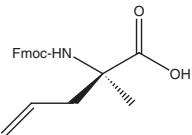
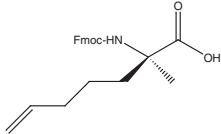
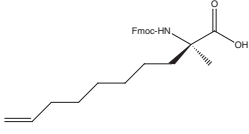
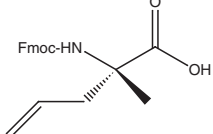
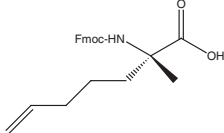
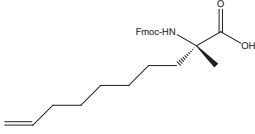
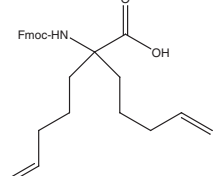


Stapled Peptide & Click Chemistry

AnaSpec, Inc is a leading provider of Combinatorial Chemistry reagents, and carries the industry's largest collection of unusual amino acids, ISO 9001

- Premier Experts in Innovative Amino Acids
- Choose from 1250 other amino acid products
- Custom synthesis available

Stapled Peptide building blocks

Product	Structure	Size	Catalog #
(R)-N-Fmoc-2-(2'-propenyl) alanine		100 mg 500 mg 1 g	64171-100 64171-500 64171-1000
(R)-N-Fmoc-2-(4'-pentenyl) alanine		100 mg 500 mg 1 g	64173-100 64173-500 64173-1000
(R)-N-Fmoc-2-(7'-octenyl) alanine		100 mg 500 mg 1 g	64175-100 64175-500 64175-1000
(S)-N-Fmoc-2-(2'-propenyl) alanine		100 mg 500 mg 1 g	64170-100 64170-500 64170-1000
(S)-N-Fmoc-2-(4'-pentenyl) alanine		100 mg 500 mg 1 g	64172-100 64172-500 64172-1000
(S)-N-Fmoc-2-(7'-octenyl) alanine		100 mg 500 mg 1 g	64174-100 64174-500 64174-1000
Fmoc-2, 2-Bis(4'-pentenyl) glycine (B5)		100 mg 500 mg 1 g	64177-100 64177-500 64177-1000



ISO 9001 Certified

Custom Synthesis & Bulk Pricing Available

©AnaSpec, Inc. | 34801 Campus Dr. | Fremont, CA 94555 | 510.791.9560 | service@anaspec.com | www.anaspec.com

Click Chemistry building blocks

Product	Structure	Size	Catalog #
2-Amino-3 mercapto-N-(prop-2-ynyl)propionamide		250 mg 1 g	64168-025 64168-1
2-Amino-N-(3-azidopropyl)-3-mercaptopropionamide		250 mg 1 g	64169-025 64169-1
Azidohomoalanine		100 mg	63669
Fmoc-Azidohomoalanine		100 mg	63670
Boc-Azidohomoalanine		100 mg	64167
R=Fmoc or Boc			
Fmoc-4-Azidophenylalanine		250 mg 1 g	64347-250 64347-1000
Fmoc-Lys(N ₃)-OH		250 mg 1 g	53100-F025 53100-F1
Boc-Lys(N ₃)-OH		250 mg 1 g	53100-B025 53100-B1
R=Fmoc or Boc			
Fmoc-L-propargylglycine		1 g 5 g	26264-F1 26264-F5
Boc-L-propargylglycine		1 g 5 g	26264-B1 26264-B5
R=Fmoc or Boc			
Fmoc-D-propargylglycine		1g 5 g	26265-F1 26265-F5
Boc-D-Propargylglycine		1g 5g	26265-B1 26265-B5
R=Fmoc or Boc			
TBTA, Tris[(1-benzyl-1H-1,2,3-triazol-4-yl)methyl]amine		50 mg 500 mg	63360-50 63360-500



ISO 9001 Certified

Custom Synthesis & Bulk Pricing Available

©AnaSpec, Inc. | 34801 Campus Dr. | Fremont, CA 94555 | 510.791.9560 | service@anaspec.com | www.anaspec.com