mg

(R)-3-(Boc-amino)-5-phenylpentanoic acid, ≥97.0% (GC)

Boc-5-phenyl-D-β-norvaline
[218608-83-4]
C₁₆H₂₃NO₄
FW 293.36

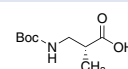


78018-500MG-F

500 mg

(R)-3-(Boc-amino)-2-methylpropionic acid, ≥98.0% (TLC)

(R)-Boc-β²-Homoala-OH
[132696-45-8]
C₉H₁₇NO₄
FW 203.24

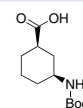


78953-500MG-F

500 mg

cis-3-(Boc-amino)cyclohexanecarboxylic acid, ≥98.0% (TLC)

[222530-33-8]
C₁₂H₂₁NO₄
FW 243.30

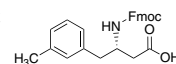


80393-500MG-F

500 mg

(S)-Fmoc-3-methyl-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Fmoc-amino)-4-(3-methylphenyl)butyric acid;
Fmoc-3-methyl-L-β-homophenylalanine
[270062-94-7]
C₂₆H₂₅NO₄
FW 415.48

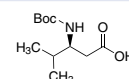


80587-500MG-F

500 mg

Boc-D-β-Leu-OH, ≥98.0% (TLC)

(S)-3-(Boc-amino)-4-methylpentanoic acid;
Boc-D-β-Homoval-OH; Boc-D-β-Leucine
[179412-79-4]
C₁₁H₂₁NO₄
FW 231.29

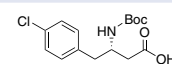


80674-250MG

250 mg

(S)-Boc-4-chloro-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-4-(4-chlorophenyl)butyric acid;
Boc-4-chloro-L-β-homophenylalanine
[270596-42-4]
C₁₅H₂₀ClNO₄
FW 313.78

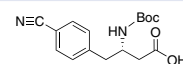


81771-500MG-F

500 mg

(S)-Boc-4-cyano-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-4-(4-cyanophenyl)butyric acid;
Boc-4-cyano-L-β-homophenylalanine
[270065-89-9]
C₁₆H₂₀N₂O₄
FW 304.34

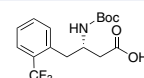


81834-500MG-F

500 mg

(S)-Boc-2-(trifluoromethyl)-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-4-[2-(trifluoromethyl)phenyl]butyric acid;
Boc-2-(trifluoromethyl)-L-β-homophenylalanine
[270065-74-2]
C₁₆H₂₀F₃NO₄
FW 347.33

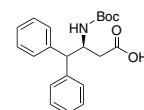


81998-500MG-F

500 mg

(S)-Boc-γ,γ-diphenyl-β-Homoala-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-4,4-diphenylbutyric acid;
Boc-γ,γ-diphenyl-D-β-homoalanine
[190190-50-2]
C₂₁H₂₅NO₄
FW 355.43

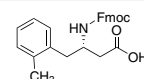


82007-500MG-F

500 mg

(S)-Fmoc-2-methyl-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Fmoc-amino)-4-(2-methylphenyl)butyric acid;
Fmoc-2-methyl-L-β-homophenylalanine
[270062-91-4]
C₂₆H₂₅NO₄
FW 415.48

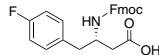


85497-500MG-F

500 mg

(R)-Fmoc-4-fluoro-β-Homophe-OH, ≥98.0% (HPLC)

(R)-3-(Fmoc-amino)-4-(4-fluorophenyl)butyric acid;
Fmoc-4-fluoro-D-β-homophenylalanine
[331763-70-3]
C₂₅H₂₂FNO₄
FW 419.44

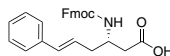


87084-500MG-F

500 mg

(S)-3-(Fmoc-amino)-6-phenyl-5-hexenoic acid, ≥97.0% (HPLC)

Fmoc-4-styryl-L-β-homoalanine
[270596-45-7]
C₂₇H₂₅NO₄
FW 427.49

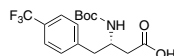


87901-500MG-F

500 mg

(S)-Boc-2-(trifluoromethyl)-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-4-[4-(trifluoromethyl)phenyl]butyric acid; Boc-4-(trifluoromethyl)-L-β-homophenylalanine
[270065-80-0]
C₁₆H₂₀F₃NO₄
FW 347.33

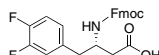


87943-500MG-F

500 mg

(S)-Fmoc-3,4-difluoro-β-Homophe-OH, ≥98.0% (HPLC)

Fmoc-3,4-difluoro-L-β-homophenylalanine;
(S)-3-(Fmoc-amino)-4-(3,4-difluorophenyl)butyric acid
[270063-55-3]
C₂₅H₂₁F₂NO₄
FW 437.44

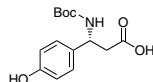


88574-500MG-F

500 mg

(R)-Boc-β-Tyr-OH, ≥97.0% (HPLC)

(R)-3-(Boc-amino)-3-(4-hydroxyphenyl)propionic acid;
Boc-L-β-tyrosine
[329013-12-9]
C₁₄H₁₅NO₅
FW 281.30

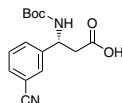


89090-500MG-F

500 mg

(R)-Boc-3-cyano-β-Phe-OH, ≥98.0% (HPLC)

(R)-3-(Boc-amino)-3-(3-cyanophenyl)propionic acid;
Boc-3-cyano-L-β-phenylalanine
[501015-21-0]
C₁₅H₁₃N₂O₄
FW 290.31

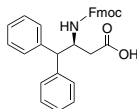


89178-500MG-F

500 mg

(S)-Fmoc-γ,γ-diphenyl-β-Homoala-OH, ≥98.0% (HPLC)

(S)-3-(Fmoc-amino)-4,4-diphenylbutyric acid;
Fmoc-γ,γ-diphenyl-D-β-homoalanine
[332062-08-5]
C₃₁H₂₇NO₄
FW 477.55

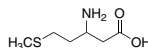


89846-500MG-F

500 mg

D,L-β-Homomethionine, ≥99.0% (TLC)

(±)-3-Amino-5-(methylthio)pentanoic acid
[158570-14-0]
C₆H₁₃NO₂S
FW 163.24

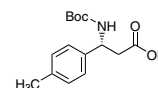


90205-1G-F

1 g

(R)-Boc-4-methyl-β-Phe-OH, ≥98.0% (HPLC)

(R)-3-(Boc-amino)-3-(4-methylphenyl)propionic acid;
Boc-4-methyl-L-β-phenylalanine
[479064-97-6]
C₁₅H₂₁NO₄
FW 279.33

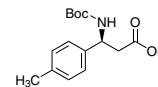


90297-500MG-F

500 mg

(S)-Boc-4-methyl-β-Phe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-3-(4-methylphenyl)propionic acid;
Boc-4-methyl-D-β-phenylalanine
[479064-96-5]
C₁₅H₂₁NO₄
FW 279.33

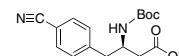


90298-500MG-F

500 mg

(R)-Boc-4-cyano-β-Homophe-OH, ≥98.0% (HPLC)

(R)-3-(Boc-amino)-4-(4-cyanophenyl)butyric acid;
Boc-4-cyano-D-β-homophenylalanine
[269726-86-5]
C₁₆H₂₀N₂O₄
FW 304.34

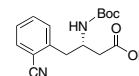


90368-500MG-F

500 mg

(S)-Boc-2-cyano-β-Homophe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-4-(2-cyanophenyl)butyric acid;
Boc-2-cyano-L-β-homophenylalanine
[270065-83-3]
C₁₆H₂₀N₂O₄
FW 304.34

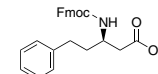


91919-500MG-F

500 mg

(R)-3-(Fmoc-amino)-5-phenyl-pentanoic acid, ≥95.0% (HPLC)

Fmoc-5-phenyl-D-β-norvaline
[269398-87-0]
C₂₆H₂₅NO₄
FW 415.48

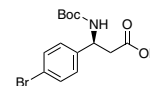


92282-500MG-F

500 mg

(S)-Boc-4-bromo-β-Phe-OH, ≥98.0% (HPLC)

(S)-3-(Boc-amino)-3-(4-bromophenyl)propionic acid;
Boc-4-bromo-D-β-phenylalanine
[261165-06-4]
C₁₄H₁₃BrNO₄
FW 344.20

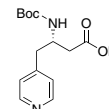


93123-500MG-F

500 mg

(S)-Boc-4-(4-pyridyl)-β-Homoala-OH, ≥97.0% (HPLC)

(S)-3-(Boc-amino)-4-(4-pyridyl)butyric acid;
Boc-4-(4-pyridyl)-L-β-homoalanine
[219297-13-9]
C₁₄H₂₀N₂O₄
FW 280.32

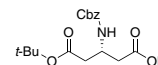


93647-500MG-F

500 mg

Z-β-Glu(OtBu)-OH, ≥94% (HPLC)

Z-L-β-Glutamic acid 5-tert-butyl ester;
Z-D-β-Glutamic acid 5-tert-butyl ester;
Z-L-β-Homoaspartic acid 5-tert-butyl ester
[118247-88-4]
C₁₇H₂₃NO₆
FW 337.37

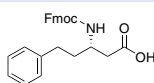


94064-500MG-F

500 mg

(S)-3-(Fmoc-amino)-5-phenyl-pentanoic acid, ≥95.0% (HPLC)

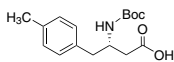
Fmoc-5-phenyl-L-β-norvaline
[219967-74-5]
C₂₆H₂₅NO₄
FW 415.48



94493-500MG-F 500 mg

(S)-Boc-4-methyl-β-Homophe-OH, ≥97.0% (HPLC)

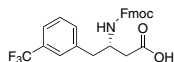
Boc-4-methyl-L-β-homophenylalanine;
(S)-3-(Boc-amino)-4-(4-methylphenyl)butyric acid
[270062-96-9]
C₁₆H₂₃NO₄
FW 293.36



94557-500MG-F 500 mg

(S)-Fmoc-3-(trifluoromethyl)-β-Homophe-OH, ≥97.0% (HPLC)

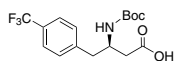
(S)-3-(Fmoc-amino)-4-[3-(trifluoromethyl)phenyl]
butyric acid; Fmoc-3-(trifluoromethyl)-L-β-homo-
phenylalanine
[270065-78-6]
C₂₆H₂₂F₃NO₄
FW 469.45



94591-500MG-F 500 mg

(R)-Boc-4-(trifluoromethyl)-β-Homophe-OH, ≥98.0% (HPLC)

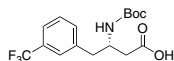
(R)-3-(Boc-amino)-4-[4-(trifluoromethyl)phenyl]
butyric acid; Boc-4-(trifluoromethyl)-D-β-
homophenylalanine
[269726-77-4]
C₁₆H₂₀F₃NO₄
FW 347.33



94592-500MG-F 500 mg

(S)-Boc-3-(trifluoromethyl)-β-Homophe-OH, ≥98.0% (HPLC)

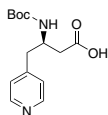
(S)-3-(Boc-amino)-4-[3-(trifluoromethyl)phenyl]
butyric acid; Boc-3-(trifluoromethyl)-L-β-
homophenylalanine
[270065-77-5]
C₁₆H₂₀F₃NO₄
FW 347.33



94603-500MG-F 500 mg

(R)-Boc-4-(4-pyridyl)-β-Homoala-OH, ≥98.0% (HPLC)

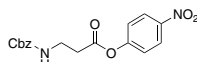
(R)-3-(Boc-amino)-4-(4-pyridyl)butyric acid;
Boc-4-(4-pyridyl)-D-β-homoalanine
[269396-68-1]
C₁₄H₂₀N₂O₄
FW 280.32



95336-500MG-F 500 mg

Z-β-Ala-ONp, ≥98.0% (HPLC)

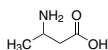
Z-β-alanine 4-nitrophenyl ester
[3642-91-9]
C₁₇H₁₆N₂O₆
FW 344.32



95875-1G 1 g

3-Aminobutanoic acid, 97%

(±)-3-Aminobutyric acid; DL-3-Aminobutyric acid
[541-48-0]
C₄H₉NO₂
FW 103.12

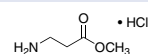


A44207-5G 5 g

A44207-25G 25 g

β-Alanine methyl ester hydrochloride, ≥98.0% (AT)

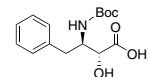
Methyl 3-aminopropionate hydrochloride
[3196-73-4]
C₄H₉NO₂ · HCl
FW 139.58



05210-10G 10 g

(2R,3R)-3-(Boc-amino)-2-hydroxy-4-phenylbutyric acid, ≥98.0% (TLC)

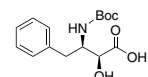
Boc-(2R,3R)-3-amino-2-hydroxy-4-phenylbutyric acid
[77171-41-6]
C₁₅H₂₁NO₅
FW 295.33



15058-25MG 25 mg

(2S,3R)-3-(Boc-amino)-2-hydroxy-4-phenylbutyric acid, ≥98.0% (TLC)

Boc-(2S,3R)-3-amino-2-hydroxy-4-phenylbutyric acid
[62023-65-8]
C₁₅H₂₁NO₅
FW 295.33

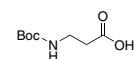


15059-25MG 25 mg

15059-100MG 100 mg

Boc-β-Ala-OH, ≥99.0% (TLC)

Boc-β-alanine
[3303-84-2]
C₈H₁₅NO₄
FW 189.21



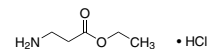
15382-5G 5 g

15382-25G 25 g

15382-100G 100 g

β-Alanine ethyl ester hydrochloride, 98%

Ethyl 3-aminopropionate hydrochloride
[4244-84-2]
C₅H₁₁NO₂ · HCl
FW 153.61

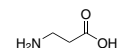


306142-10G 10 g

306142-50G 50 g

β-Alanine

β-Ala; 3-Aminopropionic acid
[107-95-9]
C₃H₇NO₂
FW 89.09

**► 99%**

146064-25G 25 g

146064-500G 500 g

146064-2KG 2 kg

► ≥99.0% (NT)

05160-50G 50 g

05160-250G 250 g

05160-1KG 1 kg

Relax and Enjoy Your Research



Tired of browsing through various catalogs and databases for the research products you need? Relax!

Sigma-Aldrich® introduces Chem Product Central, featuring the industry-leading range of 40,000+ products including cutting-edge catalysts and ligands, novel and classical synthetic reagents, and diverse libraries of building blocks. Chem Product Central uniquely categorized all chemical products by compound classes and application areas—making it easy to quickly find what you need. Regular updates ensure that you have the most innovative products for your research.

Check out Chem Product Central at sigma-aldrich.com/gochem

- 800+ Boronic Acids and Derivatives
 - 200+ Pd Catalysts
 - 400+ Privileged Ligands
 - 5,000+ Heterocyclic Building Blocks
- and so much more!*

Natural Amino Acid Building Blocks for Peptide Synthesis

With a growing peptide drug market the fast, reliable and uncomplicated synthesis of peptides is of paramount importance.⁸ Today, the most common synthetic approaches to medium and even large peptides follow the Boc/Bzl or Fmoc/tBu solid phase peptide synthesis strategies.⁹ A convenient list of the most popular Boc- and Fmoc-protected amino acid building blocks for these two synthesis strategies can be found in this issue. The Boc-protected amino acids are also frequently used in solution phase peptide synthesis procedures.

References: (8) Marx, V. *C&E News* **2005**, 83, 17. (9) Houben-Weyl, *Methods of Organic Chemistry, "Synthesis of Peptides and Peptidomimetics"*, Thieme **2001**, add. and suppl. vol. to the 4th ed. Vol E 22 a.

BOC-protected Amino Acids

Boc-Ala-OH, ≥99.0% (TLC)

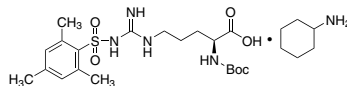
Boc-L-alanine; *N*-(*tert*-Butoxycarbonyl)-L-alanine
[15761-38-3]
C₈H₁₅NO₄
FW 189.21



15380-5G	5 g
15380-25G	25 g
15380-100G	100 g

Boc-Arg(Mts)-OH cyclohexylammonium salt, ≥99.0% (TLC)

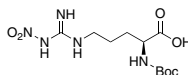
*N*_α-Boc-*N*_ω-(mesitylene-2-sulfonyl)-L-arginine cyclohexylammonium salt
[68262-72-6]
C₂₀H₃₂N₄O₆S · C₆H₁₃N
FW 555.73



15545-2.5G	2.5 g
------------	-------

Boc-Arg(NO₂)-OH, ≥98.5% (T, calc. based on dry substance)

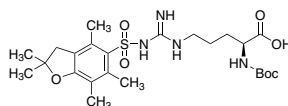
*N*_α-Boc-*N*_ω-nitro-L-arginine
[2188-18-3]
C₁₁H₂₁N₅O₆
FW 319.31



15470-5G	5 g
15470-25G	25 g

Boc-Arg(Pbf)-OH, ≥98.0% (TLC)

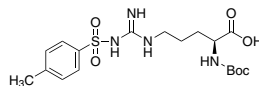
*N*_α-Boc-*N*_ω-(2,2,4,6,7-pentamethyl-dihydrobenzofuran-5-sulfonyl)-L-arginine; *N*_α-Boc-*N*_ω-Pbf-L-arginine
[200124-22-7]
C₂₄H₃₈N₄O₇S
FW 526.65



15038-5G	5 g
----------	-----

Boc-Arg(Tos)-OH

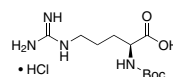
*N*_α-Boc-*N*_ω-tosyl-L-arginine
[13836-37-8]
C₁₈H₂₈N₄O₆S
FW 428.50



15506-5G	5 g
----------	-----

Boc-Arg-OH hydrochloride, ≥98.0% (TLC)

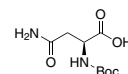
*N*_α-Boc-L-arginine hydrochloride
[35897-34-8]
C₁₁H₂₂N₄O₄ · HCl
FW 310.78



15063-5G	5 g
----------	-----

Boc-Asn-OH, ≥98.5% (T)

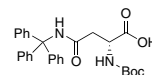
Boc-L-asparagine; *N*_α-(*tert*-Butoxycarbonyl)-L-asparagine
[7536-55-2]
C₉H₁₆N₂O₅
FW 232.23



15381-5G	5 g
15381-25G	25 g

Boc-Asn(Trt)-OH, ≥99.0% (TLC)

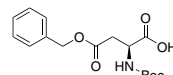
*N*_α-Boc-*N*_γ-trityl-L-asparagine
[132388-68-2]
C₂₈H₃₀N₂O₅
FW 474.55



15562-1G	1 g
----------	-----

Boc-Asp(OBzl)-OH, ≥99.0% (HPLC, sum of enantiomers)

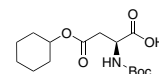
Boc-L-aspartic acid 4-benzyl ester;
N-(*tert*-Butoxycarbonyl)-L-aspartic acid 4-benzyl ester
[7536-58-5]
C₁₆H₂₁NO₆
FW 323.34



15386-5G	5 g
15386-25G	25 g

Boc-Asp(OcHx)-OH, ≥98.0% (T)

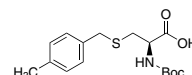
Boc-L-aspartic acid 4-cyclohexyl ester
[73821-95-1]
C₁₅H₂₅NO₆
FW 315.36



15398-5G	5 g
----------	-----

Boc-Cys(Mbz)-OH, ≥96.0% (HPLC)

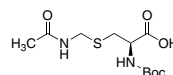
Boc-S-(4-methylbenzyl)-L-cysteine
[61925-77-7]
C₁₆H₂₃NO₄S
FW 325.42



15444-5G	5 g
----------	-----

Boc-Cys(Acm)-OH, ≥96.0% (T)

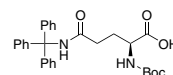
Boc-S-acetamidomethyl-L-cysteine
[19746-37-3]
C₁₁H₂₀N₂O₅S
FW 292.35



15376-5G	5 g
----------	-----

Boc-Gln(Trt)-OH, ≥98.0% (TLC)

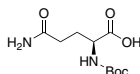
*N*_α-Boc-*N*_δ-trityl-L-glutamine
[132388-69-3]
C₂₉H₃₂N₂O₅
FW 488.57



15563-5G	5 g
----------	-----

Boc-Gln-OH, 98%

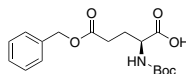
Boc-L-glutamine; *N*_α-(*tert*-Butoxycarbonyl)-L-glutamine
[13726-85-7]
C₁₀H₁₈N₂O₅
FW 246.26



408441-5G	5 g
408441-25G	25 g

Boc-Glu(OBzl)-OH, ≥98.0% (T)

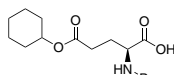
Boc-L-glutamic acid 5-benzyl ester
[13574-13-5]
C₁₇H₂₃NO₆
FW 337.37



15418-5G	5 g
15418-25G	25 g

Boc-Glu(Ochx)-OH, ≥98.0% (TLC, sum of enantiomers)

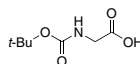
Boc-L-glutamic acid γ-cyclohexyl ester;
Boc-L-glutamic acid 5-cyclohexyl ester
[73821-97-3]
C₁₆H₂₇NO₆
FW 329.39



15437-5G	5 g
----------	-----

Boc-Gly-OH, ≥99.0% (T)

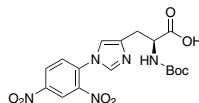
Boc-glycine; *N*-(*tert*-Butoxycarbonyl)glycine
[4530-20-5]
C₇H₁₃NO₄
FW 175.18



15420-5G	5 g
15420-25G	25 g
15420-100G	100 g

Boc-His(Dnp)-OH, ≥98.0% (TLC)

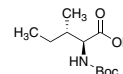
*N*_α-Boc-*N*_ε-(2,4-dinitrophenyl)-L-histidine
[25024-53-7]
C₁₇H₁₉N₅O₈ · C₃H₈O
FW 481.46



15287-5G	5 g
----------	-----

Boc-Ile-OH, 98%

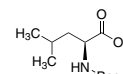
Boc-L-isoleucine; *N*-(*tert*-Butoxycarbonyl)-L-isoleucine
[13139-16-7]
C₁₁H₂₁NO₄
FW 231.29



359653-25G	25 g
------------	------

Boc-Leu-OH, ≥99.0% (HPLC)

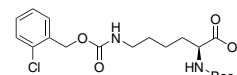
Boc-L-leucine
[13139-15-6]
C₁₁H₂₁NO₄
FW 231.29



15450-5G-F	5 g
15450-25G-F	25 g
15450-100G-F	100 g

Boc-Lys(2-Cl-Z)-OH, ≥98.0% (T)

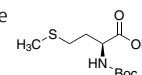
*N*_α-Boc-*N*_ε-(2-chloro-Z)-L-lysine
[54613-99-9]
C₁₉H₂₇ClN₂O₆
FW 414.88



15399-5G	5 g
----------	-----

Boc-Met-OH, 99%

N-(*tert*-Butoxycarbonyl)-L-methionine; Boc-L-methionine
[2488-15-5]
C₁₀H₁₉NO₄S
FW 249.33



408425-5G	5 g
408425-25G	25 g

Got ChemNews?

A Monthly Chemistry E-Newsletter
from Sigma-Aldrich

Visit sigma-aldrich.com/chemnews

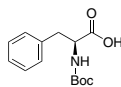


sigma-aldrich.com

SIGMA-ALDRICH®

Boc-Phe-OH, ≥99.0% (T)

Boc-L-phenylalanine; *N*-(*tert*-Butoxycarbonyl)-L-phenylalanine
[13734-34-4]
C₁₄H₁₉NO₄
FW 265.30



15480-5G	5 g
15480-25G	25 g
15480-100G	100 g

Boc-Pro-OH, ≥99.0% (T)

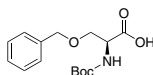
Boc-L-proline; *N*-(*tert*-Butoxycarbonyl)-L-proline
[15761-39-4]
C₁₀H₁₇NO₄
FW 215.25



15490-5G	5 g
15490-25G	25 g
15490-100G	100 g

Boc-Ser(Bzl)-OH, ≥99.0% (T)

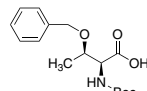
Boc-O-benzyl-L-serine; *N*-(*tert*-Butoxycarbonyl)-O-benzyl-L-serine
[23680-31-1]
C₁₅H₂₁NO₅
FW 295.33



15390-5G	5 g
15390-25G	25 g

Boc-Thr(Bzl)-OH, ≥99.0% (T)

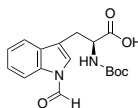
Boc-O-benzyl-L-threonine; *N*-(*tert*-Butoxycarbonyl)-O-benzyl-L-threonine
[15260-10-3]
C₁₆H₂₃NO₅
FW 309.36



15405-5G	5 g
----------	-----

Boc-Trp(For)-OH, ≥97.0% (HPLC)

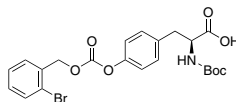
*N*_α-Boc-*N*_(in)-formyl-L-tryptophan
[47355-10-2]
C₁₇H₂₀N₂O₅
FW 332.35



15349-5G-F	5 g
------------	-----

Boc-Tyr(2-Br-Z)-OH, ≥99.0% (TLC, sum of enantiomers)

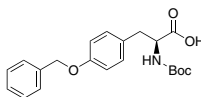
Boc-O-(2-bromo-Z)-L-tyrosine
[47689-67-8]
C₂₂H₂₄BrNO₇
FW 494.33



15431-5G	5 g
15431-25G	25 g

Boc-Tyr(Bzl)-OH, ≥99.0% (T)

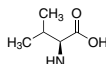
Boc-O-benzyl-L-tyrosine
[2130-96-3]
C₂₁H₂₅NO₅
FW 371.43



15410-5G	5 g
15410-25G	25 g

Boc-Val-OH, ≥99.0% (T)

(*S*)-2-(Boc-amino)-3-methylbutyric acid;
Boc-L-valine; *N*-(*tert*-Butoxycarbonyl)-L-valine
[13734-41-3]
C₁₀H₁₉NO₄
FW 217.26



15528-25G	25 g
15528-100G	100 g

FMOC-protected Amino Acids

Fmoc-Ala-OH, 95%

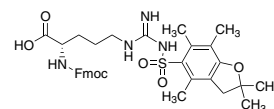
Fmoc-L-alanine; *N*-(9-Fluorenylmethoxycarbonyl)-L-alanine
[35661-39-3]
C₁₈H₁₇NO₄
FW 311.33



531480-5G	5 g
531480-25G	25 g

Fmoc-Arg(Pbf)-OH, ≥98.0% (HPLC)

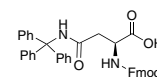
*N*_α-Fmoc-*N*_ω-(2,2,4,6,7-pentamethyl-dihydrobenzofuran-5-sulfonyl)-L-arginine; *N*_α-Fmoc-*N*_ω-Pbf-L-arginine
[154445-77-9]
C₃₄H₄₀N₄O₇S
FW 648.77



47349-1G	1 g
47349-5G	5 g
47349-25G	25 g

Fmoc-Asn(Trt)-OH, ≥97.0% (HPLC, sum of enantiomers)

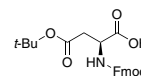
*N*_α-Fmoc-*N*_γ-trityl-L-asparagine; *N*_α-(9-Fluorenylmethoxycarbonyl)-*N*_γ-trityl-L-asparagine
[132388-59-1]
C₃₈H₃₂N₂O₅
FW 596.67



47672-1G-F	1 g
47672-5G-F	5 g
47672-25G-F	25 g

Fmoc-Asp(OtBu)-OH, ≥98.0% (HPLC)

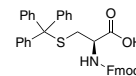
Fmoc-L-aspartic acid 4-*tert*-butyl ester
[71989-14-5]
C₂₃H₂₅NO₆
FW 411.45



47618-1G-F	1 g
47618-10G-F	10 g
47618-25G-F	25 g

Fmoc-Cys(Trt)-OH, ≥95.0% (HPLC, sum of enantiomers)

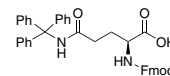
*N*_α-Fmoc-*S*-trityl-L-cysteine; *N*-(9-Fluorenylmethoxycarbonyl)-*S*-trityl-L-cysteine
[103213-32-7]
C₃₇H₃₁NO₄S
FW 585.71



47695-5G	5 g
----------	-----

Fmoc-Gln(Trt)-OH, ≥98.0% (HPLC)

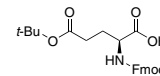
*N*_α-Fmoc-*N*_δ-trityl-L-glutamine
[132327-80-1]
C₃₉H₃₄N₂O₅
FW 610.70



47674-1G-F	1 g
47674-5G-F	5 g
47674-25G-F	25 g

Fmoc-Glu(OtBu)-OH, ≥98.0% (HPLC)

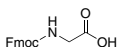
Fmoc-L-glutamic acid 5-*tert*-butyl ester
[71989-18-9]
C₂₄H₂₇NO₆
FW 425.47



47625-1G	1 g
47625-10G	10 g
47625-25G	25 g

Fmoc-Gly-OH, ≥98.0% (T)

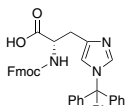
Fmoc-glycine
[29022-11-5]
C₁₇H₁₅NO₄
FW 297.31



47627-5G-F	5 g
47627-50G-F	50 g

Fmoc-His(Trt)-OH, ≥98.0% (HPLC, sum of enantiomers)

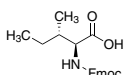
N_α-Fmoc-N_(trt)-trityl-L-histidine
[109425-51-6]
C₄₀H₃₃N₃O₄
FW 619.71



47639-5G-F	5 g
47639-25G-F	25 g

Fmoc-Ile-OH, ≥98.0% (T)

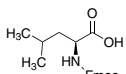
Fmoc-L-isoleucine; N-(9-Fluorenylmethoxycarbonyl)-L-isoleucine
[71989-23-6]
C₂₁H₂₃NO₄
FW 353.41



47628-5G-F	5 g
47628-50G-F	50 g

Fmoc-Leu-OH, ≥97.0% (HPLC)

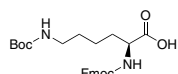
Fmoc-L-leucine; N-(9-Fluorenylmethoxycarbonyl)-L-leucine
[35661-60-0]
C₂₁H₂₃NO₄
FW 353.41



47633-5G-F	5 g
47633-50G-F	50 g

Fmoc-Lys(Boc)-OH, ≥98.0% (HPLC)

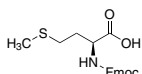
N_α-Fmoc-N_ε-Boc-L-lysine
[71989-26-9]
C₂₆H₃₂N₂O₆
FW 468.54



47624-1G-F	1 g
47624-10G-F	10 g
47624-25G-F	25 g

Fmoc-Met-OH, ≥98.0% (HPLC)

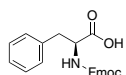
Fmoc-L-methionine
[71989-28-1]
C₂₀H₂₁NO₄S
FW 371.45



47634-5G-F	5 g
------------	-----

Fmoc-Phe-OH, 98%

Fmoc-L-phenylalanine; N-(9-Fluorenylmethoxycarbonyl)-L-phenylalanine
[35661-40-6]
C₂₄H₂₁NO₄
FW 387.43



338338-5G	5 g
338338-25G	25 g

Fmoc-Pro-OH, ≥99.0% (HPLC)

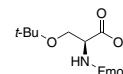
Fmoc-L-proline; N-(9-Fluorenylmethoxycarbonyl)-L-proline
[71989-31-6]
C₂₀H₁₉NO₄
FW 337.37



47636-5G-F	5 g
47636-50G-F	50 g

Fmoc-Ser(tBu)-OH, ≥98.0% (HPLC)

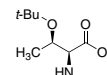
Fmoc-O-tert-butyl-L-serine
[71989-33-8]
C₂₂H₂₅NO₅
FW 383.44



47619-1G-F	1 g
47619-10G-F	10 g
47619-25G-F	25 g

Fmoc-Thr(tBu)-OH, ≥98.0% (HPLC)

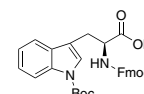
Fmoc-O-tert-butyl-L-threonine
[71989-35-0]
C₂₃H₂₇NO₅
FW 397.46



47622-1G-F	1 g
47622-10G-F	10 g

Fmoc-Trp(Boc)-OH, ≥97.0% (HPLC)

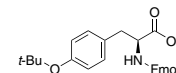
N_α-Fmoc-N_(trt)-Boc-L-tryptophan;
N_(trt)-Boc-N_α-Fmoc-L-tryptophan
[143824-78-6]
C₃₁H₃₀N₂O₆
FW 526.58



47561-1G-F	1 g
47561-5G-F	5 g
47561-25G-F	25 g

Fmoc-Tyr(tBu)-OH, ≥98.0% (HPLC)

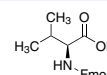
Fmoc-O-tert-butyl-L-tyrosine; N-(9-Fluorenylmethoxycarbonyl)-O-tert-butyl-L-tyrosine
[71989-38-3]
C₂₈H₂₉NO₅
FW 459.53



47623-1G-F	1 g
47623-10G-F	10 g
47623-25G-F	25 g

Fmoc-Val-OH, ≥98.0% (HPLC)

Fmoc-L-valine; N-(9-Fluorenylmethoxycarbonyl)-L-valine
[68858-20-8]
C₂₀H₂₁NO₄
FW 339.39



47638-50G-F	50 g
-------------	------

New Tools for PEGylation

PEG polymers offer numerous favorable characteristics including high water solubility, high mobility in solution, lack of toxicity and immunogenicity, and ready clearance from the body. Thus, the chemical modification of biologically active compounds, such as peptides, antibody fragments, enzymes, or small molecules with polyethylene glycol chains (referred to as "PEGylation"), is an effective method to tailor molecular properties to particular applications (**Figure 1**).

Drugs can especially benefit from PEGylation as has been proved by a continuously growing number of PEG conjugated drugs on the market. By increasing the molecular mass of proteins and peptides and shielding them from proteolytic enzymes, PEGylation improves pharmacokinetics.¹⁰ The judicious choice of PEG size and linkers enhance a potential drug's solubility and its distribution within an organism.¹¹

Numerous review articles concerning the application of PEG polymers in drug delivery are available.¹² In some recent reports PEGylation is discussed as a tool to facilitate the drug delivery of nanocarrier systems and microparticles.^{13,14}

In this issue of ChemFiles, Sigma-Aldrich® is proud to further expand its offer of tools for PEGylation with new bifunctional, small and medium-sized ethylene glycol building blocks. For a complete list of PEGylation products, please visit our online product catalog at sigma-aldrich.com/chemprod and browse through the peptide synthesis category.

References: (10) Harris, J.M.; Chess, R.B. *Nature Rev. Drug. Disc.* **2003**, 2, 214. (11) Fee, C. *Biotechn. and Bioeng.* **2007**, 98, 725. (12)(a) "Peptide and Protein PEGylation, a Review of Problems and Solutions", Veronese, F.M. *Biomaterials* **2001**, 5, 405. (b) "Chemistry for Peptide and Protein PEGylation", Roberts, M.J.; Bentley, M.D.; Jarris, J.M. *Adv. Drug Del. Rev.* **2002**, 4, 459. (c) "PEGylation, Successful Approach to Drug Delivery", Veronese, F.M.; Pasut, G. *Drug Disc. Today* **2005**, 10, 1451. (d) "Advances in PEGylation of Important Biotech Molecules: Delivery Aspects", Ryan, S.M.; Mantovani, G.; Wang, X.; Haddleton, D.M.; Brayden, D.J. *Exp. Op. on Drug Del.* **2008**, 5, 371. (13) Howard, M.D.; Jay, M.; Dziubla, T.D.; Lu, X. *Biomed. Nanotechn.* **2008**, 4, 133. (14) Wattendorf, U.; Merkle, H.P. *J. Pharm. Sci.* **2008**, 97, 4655.

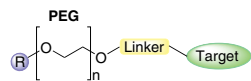
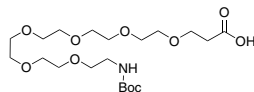


Figure 1

New PEG derivatives

21-(Boc-amino)-4,7,10,13,16,19-hexaoxaheneicosanoic acid, ≥97.0% (H-NMR)

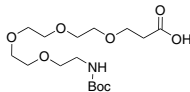
C₂₀H₃₉NO₁₀
FW 453.52



38263-100MG 100 mg
38263-500MG 500 mg

15-(Boc-amino)-4,7,10,13-tetraoxapentadecanoic acid, ≥97.0% (H-NMR)

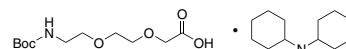
C₁₆H₃₁NO₈
FW 365.42



30953-100MG 100 mg
30953-500MG 500 mg

2-[2-(Boc-amino)ethoxy]ethoxyacetic acid (dicyclohexylammonium) salt, ≥98.0% (TLC)

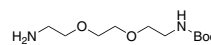
Boc-8-amino-3,6-dioxaoctanoic acid (dicyclohexylammonium) salt
C₁₁H₂₁NO₆ · C₁₂H₂₃N
FW 444.61



14766-500MG-F 500 mg

N-Boc-2,2'-(ethylenedioxy)diethylamine, ≥95.0% (NT)

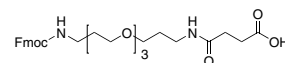
N-Boc-3,6-dioxo-1,8-octanediamine; *tert*-Butyl 2-[2-(2-aminoethoxy)ethoxy]ethylcarbamate
[153086-78-3]
C₁₁H₂₄N₂O₄
FW 248.32



89761-1G-F 1 g
89761-5G-F 5 g

N-Fmoc-N'-succinyl-4,7,10-trioxa-1,13-tridecanediamine, 95% (HPLC)

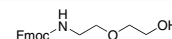
17-Oxo-6,9,12-trioxa-2,16-diazaeicosanedioic acid 1-(9H-fluoren-9-ylmethyl)ester
[172089-14-4]
C₂₉H₃₈N₂O₈
FW 542.62



671517-1G 1 g
671517-5G 5 g

2-[2-(Fmoc-amino)ethoxy]ethanol, 99%

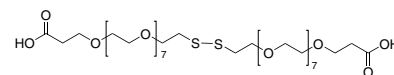
[2-(2-Hydroxy-ethoxy)ethyl]carbamic acid 9H-fluoren-9-ylmethyl ester
[299430-87-8]
C₁₉H₂₁NO₄
FW 327.37
≥98.5% (HPLC)



670529-500MG 500 mg

4,7,10,13,16,19,22,25,32,35,38,41,44,47,50,53-Hexadeca-28,29-dithiahexapentacontanedioic acid

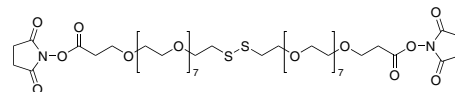
PEG acid disulfide (n=7)
[873013-93-5]
C₃₈H₇₄O₂₀S₂
FW 915.11



671525-100MG 100 mg
671525-500MG 500 mg

4,7,10,13,16,19,22,25,32,35,38,41,44,47,50,53-Hexadeca-28,29-dithiahexapentacontanedioic acid di-N-succinimidyl ester

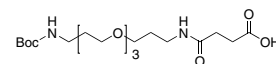
PEG NHS ester disulfide (n=7)
[947601-98-1]
C₄₆H₈₀N₂O₂₄S₂
FW 1109.26



671630-100MG 100 mg
671630-500MG 500 mg

N-Boc-N'-succinyl-4,7,10-trioxa-1,13-tridecanediamine, 95% (HPLC)

17-Oxo-6,9,12-trioxa-2,16-diazaeicosanedioic acid 1-(1,1-dimethylethyl)ester
[250612-31-8]
C₁₉H₃₆N₂O₈
FW 420.50

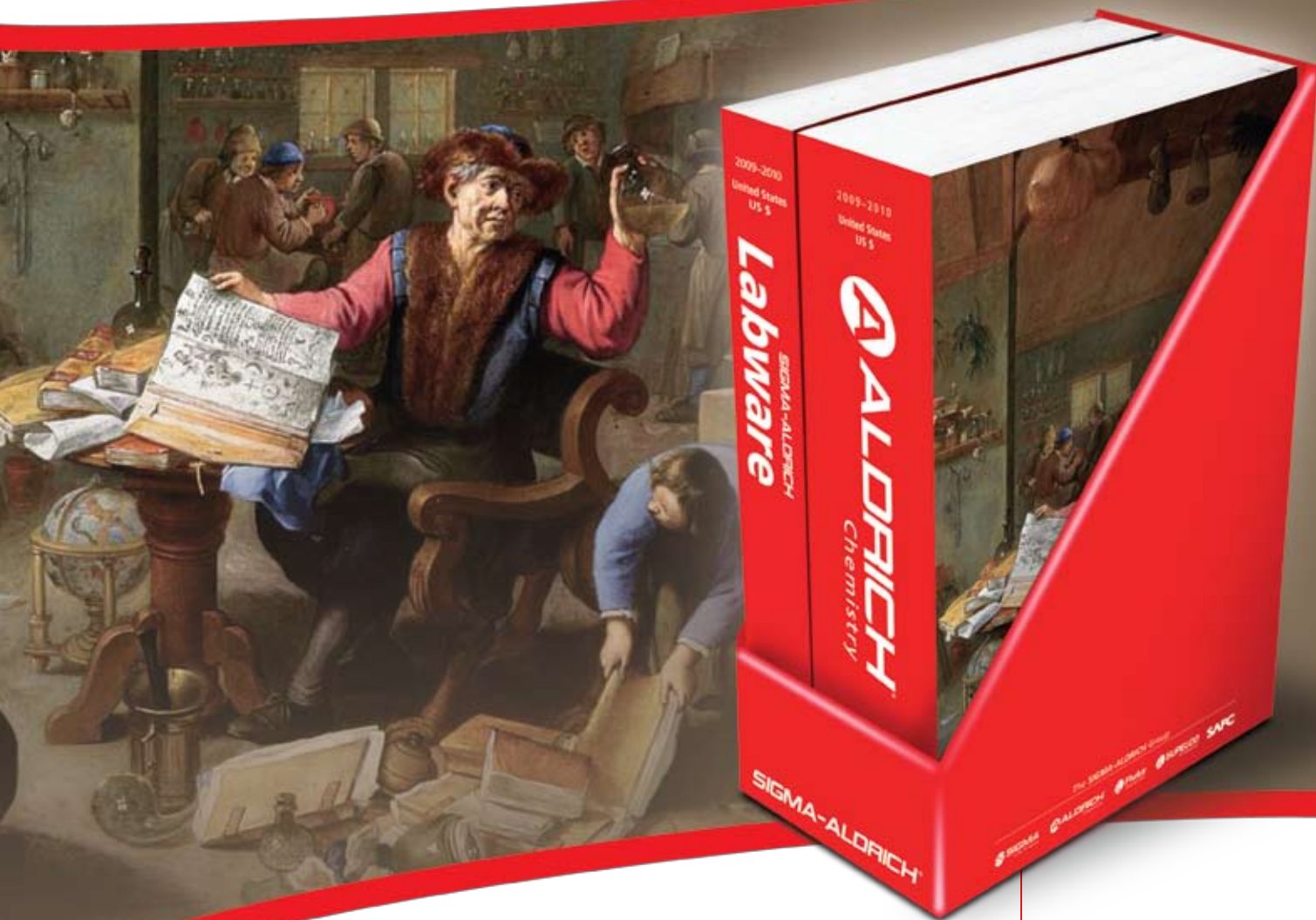


671401-1G 1 g
671401-5G 5 g

New Aldrich® Handbook!



Over 6,000 Innovative New Products



Reserve your copy of the new 2009-2010
Aldrich Handbook set today.

2009-2010 Aldrich Handbook

- 10,000 chemical structures
- 8,500 updated literature citations
- Extensive chemical & physical data

Labware Catalog

- Innovative new products
- Improved product images
- New technical information index

To request your
complimentary Aldrich
Handbook set, visit
sigma-aldrich.com/aldrichcat

Argentina

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

Australia

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

Austria

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

Belgium

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

Brazil

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

Canada

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

China

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

Czech Republic

SIGMA-ALDRICH spol. s r. o.
Tel: (+420) 246 003 200
Fax: (+420) 246 003 291

Denmark

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

Finland

SIGMA-ALDRICH FINLAND OY
Tel: (+358) 9 350 9250
Fax: (+358) 9 350 92555

France

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

Germany

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

Greece

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

Hungary

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

India

SIGMA-ALDRICH CHEMICALS
PRIVATE LIMITED
Telephone
Bangalore: (+91) 80 6621 9600
New Delhi: (+91) 11 4358 8000
Mumbai: (+91) 22 2570 2364
Hyderabad: (+91) 40 4015 5488
Fax
Bangalore: (+91) 80 6621 9650
New Delhi: (+91) 11 4358 8001
Mumbai: (+91) 22 2579 7589
Hyderabad: (+91) 40 4015 5466

Ireland

SIGMA-ALDRICH IRELAND LTD.
Free Tel: 1800 200 888
Free Fax: 1800 600 222
Tel: +353 (0) 402 20370
Fax: + 353 (0) 402 20375

Israel

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

Italy

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

Japan

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

Korea

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

Malaysia

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

Mexico

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

The Netherlands

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

New Zealand

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

Norway

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

Poland

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

Portugal

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

Russia

SIGMA-ALDRICH RUS, LLC
Tel: +7 (495) 621 6037
+7 (495) 621 5828
Fax: +7 (495) 621 5923

Singapore

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

Slovakia

SIGMA-ALDRICH spol. s r. o.
Tel: (+421) 255 571 562
Fax: (+421) 255 571 564

South Africa

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

Spain

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

Sweden

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

Switzerland

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

United Kingdom

SIGMA-ALDRICH COMPANY LTD.
Free Tel: 0800 717 181
Free Fax: 0800 378 785
Tel: (+44) 1747 833 000
Fax: (+44) 1747 833 313
SAFC (UK) Tel: 01202 712305

United States

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

Internet

sigma-aldrich.com

World Headquarters

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

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

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

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

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