

GO™ Peptides

β-Amyloid Peptides

FRET Peptides

Phosphopeptides

Dye-labeled Peptides

Enzyme Substrates & Inhibitors

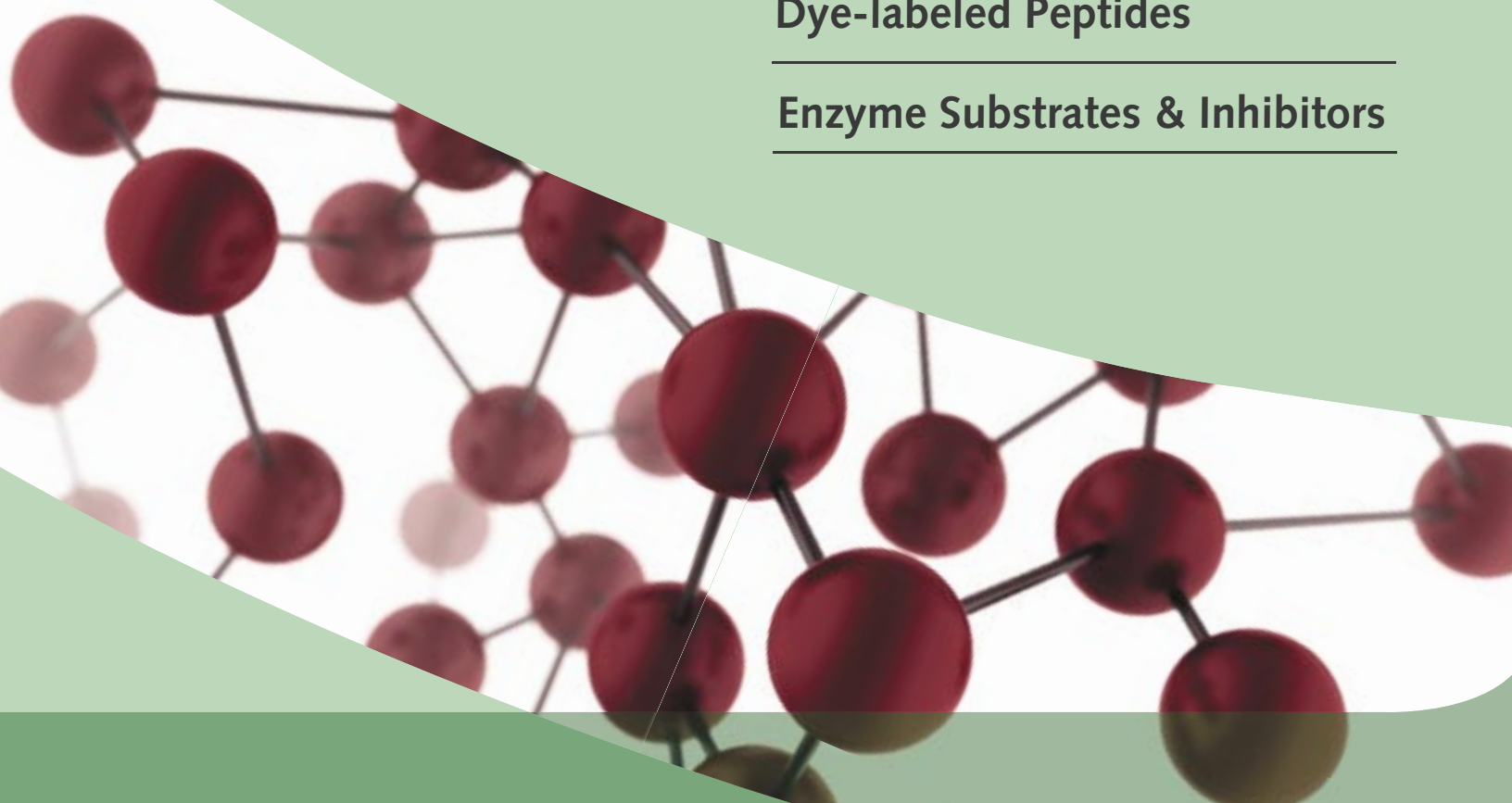


Table of Contents

A	Amylin Peptides	1	Influenza Viral Peptides.....	24	
	Amyloid Peptides	1	K	Kinases / Phosphatase Substrates	24
	Angiotensins and Related Peptides	8	L	LL-37 / CRAMP Peptides	27
	Antimicrobial and Related Peptides	9		Lymphocytic Choremeningitis Virus (LCMV) Peptides	28
	Apelin Peptides	10	M	MAGE Peptides.....	28
	Apoptosis Peptides	10		Matrix Metalloproteinase Substrates	
	Aquaporin Peptides.....	11		and Related Peptides.....	28
	Arg-Phe-Amide (RFamide) Related Peptides	11		Melan-A and Mucin Related Peptides	29
	Autoimmunity Related Peptides	11		Melanocyte Stimulating Hormones and Analogs.....	29
B	Bacterial Peptides	11		MHC Related Peptides.....	29
	Bcl / BAD Peptides.....	11		Microbial Peptides.....	30
	Bradykinins.....	13		Myelin Oligodendrocyte Glycoprotein (MOG) /	
C	Cancer Research Peptides	13		Myelin Basic Protein (MBP) Fragments	30
	Cardiac Peptides.....	15		Myosin Peptides.....	31
	CEF Control Peptide Pools	15	N	Neuropeptides	31
	Cell Adhesion Peptides	15		NF-kB/Transcription Factors Related Peptides	33
	Cell Permeable / Drug Delivery Peptides.....	15		Nuclear Export Signal Peptides	33
	Chaperone Peptides	16	O	Opioid Related Peptides	33
	Chemotaxis Peptides	16		OVA Peptides	33
	ClearPoint™-Heavy Isotope Labeled Peptides.....	17		Oxytocins, Vasopressin and Related Peptides	34
	Cytochromes and Related Peptides.....	17	P	p53 Peptides	34
D	Dengue Viral Peptides.....	17		Pancreatic Polypeptides.....	34
	Dynorphins	17		Parathyroid Hormone Related Peptides.....	34
E	Endorphins	17		Peptide Standards.....	34
	Endothelins	17		Phytochelatin Peptides.....	34
	Enkephalins.....	18		Prion Protein (PrP) Fragments	35
	Enzyme Substrates and Inhibitors	18		Protease Inhibitors	35
	Exendins	18		Protein Structure / Folding Related Peptides	35
F	Fibrinogen and Related Peptides.....	19		Protein Phosphorylation Related Peptides.....	35
	Fibronectin Fragments	19		Proteolipid Protein (PLP) Fragments	36
	FRET Peptides	19	S	Salusin Peptides	36
G	Galanins	20		Signal Transduction Peptides.....	36
	Gastrins.....	20		Somatostatins	37
	Ghrelin and Obestatins.....	21		Steroid Receptor Coactivators (SRCs)	38
	Glucagon-Like Peptides	21		Substance P and Analogs	38
	Growth Factor Fragments	21	T	Tachykinin and Related Peptides	38
	Guanylin	21		Tag Peptides	38
H	Heat Shock Protein Fragments.....	21		Thrombin Related Peptides.....	38
	Hepatitis C Virus (HCV) Related Peptides	22		TNF Peptides	39
	Histone Peptides	22		Toxins	39
	HIV Related Peptides.....	23		Tyrosinase Peptides.....	40
	Hydrocarbon Stapled Peptides	23	V	Viral Peptides	40
I	Ion Channel Peptides.....	23			

Please see our website for a complete listing of our peptides
www.anaspec.com

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
Amylin Peptides				
62763	Amylin (20-29), Islet Amyloid Polypeptide, IAPP (20-29), cat	SNNFGAILS	1019.1	1 mg
62764	Amylin (20-29), Islet Amyloid Polypeptide, IAPP (20-29), hamster	NNNLGPVLSP	1024.2	1 mg
62765	Amylin (20-29), Islet Amyloid Polypeptide, IAPP (20-29), rat	SNNLGPVLPP	1007.2	1 mg
62900	hIAPP, Human Islet Amyloid Polypeptide (8-20)	ATQLRANFLVHSS	1443.6	1 mg
62898	hIAPP, Human Islet Amyloid Polypeptide (22-27)	NFGAIL	633.8	1 mg
61937	[NMeG ²⁴ , NMeI ²⁶]-hIAPP, Human Islet Amyloid Polypeptide (22-27)	NF-(NMe-G)-A-(NMe-I)-L	661.8	1 mg
62899	hIAPP, Human Islet Amyloid Polypeptide (23-27)	FGAIL	519.7	1 mg
Amyloid Peptides				
61929	Activity-Dependent Neurotrophic Factor (6-14), ADNF (6-14)	SALLRSIPA	927.1	1 mg
62132	β-Amyloid (1-8)-Cys	DAEFRHDS	1079.1	1 mg
61970	β-Amyloid (1-9)	DAEFRHDSG	1033.0	1 mg
62131	β-Amyloid (1-9)-Ala-NH ₂	DAEFRHDSGA-NH ₂	1103.1	1 mg
62138	β-Amyloid (1-9)-Gly-Gly	DAEFRHDSGGG	1147.1	1 mg
63311	β-Amyloid (1-9), HiLyte Fluor™ 555-labeled	HiLyte Fluor™ 555-DAEFRHDSG	1501.6	0.1 mg
61973	β-Amyloid (1-9)-Lys(Biotin-LC)-NH ₂	DAEFRHDSG-K(Biotin-LC)-NH ₂	1499.7	1 mg
62144	β-Amyloid (1-10)-Cys	DAEFRHDSGYC	1299.4	1 mg
62135	β-Amyloid (1-10)-Lys(Biotin)	DAEFRHDSGY-K(Biotin)	1550.7	1 mg
62137	β-Amyloid (1-10)-Lys-Lys	DAEFRHDSGYKK	1452.6	1 mg
62134	β-Amyloid (1-10)-Gly-Gly-Lys(Biotin)	DAEFRHDSGYGG-K(Biotin)	1664.8	1 mg
62969	β-Amyloid (1-10), mouse, rat	DAEFGHDSGF	1081.1	1 mg
63331	β-Amyloid (1-11), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-DAEFRHDSGYE	1681.7	0.1 mg
63334	β-Amyloid (1-11), HiLyte Fluor™ 555-labeled	HiLyte Fluor™ 555-DAEFRHDSGYE	1793.9	0.1 mg
62967	β-Amyloid (1-12)	DAEFRHDSGYEV	1424.5	1 mg
63333	β-Amyloid (1-12)-Cys-NH ₂	DAEFRHDSGYEVC-NH ₂	1527.6	1 mg
62973	β-Amyloid (1-12)-Lys(Biotin)-NH ₂	DAEFRHDSGYEV-K(Biotin)-NH ₂	1778.0	1 mg
62978	β-Amyloid (1-13), mouse, rat	DAEFGHDSGFV	1465.5	1 mg
61798	β-Amyloid (1-15)	DAEFRHDSGYEVHHQ	1826.9	1 mg
62002-05	β-Amyloid (1-15)-Lys(HiLyte Fluor™ 488)	DAEFRHDSGYEVHHQ-K(HiLyte Fluor™ 488)	2311.4	0.5 mg
62461	Biotin-β-Amyloid (1-15)	Biotin-DAEFRHDSGYEVHHQ	2053.2	1 mg
62970	Biotin-LC-β-Amyloid (1-15)	Biotin-LC-DAEFRHDSGYEVHHQ	2166.4	1 mg
62474	β-Amyloid (1-16)-8 branched MAPS	DAEFRHDSGYEVHHQK-on 8 branched MAPS		1 mg
63332	β-Amyloid (1-16), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-DAEFRHDSGYEVHHQK	2311.4	0.1 mg
63330	β-Amyloid (1-16), HiLyte Fluor™ 555-labeled	HiLyte Fluor™ 555-DAEFRHDSGYEVHHQK	2423.6	0.1 mg
62136	β-Amyloid (1-16)-Lys(Biotin-LC)-NH ₂	DAEFRHDSGYEVHHQK-K(Biotin-LC)-NH ₂	2421.7	1 mg
62457	Biotin-β-Amyloid (1-16)	Biotin-DAEFRHDSGYEVHHQK	2181.3	1 mg
62458	Biotin-LC-β-Amyloid (1-16)	Biotin-LC-DAEFRHDSGYEVHHQK	2294.5	1 mg
62464	β-Amyloid (1-16)-Lys(Biotin-LC)-NH ₂	DAEFRHDSGYEVHHQK-K(Biotin-LC)-NH ₂	2421.7	1 mg
61955	β-Amyloid (1-17)	DAEFRHDSGYEVHHQKL	2068.2	1 mg
62473	β-Amyloid (1-17)-Cys	DAEFRHDSGYEVHHQKLC	2171.3	1 mg
62952	β-Amyloid (1-17), FAM-labeled	5-FAM-DAEFRHDSGYEVHHQKL	2426.5	0.1 mg
62953	β-Amyloid (1-17), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-DAEFRHDSGYEVHHQKL	2424.6	0.1 mg
62977	β-Amyloid (1-17), mouse, rat	DAEFGHDSGFVHHQKL	1972.1	1 mg
62456	β-Amyloid (1-20)	DAEFRHDSGYEVHHQKLVFF	2461.7	1 mg
62966	β-Amyloid (1-22)	DAEFRHDSGYEVHHQKLVFFAE	2661.9	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62141	β-Amyloid (1-23)-Gly	DAEFRHDSGYEVHHQKLVFFAEDG	2834.0	1 mg
62475	β-Amyloid (1-24)-Cys	DAEFRHDSGYEVHHQKLVFFAEDVC	2979.3	1 mg
62465	β-Amyloid (1-28)-Lys(Biotin)-NH ₂	DAEFRHDSGYEVHHQKLVFFAEDVGSNK-K(Biotin)-NH ₂	3616.0	0.5 mg
62995	β-Amyloid (1-28), mouse, rat	DAEFGHDSGFVEVRHQKLVFFAEDVGSNK	3166.4	1 mg
62890	[Ala ^{16,17,20}]-β-Amyloid (1-28)	DAEFRHDSGYEVHHQAAVFAEDVGSNK	3087.2	1 mg
62441	Bromoacetyl-β-Amyloid (1-28)	Bromoacetyl-DAEFRHDSGYEVHHQKLVFFAEDVGSNK	3383.5	1 mg
62140	β-Amyloid (1-33)-Ala	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAI-IGA	3745.1	1 mg
61799	β-Amyloid (1-34)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGL	3787.2	1 mg
62998	β-Amyloid (1-34), mouse, rat	DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGL	3691.1	1 mg
63748	[Leu ³²]-β-Amyloid(1-36)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAILGLMV	4017.5	0.5 mg
63329	β-Amyloid (1-37)-Lys(Biotin), FAM-labeled	5-FAM-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVG-K(Biotin)	4787.3	0.1 mg
62476	β-Amyloid (1-38), mouse, rat	DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGLMVGG	4035.5	1 mg
63000	β-Amyloid (1-39), mouse, rat	DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGLMVGGV	4134.7	0.5 mg
63318	β-Amyloid (1-40), 5-ROX-labeled	5-ROX-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	4683.5	0.1 mg
62426	β-Amyloid (1-40) Binding Peptide	DWGKGGWRWRLWPGASGKTEA	2215.5	1 mg
62427	β-Amyloid (1-40) Binding Peptide, Biotin-labeled	Biotin-DWGKGGWRWRLWPGASGKTEA	2441.8	1 mg
62943	β-Amyloid (1-40) Binding Peptide, FAM-labeled	5-FAM-DWGKGGWRWRLWPGASGKTEA	2573.8	0.1 mg
62233	β-Amyloid (1-40)-Cys	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVC	4433	0.1 mg
61949-01	β-Amyloid (1-40), DEAC-labeled	DEAC-DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV, DEAC=(7-Diethylaminocoumarin-3-yl)carboxylic acid	4573.1	0.1 mg
61483-01	β-Amyloid (1-40)-Lys(Biotin)-NH ₂	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂	4684.4	0.1 mg
61483-05	β-Amyloid (1-40)-Lys(Biotin)-NH ₂	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂	4684.4	0.5 mg
63356	β-Amyloid (1-40)-Lys(Biotin)-NH ₂ , mouse, rat	DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂	4587.3	0.1 mg
61968-05	[Ala ²⁸]-β-Amyloid (1-40)	DAEFRHDSGYEVHHQALVFFAEDVGSNKGAIIGLMVGGVV	4215.7	0.5 mg
63320	[Asn ⁷]-β-Amyloid (1-40), Tottori-Japanese Mutation	DAEFRHNSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	4328.9	0.5 mg
62145	[Asn ²³]-β-Amyloid (1-40), Iowa Mutation	DAEFRHDSGYEVHHQKLVFFAENVGSNKGAIIGLMVGGVV	4328.9	0.5 mg
61717-01	Biotin-LC-β-Amyloid (1-40), mouse, rat	Biotin-LC-DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGLMVGGVV	4573.3	0.1 mg
61717-05	Biotin-LC-β-Amyloid (1-40), mouse, rat	Biotin-LC-DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGLMVGGVV	4573.3	0.5 mg
61717-1	Biotin-LC-β-Amyloid (1-40), mouse, rat	Biotin-LC-DAEFGHDSGFVEVRHQKLVFFAEDVGSNKGAIIGLMVGGVV	4573.3	1 mg
63321	[Arg ⁶]-β-Amyloid (1-40), English Mutation	DAEFRDSDGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	4348.9	0.5 mg
62815-1	[Cys ⁷]-β-Amyloid (1-40)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	4317.9	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62815-05	[Cys ⁷]-β-Amyloid (1-40)	DAEFRHCSGYEVHHQKLVFFAEDVGSNKGAI-IGLMVGGVV	4317.9	0.5 mg
63895	[Cys ²⁰]-β-Amyloid (1-40)	DAEFRHDSGYEVHHQKLVFCAEDVG-SNKGAIIGLMVGGVV	4285.8	0.5 mg
61962-01	β-Amyloid (1-40)-Lys(LC-biotin)-NH ₂ , FAM-labeled	5-FAM-DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVV-K(LC-Biotin)-NH ₂	5154.9	0.1 mg
61948-05	[Gln ⁹]-β-Amyloid (1-40)	DAEFRHDSQYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVV	4401.0	0.5 mg
62150	[Gly ²¹]-β-Amyloid (1-40), A21G, Flemish Mutation	DAEFRHDSGYEVHHQKLVFFGEDVG-SNKGAIIGLMVGGVV	4315.9	0.5 mg
61261	[Gln ²²]-β-Amyloid (1-40), Dutch Mutation	DAEFRHDSGYEVHHQKLVFFAQDVGSNKGAI-IGLMVGGVV	4328.9	0.5 mg
62146	[Gln ²² , Asn ²³]-β-Amyloid (1-40), E22Q/D23N Dutch/Iowa double mutation	DAEFRHDSGYEVHHQKLVFFAQNVGSNKGAI-IGLMVGGVV	4327.9	0.5 mg
62830-01	[Leu ³³]-β-Amyloid (1-40)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAI-ILLMVGCVV	4386	0.1 mg
62830-05	[Leu ³³]-β-Amyloid (1-40)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAI-ILLMVGCVV	4386	0.5 mg
62147	[Lys ²²]-β-Amyloid (1-40), Italian Mutation	DAEFRHDSGYEVHHQKLVFFAKDVG-SNKGAIIGLMVGGVV	4328.9	0.5 mg
63671-05	[Cys ²⁶]-β-Amyloid (1-40), S26C β-Amyloid (1-40)	DAEFRHDSGYEVHHQKLVFFAED-VGCNKGAIIGLMVGGVV	4345.9	0.5 mg
63671-1	[Cys ²⁶]-β-Amyloid (1-40), S26C β-Amyloid (1-40)	DAEFRHDSGYEVHHQKLVFFAED-VGCNKGAIIGLMVGGVV	4345.9	1 mg
62428	β-Amyloid (1-40)-Binding Peptide	AECDWKGKGRWRLWPGASGKTEACGP	2776.1	1 mg
62429	Cys-containing β-Amyloid (1-40) Binding Peptide, Biotin-labeled	Biotin-AECDWKGKGRWRLWPGASGKTEAC-GP	3002.4	1 mg
61484-01	β-Amyloid (1-42)-Lys(Biotin)-NH ₂	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA-K(Biotin)-NH ₂	4867.6	0.1 mg
61484-05	β-Amyloid (1-42)-Lys(Biotin)-NH ₂	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA-K(Biotin)-NH ₂	4867.6	0.5 mg
63319	β-Amyloid (1-42), 5-ROX-labeled	5-ROX-DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	5048.7	0.1 mg
62446	[Ala ²⁰]-β-Amyloid(1-42)	DAEFRHDSGYEVHHQKLVFAEDVG-SNKGAIIGLMVGGVVIA	4438.0	0.5 mg
63323	[Arg ⁶]-β-Amyloid (1-42), English Mutation	DAEFRDSDGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	4533.2	0.5 mg
61958-01	[Asn ^{6,13,14}]-β-Amyloid (1-42)	DAEFRNDSGYEVNNQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	4445.0	0.1 mg
63324	[Asn ⁷]-β-Amyloid (1-42), Tottori-Japanese Mutation	DAEFRHNSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	4513.1	0.5 mg
63705	[Asn ²³]-β-amyloid (1-42), Iowa Mutation	DAEFRHDSGYEVHHQKLVFFAENVG-SNKGAIIGLMVGGVVIA	4513.1	0.5 mg
62444	[Asp ³⁷]-β-Amyloid(1-42)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	4573.2	0.5 mg
61718-01	Biotin-LC-β-Amyloid (1-42), mouse, rat	Biotin-LC-DAEFGHDSGFVRRHQKLVFFAED-VGSNKGAIIGLMVGGVVIA	4757.5	0.1 mg
61718-05	Biotin-LC-β-Amyloid (1-42), mouse, rat	Biotin-LC-DAEFGHDSGFVRRHQKLVFFAED-VGSNKGAIIGLMVGGVVIA	4757.5	0.5 mg
62142	[Gln ²²]-β-Amyloid (1-42), E22Q Dutch Mutation	DAEFRHDSGYEVHHQKLVFFAQDVG-SNKGAIIGLMVGGVVIA	4513.1	0.5 mg
63704	[Gly ²¹]-β-Amyloid (1-42), A21G Flemish Mutation	DAEFRHDSGYEVHHQKLVFFGEDVG-SNKGAIIGLMVGGVVIA	4500.1	0.5 mg
61967-05	[Gly ²²]-β-Amyloid (1-42), Arctic Mutation	DAEFRHDSGYEVHHQKLVFFAGDVG-SNKGAIIGLMVGGVVIA	4442.1	0.5 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62832-01	[Gly ³⁵ , Asp ³⁷]-β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAI-IGLGVDGVVIA	4498.0	0.1 mg
62832-05	[Gly ³⁵ , Asp ³⁷]-β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAI-IGLGVDGVVIA	4498.0	0.5 mg
62143	[Leu ³⁵]-β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLLVGGVVIA	4496.1	0.5 mg
62148	[Lys ²²]-β-Amyloid (1-42), Italian Mutation	DAEFRHDSGYEVHHQKLVFFAKDVG-SNKGAIIGLMVGGVVIA	4513.2	0.5 mg
62478	[Met(O) ³⁵]-β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGL-M(O)-VGGVVIA	4530.1	0.5 mg
62443	[Pro ³¹]-β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAPIGLMVGGVVIA	4498.1	0.5 mg
62442	[Val ³³]-β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAI-IVLMVGGVVIA	4556.2	0.5 mg
63672-05	[Cys ²⁶]-β-Amyloid (1-42), S26C β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAED-VGCNKGAIIGLMVGGVVIA	4530.2	0.5 mg
63672-1	[Cys ²⁶]-β-Amyloid (1-42), S26C β-Amyloid (1-42)	DAEFRHDSGYEVHHQKLVFFAED-VGCNKGAIIGLMVGGVVIA	4530.2	1 mg
62479	β-Amyloid (1-43), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIAT	4519.1	1 mg
61966-01	β-Amyloid (1-44)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATV	4714.4	0.1 mg
61956-01	β-Amyloid (1-45)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVI	4827.5	0.1 mg
62991	β-Amyloid (1-45), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVI	4731.4	0.1 mg
62076-01	β-Amyloid (1-46)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIV	4926.6	0.1 mg
62992	β-Amyloid (1-46), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIV	4830.6	0.1 mg
61964-01	β-Amyloid (1-47)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVI	5039.8	0.1 mg
62993	β-Amyloid (1-47), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVI	4943.7	0.1 mg
61965-01	β-Amyloid (1-48)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVIT	5140.9	0.1 mg
62994	β-Amyloid (1-48), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVIT	5044.8	0.1 mg
61963-01	β-Amyloid (1-49)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITL	5254.1	0.1 mg
62979	β-Amyloid (1-49), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITL	5158	0.1 mg
63347	β-Amyloid (1-50), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITLV	5257.1	0.1 mg
63346	β-Amyloid (1-51), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITLVM	5588.3	0.1 mg
63345	β-Amyloid (1-52), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITLVML	5501.5	0.1mg
63344	β-Amyloid (1-53), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITLVMLK	5629.6	0.1 mg
63335	β-Amyloid (1-55)	DAEFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITLVMLKKK	5982.1	0.1 mg
63354	β-Amyloid (1-55), mouse, rat	DAEFGHDSGFVVRHQKLVFFAEDVG-SNKGAIIGLMVGGVVIATVIVITLVMLKKK	5886	0.1 mg
61029	β-Amyloid (3-40)	EFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVV	4143.7	0.1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61029-1	β-Amyloid (3-40)	EFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVV	4143.7	1 mg
63715-01	β-Amyloid (3-42)	EFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	4327.9	0.1 mg
63715-1	β-Amyloid (3-42)	EFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	4327.9	1 mg
61959-01	β-Amyloid (3-42)-Lys(Biotin)-NH ₂	EFRHDSGYEVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA-K(Biotin)-NH ₂	4682.4	0.1 mg
61979	Cys-β-Amyloid (4-10)	CFRHDSGY	984.1	1 mg
62954	β-Amyloid (4-13)	FRHDSGYEVH	1246.3	1 mg
62961	β-Amyloid (4-24)	FRHDSGYEVHHQKLVFFAEDV	2560.8	1 mg
63326	β-Amyloid (4-24)-Cys	FRHDSGYEVHHQKLVFFAEDVC	2664	1 mg
62965	β-Amyloid (5-14)	RHDSGYEVHH	1236.3	1 mg
62960	β-Amyloid (5-24)-Cys	RHDSGYEVHHQKLVFFAEDVC	2516.8	1 mg
63338	β-Amyloid (5-34), mouse, rat	GHDSGFVVRHQKLVFFAEDVGSNKGAIIGL	3228.6	1 mg
63341	β-Amyloid (5-36), mouse, rat	GHDSGFVVRHQKLVFFAEDVGSNKGAIIGLMV	3459	1 mg
62964	β-Amyloid (5-40)	RHDSGYEVHHQKLVFFAEDVGSNKGAIIGL-MVGGVV	3867.4	1 mg
63343	β-Amyloid (5-55), mouse, rat	GHDSGFVVRHQKLVFFAEDVGSNKGAIIGL-MVGGVVIATVIVITLVMLKKK	5423.5	0.1 mg
62888	β-Amyloid (6-30)	HDSGYEVHHQKLVFFAEDVGSNKG	2772.0	1 mg
61803	β-Amyloid (7-22)	DSGYEVHHQKLVFFAE	1906.1	1 mg
62889	β-Amyloid (7-29)	DSGYEVHHQKLVFFAEDVGSNKG	2563.8	1 mg
62887	β-Amyloid (8-28)	SGYEVHHQKLVFFAEDVGSNK	2391.6	1 mg
61879	β-Amyloid (8-38)	SGYEVHHQKLVFFAEDVGSNKGAIIGLMVGG	3260.7	1 mg
61975	β-Amyloid (8-40)	SGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVV	3459.0	1 mg
62996	β-Amyloid (8-40), mouse, rat	SGFEVRHQKLVFFAEDVGSNKGAIIGL-MVGGVV	3462	1 mg
62886	β-Amyloid (9-27)	GYEVHHQKLVFFAEDVGSN	2176.4	1 mg
62462	β-Amyloid (9-42)-Lys(Biotin)-NH ₂	GYEVHHQKLVFFAEDVGSNKGAIIGLMVGGV-VIA-K(Biotin)-NH ₂	3909.7	0.5 mg
62963	β-Amyloid (10-19)	YEVHHQKLVF	1299.5	1 mg
62972	Biotin-LC-β-Amyloid (10-20)	Biotin-LC-YEVHHQKLVFF	1786.2	1 mg
62885	β-Amyloid (10-26)	YEVHHQKLVFFAEDVGS	2005.2	1 mg
61931	β-Amyloid (10-35)	YEVHHQKLVFFAEDVGSNKGAIIGLM	2903.4	1 mg
63306	β-Amyloid (10-35), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-YEVHHQKLVFFAEDVG-SNKGAIIGLM	3259.8	0.1 mg
63307	β-Amyloid (10-35), HiLyte Fluor™ 555-labeled	HiLyte Fluor™ 555-YEVHHQKLVFFAEDVG-SNKGAIIGLM	3372	0.1 mg
62884	β-Amyloid (11-25)	EVHHQKLVFFAEDVG	1755.0	1 mg
62472	β-Amyloid (11-28)-Lys(Biotin)-NH ₂	EVHHQKLVFFAEDVGSNK-K(Biotin)-NH ₂	2437.8	0.5 mg
63339	β-Amyloid (11-28), mouse, rat	EVRHQKLVFFAEDVGSNK	2103.4	1 mg
62950	β-Amyloid (11-40), FAM-labeled	5-FAM-EVHHQKLVFFAEDVGSNKGAIIGLM-VGGVV	3510	0.1 mg
62951	β-Amyloid (11-40), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-EVHHQKLVFFAEDVG-SNKGAIIGLMVGGVV	3508.1	0.1 mg
62471	β-Amyloid (11-40)-Lys(Biotin)-NH ₂	EVHHQKLVFFAEDVGSNKGAIIGLMVGGVV-K(Biotin)-NH ₂	3505.2	0.5 mg
62997	β-Amyloid (11-40), mouse, rat	EVRHQKLVFFAEDVGSNKGAIIGLMVGGVV	3170.7	1 mg
63317	β-Amyloid (11-42)	EVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	3335.9	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
63328	β-Amyloid (11-42), FAM-labeled	5-FAM-EVHHQKLVFFAEDVGSNKGAIIGLM-VGGVVIA	3694.2	0.1 mg
63327	β-Amyloid (11-42), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-EVHHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA	3692.3	0.1 mg
62470	β-Amyloid (11-42)-Lys(Biotin)-NH ₂	EVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA-K(Biotin)-NH ₂	3689.4	0.5 mg
62939	β-Amyloid (12-22)	VHHQKLVFFAE	1354.6	1 mg
62883	β-Amyloid (12-24)	VHHQKLVFFAEDV	1568.8	1 mg
63342	β-Amyloid (12-33), mouse, rat	VRHQKLVFFAEDVGSNKGAIIG	2385.8	1 mg
63340	β-Amyloid (12-38), mouse, rat	VRHQKLVFFAEDVGSNKGAIIGLMVGG	2843.4	1 mg
62452	β-Amyloid (12-42)	VHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA	3206.8	1 mg
62894	β-Amyloid (13-22)	HHQKLVFFAE	1255.5	1 mg
62882	β-Amyloid (13-23)	HHQKLVFFAED	1370.5	1 mg
61802	β-Amyloid (13-28)	HHQKLVFFAEDVGSNK	1856.1	1 mg
61800	β-Amyloid (13-29)-Gly-Cys	HHQKLVFFAEDVGSNKGGC	2073.3	1 mg
62962	β-Amyloid (13-40)	HHQKLVFFAEDVGSNKGAIIGLMVGGVV	2923.4	1 mg
62893	β-Amyloid (14-21)	HQKLVFFA	989.2	1 mg
62881	β-Amyloid (14-22)	HQKLVFFAE	1118.3	1 mg
62024	β-Amyloid (14-23)	HQKLVFFAED	1233.4	1 mg
62891	[Gln ²²]-β-Amyloid (14-23)	HQKLVFFAQD	1232.4	1 mg
62760	β-Amyloid (14-25)	HQKLVFFAEDVG	1389.6	1 mg
62880	β-Amyloid (15-23)	QKLVFFAED	1096.3	1 mg
62892	[Gln ²²]-β-Amyloid (15-23)	QKLVFFAQD	1095.3	1 mg
62906	β-Amyloid (15-25)	QKLVFFAEDVG	1252.4	1 mg
62907	Biotin-LC-β-Amyloid (15-25)	Biotin-LC-QKLVFFAEDVG	1591.9	1 mg
62469	β-Amyloid (15-26)-Lys(Biotin)-NH ₂	QKLVFFAEDVGS-K(Biotin)-NH ₂	1693.0	0.5 mg
62023	β-Amyloid (15-36)	QKLVFFAEDVGSNKGAIIGLMV	2336.8	1 mg
61344	[Ac-Lys(Me) ¹⁶ , N-Me-Val ¹⁸]-β-Amyloid (16-20)-NH ₂	Ac-K(Me)-L-(N-Me-V)-FF-NH ₂	721.9	1 mg
61343	β-Amyloid (16-20), all d-isomers	klvff	652.8	1 mg
63304	β-Amyloid (16-20), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-KLVFF	1009.2	1 mg
63305	β-Amyloid (16-20), HiLyte Fluor™ 555-labeled	HiLyte Fluor™ 555-KLVFF	1121.4	1 mg
62468	β-Amyloid (16-20)-Lys(Biotin)-NH ₂	KLVFF-K(Biotin)-NH ₂	1006.3	1 mg
62897	β-Amyloid (16-22)	KLVFFAE	853	1 mg
61801	β-Amyloid (16-23)	KLVFFAED	968.1	1 mg
62139	β-Amyloid (16-24)	KLVFFAEDV	1067.3	1 mg
61804	β-Amyloid (16-26)	KLVFFAEDVGS	1211.4	1 mg
62467	β-Amyloid (17-21)-Lys(Biotin)	LVFFA-K(Biotin)	950.2	1 mg
62477	Myristoyl-β-Amyloid(17-21)	Myristoyl-LVFFA	788.1	1 mg
61978	β-Amyloid (17-24)	LVFFAEDV	939.1	1 mg
61980	Biotin-β-Amyloid (17-24)	Biotin-LVFFAEDV	1165.4	1 mg
62447	β-Amyloid (17-28)	LVFFAEDVGSNK	1325.5	1 mg
62466	β-Amyloid (17-28)-Lys(Biotin)	LVFFAEDVGSNK-K(Biotin)	1680.0	0.5 mg
62975	β-Amyloid (17-40)-KKK-Lys(Biotin)-NH ₂	LVFFAEDVGSNKGAIIGLMVGGVKKK-K(Biotin)-NH ₂	3130.9	0.1 mg
62448	Biotin-β-Amyloid (18-28)	Biotin-VFFAEDVGSNK	1438.6	1 mg
62971	Biotin-LC-β-Amyloid (18-28)	Biotin-LC-VFFAEDVGSNK	1551.8	1 mg
61989	β-Amyloid (20-42)	FAEDVGSNKGAIIGLMVGGVVIA	2217.6	1 mg
62453	β-Amyloid (22-40)	EDVGSNKGAIIGLMVGGVV	1815.1	1 mg
62454	β-Amyloid (22-40)-NH ₂	EDVGSNKGAIIGLMVGGVV-NH ₂	1814.1	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62455	Biotin-LC-β-Amyloid (22-41)	Biotin-LC-EDVGSNKGAIIGLMVGGVVI	2267.8	1 mg
61972	β-Amyloid (22-42)	EDVGSNKGAIIGLMVGGVVIA	1999.4	1 mg
62761	β-Amyloid (23-37)	DVGSNKGAIIGLMVG	1430.7	1 mg
62766	β-Amyloid (23-42)	DVGSNKGAIIGLMVGGVVIA	1870.3	1 mg
63308	β-Amyloid (25-35), HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-GSNKGAIIGLM	1416.7	0.1 mg
61971	β-Amyloid (25-35), scrambled	MAKGINGISGL	1060.3	1 mg
62451	Biotin-β-Amyloid (25-35)	Biotin-GSNKGAIIGLM	1286.6	1 mg
61984	[Phe ³⁴]-β-Amyloid (25-35)	GSNKGAIIGFM	1094.3	1 mg
62920	β-Amyloid (26-37)	SNKGAIIGLMVG	1159.4	1 mg
62919	β-Amyloid (26-38)	SNKGAIIGLMVGG	1216.5	1 mg
62916	β-Amyloid (26-39)	SNKGAIIGLMVGGV	1315.6	1 mg
61985	β-Amyloid (26-40)	SNKGAIIGLMVGGVV	1414.7	1 mg
62915	β-Amyloid (26-42)	SNKGAIIGLMVGGVVIA	1599	1 mg
62917	β-Amyloid (26-43)	SNKGAIIGLMVGGVVIAT	1700.1	1 mg
61990	β-Amyloid (28-40)	KGAIIGLMVGGVV	1213.6	1 mg
61976	β-Amyloid (28-42)	KGAIIGLMVGGVVIA	1397.8	1 mg
61974	β-Amyloid (28-43)	KGAIIGLMVGGVVIAT	1498.9	1 mg
61988	β-Amyloid (28-40)-Gly-Gly-Gly-Cys	KGAIIGLMVGGVVGCGC	1487.9	1 mg
62924	β-Amyloid (29-37)	GAIIGLMVG	830.1	1 mg
62923	β-Amyloid (29-38)	GAIIGLMVGG	887.1	1 mg
62921	β-Amyloid (29-39)	GAIIGLMVGGV	986.3	1 mg
62449	Biotin-β-Amyloid (29-40)	Biotin-GAIIGLMVGGVV	1311.7	1 mg
61987	β-Amyloid (29-42)	GAIIGLMVGGVVIA	1269.6	1 mg
61986	Biotin-β-Amyloid (29-42)	Biotin-GAIIGLMVGGVVIA	1495.9	1 mg
61977	β-Amyloid (30-34)	AIIGL	485.6	1 mg
61982	β-Amyloid (33-40)	GLMVGGVV	730.9	1 mg
61981	β-Amyloid (35-42)	MVGGVVIA	745.0	1 mg
62759	β-Amyloid (35-43)	MVGGVVIAT	846.1	1 mg
63747	β-Amyloid (36-42)	VGGVVIA	613.8	1 mg
63749	β-Amyloid (36-44)	VGGVVIATV	814.0	1 mg
63750	β-Amyloid (36-45)	VGGVVIATVI	927.2	1 mg
62533	Cys-β-hairpin (BHA) Protein G B1 Domain (41-56)	CGEWYDDATKTFVTVE	1864	1 mg
61983	Amyloid Precursor Protein (APP) (44-62)	HMNQNGKWSDPSGKTC	2105.3	1 mg
62913	Amyloid Precursor-Like Protein 1 (APLP1) (594-610)	GGGSLIVLSLLLLRKKK	1795.3	1 mg
62289	[Phe ⁶⁶⁸]-Amyloid Precursor Protein (APP) (659-676)	GVVEVDAAVFPEERHLSK	1982.2	1 mg
62065	Amyloid Precursor Protein (APP) (667-676)	SEVKMDAEFR	1211.4	1 mg
62059	Amyloid Precursor Protein (APP) (667-676)-Cys	SEVKMDAEFRC	1314.5	1 mg
62074	Amyloid Precursor Protein (APP) (667-676), Dnp-labeled	SEVKMDAEFR-DAP(Dnp)-KK	1719.9	1 mg
62075	Amyloid Precursor Protein (APP) (667-676), FRET	Mca-SEVKMDAEFR-DAP(Dnp)-NH ₂	1678.8	1 mg
62060	Amyloid Precursor Protein (APP) (667-676), Swedish Mutation	SEVNLDAEFR	1179.3	1 mg
62515	[Asn ⁶⁷⁰ , Leu ⁶⁷¹]-Amyloid Precursor Protein (APP) (668-675), Swedish Mutation	EVNLDAEF	936	1 mg
62236	[Asn ⁶⁷⁰ , Leu ⁶⁷¹]-RE(EDANS)-APP(668-675)-Lys(DABCYL)-R	R-E(EDANS)-EVNLDAEF-K(DABCYL)-R	2005.3	0.5 mg
62762	Amyloid-Forming Peptide, GNNQQNY	GNNQQNY	836.8	1 mg
62895	β-Amyloid-Related Death-Inducing Protein (A-β-DIP) (768-776)	MLTRILVIREIQEIAIVANASTMN	2658.2	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62896	β-Amyloid-Related Death-Inducing Protein (A-β-DIP) (183-196)	AEIPEEHIPHQIQ	1668.8	1 mg
62834	Amyloid Toxicity Inhibitor Peptide 11	RGTFEQKF-NH ₂	940.1	1 mg
62835	Amyloid Toxicity Inhibitor Peptide 13	RGTLEGKL-NH ₂	872	1 mg
62837	Amyloid Toxicity Inhibitor Peptide 110	RGTWEGKW-NH ₂	1018.2	1 mg
62838	Amyloid Toxicity Inhibitor Peptide 111	RGTYEGKY-NH ₂	972.1	1 mg
62912	Amyloid Precursor-Like Protein 2 (APLP2) (706-721)	AIATVIVISLVMLRKR	1783.3	1 mg
63322	Amyloid Precursor Protein, (APP) (721-770)	VMLKKKQYTSIHGGVVEVDAAVTPEERHLSK-MQQNGYENPTYKFFEQMQN	5910.7	0.5 mg
62073	Amyloid Precursor Protein, (APP) (740-770)	AAVTPEERHLSKMQQNGY-ENPTYKFFEQMQN	3717.1	1 mg
63303	Amyloid Precursor Protein, (APP) (741-770)	AVTPEERHLSKMQQNGYENPTYKFFEQMQN	3646.1	1 mg
63302	Cys-Amyloid Precursor Protein, (APP) (751-770)	CKMQQNGYENPTYKFFEQMQN	2629	1 mg
63301	Amyloid Precursor Protein, (APP) (756-762)	Ac-GYENPTY-NH ₂	883.9	1 mg
61805	Anti-parallel Topology β-Amyloid Modified Peptide	CHQKLFFAEDYNGKHQKLVFFAED	3014.4	1 mg
62531	Betanova (BET)	RGWSVQNGKYTNNGKTTEGR	2253.4	1 mg
63737	ClearPoint™ β-Amyloid (1-40), ¹³ C, ¹⁵ N-labeled at Arg & Lys	DAEF-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVV [R*=R(U- ¹³ C ₆ , U- ¹⁵ N ₄) & K*=K(U- ¹³ C ₆ , U- ¹⁵ N ₂)]	4355.9	50 µg
63740	ClearPoint™ β-Amyloid (1-40), U- ¹³ C-labeled at Leu ¹⁷	DAEFRHDSGYEVHHQK-L*-VFFAEDVG-SNKGAIIGLMVGGVV [L*=L(U- ¹³ C ₆)]	4335.9	50 µg
63738	ClearPoint™ β-Amyloid (1-42), ¹³ C, ¹⁵ N-labeled at Arg & Lys	DAEF-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVVIA [R*=R(U- ¹³ C ₆ , U- ¹⁵ N ₄) & K*=K(U- ¹³ C ₆ , U- ¹⁵ N ₂)]	4540.1	50 µg
63739	ClearPoint™ β-Amyloid (4-42), ¹³ C, ¹⁵ N-labeled at Arg & Lys	F-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVVIA [R*=R(U- ¹³ C ₆ , U- ¹⁵ N ₄) & K*=K(U- ¹³ C ₆ , U- ¹⁵ N ₂)]	4224.8	50 µg
63741	ClearPoint™ β-Amyloid (4-42), U- ¹³ C, ¹⁵ N -labeled at Val ¹²	FRHDSGYE-V*-HHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA [V*=V(U- ¹³ C ₅ , U- ¹⁵ N)]	4204.8	50 µg
63742	ClearPoint™ β-Amyloid (8-42), U- ¹³ C, ¹⁵ N -labeled at Val ¹²	SGYE-V*-HHQKLVFFAEDVGSNKGAIIGLMVGGVVIA [V*=V(U- ¹³ C ₅ , U- ¹⁵ N)]	3649.2	50 µg
61930	Colivelin	SALLRSIPAPAGASRLLLLTGEIDL	2645.1	1 mg
62905	Cys-NC2-2 CLAC-P (155-169)	CKGEQGDQGPRMVFQK	1777.1	1 mg
62080	DA-A, BACE Substrate	VVEVDAAVTP	999.1	1 mg
61927	[D-Ser ¹⁴] Humanin (HN)	MAPRGFSCLLLLTsEIDLQVKKRA	2687.3	1 mg
61928	AGA-(C8R)HNG17, Humanin Derivative	PAGASRLLLLTGEIDL	1736.1	1 mg
62903	NC1 CLAC-P (14-27)	EPRSEDPTPAEQHC	1595.7	1 mg
62904	NC2-1 CLAC-P (171-185)	GCNHGFLSADQQLIK	1630.9	1 mg
62901	NC3 CLAC-P (430-443)	DYNGNLHEALQRITC	1746.9	1 mg
62902	NC4 CLAC-P (641-654)	LPGDGLPMPGCWQK	1498.8	1 mg
61807	Parallel topology β-Amyloid modified peptide	CHQKLFFAEDYNGKDEAFFVLKQHW	3200.6	1 mg
62922	Serum Amyloid P-Component	FTLCFR	786	1 mg
62959	Tau-Protein (1-16)	MAEPRQEFVEMDHAG	1876.1	1 mg
62958	Tau-Protein (323-335)	GSLGNIHHPGGG	1230.4	1 mg

Angiotensins and Related Peptides

63702	[Des-Asp ¹]-Angiotensin I, human	RVYIHPFHL	1181.4	1 mg
61182	ClearPoint™ Angiotensin I, human, ¹³ C and ¹⁵ N-labeled	DRVY-I*-HPFHL [I*=I(U- ¹³ C ₆ , ¹⁵ N)]	1303.5	0.1 mg
61901	Angiotensin II Receptor, AT2, Amino Terminal Fragment	MKDNFSFAATSRNITSS	1877.1	1 mg
61873	Angiotensin II type 1 receptor (181-187), AT1, ATE.	AFHYESQ	880.9	1 mg
61902	AT1A, Angiotensin II receptor (225-237)	AYEIQKNKPRNDD	1590.7	1 mg
63703	[Sar ¹ , Ala ⁸]-Angiotensin II	Sar-RVYIHPA	926.1	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62337	DX 600, ACE2 Inhibitor	Ac-GDYSHCSPLRYYPWWKCTYPDPEGGG-NH ₂	3076.4	0.1 mg
61872-01	Peptide Substrate for Renin 520 Assay kit		2000 - 2200	0.1 mg
61872	Peptide Substrate for Renin 520 Assay kit		2000 - 2200	1 mg
62334-01	Rat Renin FRET Substrate (5-FAM/QXL™520)		2000 - 2200	0.1 mg
62334	Rat Renin FRET Substrate (5-FAM/QXL™520)		2000 - 2200	1 mg
62022	Renin Substrate 1, FRET	R-E(EDANS)-IHPFHLVIHT-K(DABCYL)-R	2282.7	1 mg
72065	Renin Inhibitor III	RRPFH-Sta-IHK-NH ₂	1247.6	1 mg
72066	Renin Inhibitor III	RRPFH-Sta-IHK-NH ₂	1247.6	5 mg
62118	HRP Prorenin, Decoy Handle Region Peptide (31-40)	RILLKKMPVS	1184.6	1 mg
62119	HRP Prorenin Peptide, scrambled (33-42)	RIFLKRMPSI	1260.6	1 mg

Antimicrobial and Related Peptides

61891	Acetyl-Adhesin (1025-1044), amide	Ac-QLKTADLPAGRDETTSFVLV-NH ₂	2202.5	1 mg
62783	Alloferon 1	HGVSGHGQHGCVHG	1265.3	1 mg
62784	Alloferon 2	GVSFGHGQHGCVHG	1128.2	1 mg
62046	Andropin	VFIDILDKVENAIHNAAQVGIGFAKPFKELINPK	3750.4	1 mg
62045	Apidaecin IA	GNNRPVYIPQPRPPHPRI	2108.4	1 mg
62044	Apidaecin IB	GNNRPVYIPQPRPPHPRL	2108.4	1 mg
62043	Aurein 1.1	GLFDIIKKIAESI-NH ₂	1445.8	1 mg
61442	BTM-P1, <i>Bacillus thuringiensis</i> Protein P1	VAPIAKYLATALAKWALKQGFAKLKS	2788.4	1 mg
61915-05	Carnobacteriocin B2	VNYGNGVSCSKTKCSVNWGQAFQERY-TAGINSFVSGVASGAGSIGRRP (Disulfide bridge: 9-14)	4967.5	0.5 mg
61307	CAP-18, Cathelicidin-Derived Peptide-18, rabbit	GLRKRLRKFNRNIIKEKLLKIGQKIQLLP-KLAPRTDY	4433.5	1 mg
62649	κ Casein A (138-158), bovine	AVESTVATLEDSPEVIESPPE	2199.4	1 mg
63731	κ Casein (106-115), bovine	MAIPPKKNQD	1123.4	1 mg
63730	κ Casein (138-158, pSer ¹⁴⁹)	AVESTVATLEA-pS-PEVIESPPE	2235.3	1 mg
61881	DDDDDDD, Antimicrobial Anionic Peptide, Surfactant, ovine	DDDDDDD	822.7	1 mg
61913-05	Dermaseptin	ALWKTMLKKLGTMLHAGKAAL-GAAADTISQGTQ	3455.1	0.5 mg
61910	[Des Ala ²⁰ , Gln ³⁴] Dermaseptin	ALWKTMLKKGTMLHAGKALWKTMLKKGTM-ALHAGK	1985.5	1 mg
63713	Dermcidin, DCD-1L	SSLLEKGLDGAKKAVGGLGKLGKDAVEDLES-VGKGAVHDVKDVLDSVL	4818.5	0.1 mg
62600	Drosocin	GKPRPYSRPTSHRPIRV	2198.6	1 mg
61641	Elafin	AQEPVKGPVSTKPGSCPIILIRCAML-NPPNRCLKDTCPGIKKCCGSCGMACFVPQ (Disulfide bridges: 16-45, 23-49, 32-44, 38-53)	5999.3	0.1 mg
61880	GADDDDD, Antimicrobial Anionic Peptide, Surfactant, ovine	GADDDDD	721.6	1 mg
61359	Histatin-3 (H3)	DSHAKRHHGYKRKFHEKHHSHRGYRSNYL-YDN	4062.4	1 mg
62412	Human Platelet Factor IV 18, C18G	ALYKLLKLLKLSAKKLG	2043.7	1 mg
62514	Innate Defense Regulator-1 (IDR-1)	KSRIVPAIPVSL	1392.8	1 mg
62651	Lactoferricin B, Lactoferrin (17-41)	FKCRRWQWRMKKLGAPSITCVRRAF (S-S BOND)	3123.9	1 mg
61432	LEAP-1, Hecpidin, human	DTHFPICIFCCGCHRSKCGMCKCT (Disulfide bridges: 7-13, 10-23, 11-22, 14-19)	2789.5	0.1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61640	LEAP-2 (38-77), human	MTPFWRGVSLRPIGASCRDDSECITRL-CRKRRCLSLSVAQE	4585.4	1 mg
62366	Melittin, honey bee	GIGAVLKVLTTGLPALISWIKRKRQQ	2847.5	1 mg
62603	Metalnikowin I	VDKPDYRPRPRPPNM	1838.1	1 mg
62602	Metalnikowin IIA	VDKPDYRPRPWPRPN	1893.2	1 mg
61918	Mundtacin	KYYGNGVSCNKKGCSVDWGKAIGIIGNNSAANLATGGAAGWSK (Disulfide bridge: 9-14)	4287.8	0.5 mg
61917-05	Mundtacin KS	KYYGNGVSCNKKGCSVDWGKAIGIIGNNSAANLATGGAAGWKS (Disulfide bridge: 9-14)	4287.8	0.5 mg
61310	OV-1, Ovispirin-1, sheep	KNLRRRIIRKIIHIIKKYG	2262.9	1 mg
61311	OV-2, Ovispirin-2, sheep	LRRRIIRKIIHIIKK-NH ₂	1799.4	1 mg
61312	OV-3, Ovispirin-3, sheep	IRRIIRKIIHIIKK-NH ₂	1799.4	1 mg
61916-05	Piscicolin 126	KYYGNGVSCNKGCTVDWSKAIGIIGNNAANLTTGGAAGWNKG (Disulfide bridge: 9-14)	4415.9	0.5 mg
62130	<i>Plasmodium yoelii</i> Circumsporozoite Protein, (PyCSP), (57-70)	KIYNRNIVNRLLDG	1688.0	1 mg
62087	<i>Plasmodium yoelii</i> Circumsporozoite Protein, (PyCSP), (280-288)	SYVPSAEQI	993.1	1 mg
62047-01	Plectasin	GFGCNGPWDRDEDDMQCHNHCLSIKGYKG-GYCAKGGFVCKCY (With three disulfide bridges)	4658.3	0.1 mg
63690	PR-19	RRRPRPPYLPRPRPPFFFP	2399.9	1 mg
63313	Pseudin-2	GLNALKKVFQGIHEAIKLINNHVQ	2685.2	1 mg
62601	Pyrrhocoricin	VDKGSYLPRPTPRPIYNRN	2340.7	1 mg
62652	Serorphin, BSA (399-404)	YGFQNA	698.7	1 mg
61308	SMAP 29, Sheep Myeloid Antimicrobial Peptide 29	RGLRRLGRKIAHGKYGPTVLRRIIAG	3256.0	1 mg
61352	SP2, <i>C. albicans</i> Mating Pheromone	GFRLTNFGYFEPGK	1632.9	1 mg
61440	Theromacin (49-63)	QRLPNKQCRCINAR (Disulfide bridge 57-59)	1812.1	1 mg
62587	Ubiquitin (1-34)	MQIFVKLTGTITLEVEPSDTIENVKAKIQDKE	3848.5	1 mg
62586	Ubiquitin (65-76), Ub2	STLHLVLRRLGG	1321.6	1 mg

Apelin Peptides

63699	Apelin-16, human, bovine	FRRQRPRLSHGKMPF	2010.4	1 mg
63701	Apelin-36, bovine	LVQPRGPRSGPGPWQGGRRKFRQR-PRLSHKGPMPF	4178.9	1 mg

Apoptosis Peptides

62198	CGDKAKGRC-GG-klaklaklaklak, Obesity and Apoptosis Related Peptide (OAR), scrambled	CGDKAKGRC-GG-klaklaklaklak-NH ₂ (Disulfide bridge: 1-9)	2554.2	1 mg
62201	CKGGRAKDC, Obesity Related Peptide	CKGGRAKDC-NH ₂ (Disulfide bridge: 1-5)	934.1	1 mg
62206	Pro-apoptotic Peptide, klaklaklaklak, FAM-labeled	5-FAM-klaklaklaklak-NH ₂	1881.3	1 mg
62371	DQ (65-79), α Chain HLA II fragment	NIAVLKHNLNIVIKR	1745.2	1 mg
62403	Noxa BH3, Peptide 2	VPADLKDECAQLRRIGDKVNLRQKL	2879.4	1 mg
62199	(klaklak) ₂ , Proapoptotic Peptide	klaklaklaklak-NH ₂	1523.0	1 mg
62207	RGD-targeted Proapoptotic Peptide	ACDCRGDCFC-GG-klaklaklaklak-NH ₂ (S-S bonded C1-C4 & C2-C3)	2707.4	1 mg
63688-05	PR39, Anti-Apoptotic Factor	RRRPRPPYLPRPRPPFFPPRLPPRIPPGFP-PRFPPRFP	4720.7	0.5 mg
63689-05	PR39-S0, Scrambled PR39	RRRPRFPFGPIPPFPPLRPPFPFPPFR-PRPYRPRLPR	4720.7	0.5 mg
63689-1	PR39-S0, Scrambled PR39	RRRPRFPFGPIPPFPPLRPPFPFPPFR-PRPYRPRLPR	4720.7	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62404	PUMA BH3	EEQWAREIGAQLRRMADDLNAQYER	3049.3	1 mg
Aquaporin Peptides				
61324	Aquaporin-2 (254-267), human	RQSVELHSPQSLPR	1633.8	1 mg
61325	[pSER ²⁵⁶]-Aquaporin-2 (254-267), human	RQ-pS-VELHSPQSLPR	1713.8	1 mg
61328	[pSER ²⁶¹]-Aquaporin-2 (254-267), human	RQSVELH-pS-PQSLPR	1713.8	1 mg
61326	Aquaporin-2 (255-271), human	QSVELHSPQSLPRGSKA	1821.0	1 mg
61327	Aquaporin-2 (255-271), rat	QSVELHSPQSLPRGTGA	1835.1	1 mg
Arg-Phe-Amide (RFamide) Related Peptides				
61649	26Rfa, Hypothalamic RFamide Peptide, frog	VG TALGSLAEELNGYNRKKGGFSFRF-NH ₂	2347.0	1 mg
61651	26Rfa, Hypothalamic RFamide Peptide, human	TSGPLGNLAEELNGYSRKKGGFSFRF-NH ₂	2832.2	1 mg
61650	26Rfa, Hypothalamic RFamide Peptide, rat	ASGPLGTLAEELSSYSRRKGGFSFRF-NH ₂	2820.2	1 mg
61438	PQRFamide, Lamprey	SWGAPAEKFWMRAMPQRF-NH ₂	2195.6	1 mg
61851	Neuropeptide AF (hNPAF), human	AGEGLNSQFWSLAAPQRF-NH ₂	1978.2	1 mg
Autoimmunity Related Peptides				
62726	Autoimmune Lupus Erythematosus Epitope 1	PPPGIRGP	789.9	1 mg
61514	GAD65, Glutamic Acid Decarboxylase (78-97)	KPCNCPKGDVNYAFLHATDL	2206.5	1 mg
61515	GAD65, Glutamic Acid Decarboxylase (206-220)	TYEIAPVFVLLLEYVT	1757.1	1 mg
62727	GAD65, Glutamic Acid Decarboxylase (524-538)	SRLSKVAPVIKARMM	1687.2	1 mg
62758	GAD65, Glutamic Acid Decarboxylase (530-543)	APVIKARMMMEYGT	1567.9	1 mg
62089	H/K-ATPase α -chain Peptide (889-900), PLL (889-900)	PLLCVGLRPQWE	1410.7	1 mg
62088	H ⁺ /K ⁺ -ATPase α -chain Peptide (630-641), PIT (630-641)	PITAKAIAASVG	1098.3	1 mg
Bacterial Peptides				
62231	Bacterial Sortase Substrate I, FRET	DABCYL-LPETG-EDANS	1015.2	1 mg
62107	CAT Peptide, Catalytic Region of Streptococcal Glucosyltransferases (442-462)	DANFDSIRVDAVDNVDADLLQ	2305.4	1 mg
61689	CFP 10, Culture Filtrate Protein 10 (71-85)	EISTNIRQAGVQYSR	1721.9	1 mg
61810-1	CFTR, Cystic Fibrosis Transmembrane Conductance Regulator (108-117), <i>P. aeruginosa</i> Inhibitor	SYDPDNKEER	1252.3	1 mg
61810-5	CFTR, Cystic Fibrosis Transmembrane Conductance Regulator (108-117), <i>P. aeruginosa</i> Inhibitor	SYDPDNKEER	1252.3	5 mg
62399	CRGKA, Giardia Variable Surface Proteins Conserved Region	CRGKA	533.7	1 mg
62634	flg15, Flagellin Fragment	RINSAKDDAAGLQIA	1542.7	1 mg
62635	flg15- δ -4, Flagellin Fragment	RINSAKDDAAG	1117.2	1 mg
62108	GLU Peptide, Glucan-Binding Region of Streptococcal Enzymes	TGAQTIKQKLYFKANGQQVKG	2365.7	1 mg
62818	Lysosomal Integral Membrane Protein LIMP-2	CDEGTADERAPLIRT	1646.8	1 mg
62710	Malaria Circumsporozoite Protein, CSP (334-342)	YLKKIKNSL	1106.4	1 mg
62425	<i>Mycobacterium tuberculosis</i> Protein Ag85A	LTSELPGWLQANRHVKPTGS	2191.5	1 mg
62424	<i>Mycobacterium tuberculosis</i> Protein Ag85A-CD8	MPVGGQSST	863	1 mg
62550	<i>P. falciparum</i> Liver-Stage Antigen 3-NRI (LSA3-NRI)	DELFNELLNSVDVNGENILEESQ	2620.8	1 mg
62551	<i>P. falciparum</i> Liver-Stage Antigen 3-NRII (LSA3-NRII) (81-106)	LEESQVNDIFNSLVKSVQQEQQHNV	3028.3	1 mg
62549	<i>P. falciparum</i> Liver-Stage Antigen 3-RE (LSA3-RE) (183-210)	VESVAPSVEESVAPSVEESVAENVEESV	2888.1	1 mg
Bcl / BAD Peptides				
62271	BAD (CT-1)	KKGLPRPKSAGTATQMRQSSSWTC-NH ₂	2606	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62270	BAD (NT-1)	PEFEPSEQEDSSSAERC-NH ₂	1926	1 mg
62322	Bad (103-127), human, all d-isomers	nlwaaqrygrellrmsdefvdsfkk	3103.5	1 mg
62489	BAD BH3 (103-123)	NLWAAQRYGRELRRMSDEFVD	2612.9	1 mg
62483	BAD BH3 (146-159)	RYGRELRRMSDEFE	1844.1	1 mg
62482	[pSer ¹⁵⁵]-BAD BH3 (146-159)	RYGRELRRM-pS-DEFE	1924.1	1 mg
62269	BAD Peptide, Biotin-LC-labeled	Biotin-LC-AGAGRSRHSSYPAGT	1814.1	1 mg
62523	BAFF-R (160-183)	SVPVPATELGSTELVTTKTAGPEQ	2412.7	1 mg
61616	Bak-BH3	GQVGRQLAIIGDDINR	1724.9	1 mg
62492	Bak-BH3 (67-87)	PSSTMGQVGRQLAIIGDDINR	2228.5	1 mg
62272	Bak-BH3 (69-93)	STMGQVGRQLAIIGDDINRRYDSEF	2842.2	1 mg
62261	Bak-BH3 (71-89)	MGQVGRQLAIIGDDINRRY	2175.5	1 mg
62262	Antennapedia Bak BH3 (Ant-BH3) (71-105) Fusion Peptide	RQIKIWFQNRMMKWKMGQVGRQLAI-IGDDINRRY	4404.3	1 mg
62480	Bak-BH3 (73-87)	QVGRQLAIIGDDINR	1667.9	1 mg
62235	Bak-BH3, Bcl2 (72-90)	KGGGQVGRQLAIIGDDINR	1967.2	1 mg
62029	Bak-BH3, Mca-labeled	Mca-GQVGRQLAIIGDDINR	1941.1	1 mg
62030	Bak-BH3, TAMRA-labeled	GQVGRQLAIIGDDINRK(TAMRA)-NH ₂	2265.6	0.5 mg
61629	Bax-BH3	KKLSECLKRIGDELDS	1834.1	1 mg
61630	Bax-BH3 L63A	KKLSECAKRIGDELDS	1792.1	1 mg
62486	Bax-BH3 (58-71)	KLSECLKRIGDELD	1618.9	1 mg
62267	[Glu ⁶³] Bax BH3, mutant	STKKLSECEKRIGDELDSNM	2283.6	1 mg
62266	Bax BH3 Peptide (55-74), wild type	STKKLSECLKRIGDELDSNM	2267.6	1 mg
62283	Bax H2-H3 (53-86), Helix 2-3	DASTKKLSECLKRIGDELDSNMELQRMI-AAVDTD	3797.3	1 mg
62330	[Arg ⁶⁷]-Bax H2-H3 (53-86), [Arg ⁶⁷] Helix 2-3 (53-86), mutant	DASTKKLSECLKRIRDELDSNMELQRMI-AAVDTD	3896.5	1 mg
62495	Bax H3 (71-86), Helix 3 (71-86)	DSNMELQRMIAAVDTD	1809.5	1 mg
62268	Bax I	PQDASTKKLSECLKRIGDELDSNMEL	2808.2	1 mg
62274	Bcl 9-2	GSEGLSKEQLEHRERSLQTLRDIERLLLRSGETEPFLKGGPPGGAG-NH ₂	4944.6	0.5 mg
62275	Bcl 9-2, mutant	GSEGLSKEQLEHRERSFQTLRDIERLLLRSGETEPFLKGGPPGGAG-NH ₂	4978.6	0.5 mg
62494	Bcl-2 BH3 (85-105)	ALSPVPVHLLTRQAGDFSRR	2319.7	1 mg
62273	Bcl-2 BH3 Peptide II	LSPVPPVHLLRQAGDDFSRRYRG	2807.2	1 mg
62481	Bcl-XL BH3 (85-98)	AVKQALREAGDEFE	1562.7	1 mg
61516	BDC2.5(A)	GKKVAAPAWARMG	1342.6	1 mg
62439	BH3 BIM Peptide (52-71), human	Ac-MRPEIWIWAQELRRIGDEFNA	2485.9	1 mg
62321	[Cys ⁹⁶] Bid BH3, mouse	ESQEEIHHNIARHLAQIGDCMDHN	2774.1	1 mg
62317	Bid BH3 (77-100), mouse	ESQEEIHHNIARHLAQIGDEMDHN	2800	1 mg
62318	[Ala ^{90,95}] Bid BH3 (77-100), mouse	ESQEEIHHNIARHAAQIGAEMDHN	2713.9	1 mg
62320	[Cys ⁹²] Bid BH3 (77-100), mouse	ESQEEIHHNIARHLACIGDEMDHN	2775	1 mg
62319	[Cys ^{92,96}] Bid BH3 (77-100), mouse	ESQEEIHHNIARHLACIGDCMDHN	2749.1	1 mg
62488	Bid BH3 (79-99)	QEDIIRNIARHLAQVGDSDMR	2437.7	1 mg
61711	Bid BH3 (80-99), human	EDIIRNIARHLAQVGDSDMR	2309.6	1 mg
61631	Bid BH3 (84-99), human	RNIARHLAQVGDSDMR	1839.1	1 mg
61712	Bid BH3 (84-99), human, FAM-labeled	5-FAM-EDIIRNIARHLAQVGDSDMR	2667.9	0.5 mg
62485	Bid BH3 (85-98)	NIARHLAQVGDSDMD	1526.7	1 mg
61715	Bid BH3-r8	rrrrrrrr-GEDIIRNIARHLAQVGDSDMR	3616.2	0.5 mg
62057	Bid BH3-R8	RRRRRRRRGEDIIRNIARHLAQVGDSDMR	3616.2	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61713	Bid BH3-R9	RRRRRRRRRRGEDIIRNIARHLAQVGDSMDR	3772.3	1 mg
62277	Bid BH3, Peptide II, TAMRA-labeled	5-TAMRA-EDIIRNIARHLAQVGDSMDR	2722.1	1 mg
62490	Bik BH3 (50-70)	CMEGSDALALRLACIGDEMDV	2212.6	1 mg
62487	Bik BH3 (56-69)	ALALRLACIGDEMD	1490.8	1 mg
62278	Bim BH3 (81-105), human, Peptide III	DMRPEIWIAQELRRIGDEFNAYYAR	3113.5	1 mg
62279	Bim BH3 (81-106), human, Peptide IV	DMRPEIWIAQELRRIGDEFNAYYARR	3269.7	1 mg
62265	Bim BH3 (82-100), human, TAMRA-labeled	5-TAMRA-MRPEIWIAQELRRIGDEFNA	2857.3	0.5 mg
62484	Bim BH3 (87-100)	WIAQELRRIGDEFN	1747	1 mg
62263	Bim-23056	fFYwKVFnal-NH ₂	1232.5	1 mg
62281	BMf-BH3	LQHRAEVQIARKLQCIADQFHLHT	3012.5	1 mg
62282	Noxa BH3, Peptide 1	PAELEVECATQLRRFGDKLNFRQKLL	3075.6	1 mg
62403	Noxa BH3, Peptide 2	VPADLKDECAQLRRIGDKVNLRQKL	2879.4	1 mg
62323	NuBCP-9, Nur77-derived Bcl-2 Converting Peptide	FSRSLHSLL	1059.2	1 mg
62404	PUMA BH3	EEQWAREIGAQLRRMADDLNAQYER	3049.3	1 mg

Bradykinins

62162	Bradykinin, R9C	RPPGFSPFC	1007.2	1 mg
63312	Hemopressin	PVNFKLLSH	1054.3	1 mg

Cancer Research Peptides

61371	234 CM, Codon 234 Point Mutation, p53 Product	KYICNSSCM	1048.3	1 mg
61372	234 CW, Codon 234, p53 Product	KYMCNSSCM	1066.3	1 mg
61271	[-2] pPSA, Prostate-Specific Antigen, truncated	SRIVGGWECEK	1263.4	1 mg
61273	[-4] pPSA, pro-Prostate-Specific Antigen	ILSRIVGGWECEK	1489.8	1 mg
62926	53BP2 (490-498), p53-Binding Loop (CDB3)	REDEDEIEW	1220.2	1 mg
61443	Adipophilin	SVASTITGV	833.9	1 mg
61323	Amphoterin (150-183)	KLKEKEYEKDIAAYRAKGKPDAAKKGVVKAEK-SKK	3805.5	1 mg
61368	anti-Fit1, VEGF Related	GNQWFI	763.9	1 mg
61272	[APLILSR] pPSA, pro-Prostate-Specific Antigen	APLILSRIVGGWECEK	1771.1	1 mg
61502	BAGE (2-10)	AARAVFLAL	931.2	1 mg
62536	BRCA1 (610-619)	SSKKQKRSHK	1213.4	1 mg
61793	Cathepsin D and E Substrate	Mca-GKPILFFRLK(Dnp)-r-NH ₂	1756.1	1 mg
62368	Cathepsin K Substrate	Abz-HPGGPQ-EDDnp	920	0.1 mg
61861-05	[Cys(Acm) ³³]-Endostatin (6-49)	FQPVLHLVALNSPLSGGMRGIRGADFQ-C(Acm)-FQQARAVGLAGTFRAF	4777.6	0.5 mg
62616	Cys-p21 (139-154)	CGRKRRQTSMTDFYHSK	2101.4	1 mg
62013	EGFR (662-682)	RRELVEPLTPSGEAPNQALLR	2346.7	1 mg
62695	F4.2, Gastric Signet Ring Cell Carcinoma Derived Peptide	YSWMDISWVI	1303.5	1 mg
61274	G154, gp100 (154-162)	KTWGQYWQV	1195.4	1 mg
61276	G209, gp100 (209-217)	ITDQVPFSV	1005.1	1 mg
61277	G209-2M, gp100 (209-217)	IMDQVPFSV	1035.2	1 mg
61280	G280-9, gp100 (280-288)-Lys(Biotin)	YLEPGPVTA-K(Biotin)	946.1	1 mg
61278	G280-9V, gp100(280-288)-Lys(Biotin)	YLEPGPVTV-K(Biotin)	974.1	1 mg
62589	gp100 (25-33), human	KVPRNQDWL	1155.3	1 mg
61505	gp100 (177-186)	AMLGTHTMEV	1089.3	1 mg
61487	gp100 (178-186)	MLGTHTMEV	1018.2	1 mg
61466	gp100 (457-466)	LLDGTATLRL	1072.3	1 mg
61482	gp100 (476-485)	VLYRYGSFSV	1190.4	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61481	gp100 (570-579)	SLADTNSLAV	990.1	1 mg
61478	gp100 (614-622)	LIYRRRLMK	1248.6	1 mg
61480	gp100 (619-627)	RLMKQDFSV	1123.4	1 mg
61479	gp100 (639-647)	RLPRIFCSC	1094.4	1 mg
61463	GPC3, Glypican-3 (144-152)	FVGEFFTDV	1060.2	1 mg
61464	GPC3, Glypican-3 (298-306), mouse	EYILSLEEL	1108.3	1 mg
62205	GRP78 Binding Chimeric Peptide Motif	WIFPWQLGGklaklaklaklak-NH ₂	2721.5	1 mg
62511	HB-1, B-cell Leukemia-Associated Antigen (26-41)	EDDVYLRHSSSLTYRL	1954.1	1 mg
62512	HB-1 (18-41)	WKSELVEVDVYLRHSSSLTYRL	2796.1	1 mg
62011	HPV Oncoprotein, E7 (43-62)	GQAEPDRAHYNIVTFCKCD	2270.6	1 mg
62012	HPV Oncoprotein, E7 (43-77)	GQAEPDRAHYNIVTFCKCDSTLRLCVQSTH-VDIR	3980.6	1 mg
62941	HPV16 E7(86-93)	TLGIVCPI	815.1	1 mg
62401	IL-8 Inhibitor	Ac-RRWWCR-NH ₂	1003.2	1 mg
62204	IL-11R _a Binding Peptide	CGRRAGGSC (Disulfide bridge: 1-9)	864.0	1 mg
61655	iLRP, Immature Laminin Receptor Protein (58-66)	LLLAARAIIV	939.2	1 mg
61656	iLRP2, Immature Laminin Receptor Protein (60-68)	LAARAIVAI	897.1	1 mg
61654	iLRP3, Immature Laminin Receptor Protein (146-154)	ALCNTDSPL	933.1	1 mg
61657	iLRP4, Immature Laminin Receptor Protein (7-15)	VLQMKEEDV	1090.3	1 mg
61380	IP3 peptide, MDM-2 Binding Motif, Biotin-labeled	Ac-MPRFMDYWEGLNKK-Biotin	2083.5	1 mg
61379	IP3 peptide, MDM-2 Binding Motif -Lys(TAMRA)	MPRFMDYWEGLN-K(TAMRA)	2099.5	1 mg
61378	IP3 Truncated Peptide, MDM-2 Binding Motif, acetylated	Ac-FMDYWEGLN	1216.3	1 mg
61507	KM-HN-1, Cancer/Testis Antigen (65-74)	SFQALRMQTL	1194.4	1 mg
61508	KM-HN-1, Cancer/Testis Antigen (107-116)	VFGTRIEKDL	1177.4	1 mg
61506	KM-HN-1, Cancer/Testis Antigen (196-204)	NYNNFYRFL	1250.4	1 mg
61509	KM-HN-1, Cancer/Testis Antigen (335-343)	HFCRKCKKL	1162.5	1 mg
61510	KM-HN-1, Cancer/Testis Antigen (499-508)	EYSKECLKEF	1275.5	1 mg
61511	KM-HN-1, Cancer/Testis Antigen (770-778)	EYLSLSDKI	1067.2	1 mg
61455	Livin7 / ML-IAP, Melanoma Inhibitor of Apoptosis Protein	KWFPSQCFL	1268.6	1 mg
62169	LyP-1, Lymphatic Tumor Antigen, cyclized	CGNKRTRGC (Disulfide bridge: 1-9)	992.2	1 mg
62170	LyP-1, Lymphatic Tumor Antigen, linear	CGNKRTRGC	994.2	1 mg
62710	Malaria Circumsporozoite Protein, CSP (334-342)	YLKKIKNSL	1106.4	1 mg
62576	Maspin Reactive Site Loop (RSL), (330-345)	GGDSIEVPGARILQHK	1676.9	1 mg
62575	[Gln ³⁴⁰]-Maspin, Reactive Site Loop (RSL), (330-345)	GGDSIEVPGAQILQHK	1648.8	1 mg
61358	MUM-1, Melanoma Associated Antigen, mutated (261-269)	EEKLIVVLF	1089.4	1 mg
61470	GnT-V, (nt 38-47) Melanoma Antigen	VLPDVFIRCV	1160.5	1 mg
61471	MZ2-F, HLA-Cw6-Restricted Epitope	YRPRPRRY	1163.4	1 mg
62712	NY-ESO-1 (53-62)	ASGPGGGAPR	825.9	1 mg
61461	ORF2, RACE, Renal Tumor Antigen Fragment	SPSSNRIRNT	1131.2	1 mg
61462	ORF5, RACE, Renal Tumor Antigen Fragment	PASKKTDPQK	1099.3	1 mg
62614	p21 (140-147)	RKRRQTSM	1062.3	1 mg
62613	[pThr ¹⁴⁵]-p21 (140-147)	RKRRQ-pT-SM	1142.3	1 mg
62852	Pirh2-Derived Peptide (CDB62) (120-137)	LKC�LCLTTNLRGKHKCI	2058.6	1 mg
62105	Proapoptotic Chimera Prostate Vasculature Marker	SMSIARLGGklaklaklaklak-NH ₂	2396.1	1 mg
62106	Prostate Vasculature Marker, Lys(Biotin)	SMSIARLK(LC-Biotin)	1244.6	1 mg
61459	PSA 1, Prostate-Specific Antigen 1 (141-150)	FLTPKKLQCV	1176.5	1 mg
61460	PSA 2, Prostate-Specific Antigen 2 (146-154)	KLQCVDLHV	1054.3	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62851	Rad51 (175-190) (CDB55)	AERYGLSGSDVLDNVA	1665.8	1 mg
62850	Rad51 (179-190)	GLSGSDVLDNVA	1146.2	1 mg
61297	Stromal Target Antigen	SIYYRYRGL	1197.4	1 mg
61447	Survivin	ELTLGEFLKL	1162.4	1 mg
62691	Survivin (85-93)	AFLSVKKQF	1067.3	1 mg
62693	Survivin 2B (80-88)	AYACNTSTL	943.1	1 mg
62413	Telomerase Reverse Transcriptase p572Y (TERT572Y)	RLFFYRKS	1215.5	1 mg
61475	TRP-1, Tyrosinase-Related Protein-1 Fragment	MSLQRQFL	1022.2	1 mg
61474	TRP-2, Tyrosinase-Related Protein-2 Coding Region Fragment	LLPGGRPYR	1028.2	1 mg
62090	VEGFR-2/KDR I, murine	FSNSTNDILI	1123.2	1 mg
62091	VEGFR-2/KDR 2, murine	VILTNPISM	987.2	1 mg
62621	WP9QY, TNF- α Antagonist	YCWSQYLCY (Disulfide bridge: 2-8)	1226.4	1 mg

Cardiac Peptides

62911	Cardiac Troponin I (hcTnI) (39-58), human	SKISASRKLQLKTLQLQIAK	2239.8	1 mg
63775	Cardiac Troponin I (41-49)	SASRKLQLK	1030.2	1 mg

CEF Control Peptide Pools

61036-003	CEF Control Peptide Pool		N/A	1 mg
62339	Cytomegalovirus (CMV) Control Peptide Pool	CEF20,21,22,23,33	N/A	1.25 mg
62341	Epstein-Barr Virus (EBV) Control Peptide Pool	CEF10,11,12,14,15,16,17,18,19,27,28,29,30,31,32	N/A	3.75 mg
62340	Influenza Virus Control Peptide Pool	CEF1,2,3,4,5,6,7,8,9,24,25,26	N/A	3 mg

Cell Adhesion Peptides

62639	37, 40 GAP26, Connexin Mimetic	VCYDQAFPISHIR	1548.8	1 mg
62641	40Gap 27, Connexin Mimetic	SRPTEKNVFIV	1289.5	1 mg
62644	43Gap 26, Connexin Mimetic	VCYDKSFPISHVR	1550.8	1 mg
62643	43Gap 36, Connexin Mimetic	KRDPCHQVDCFLSRPTEK	2159.5	1 mg
62411	α B-Crystallin (73-92)	DRFSVNLVDVKHFSPEELKVK	2387.7	1 mg
62697	Binin (103-120), <i>Strongylocentrotus purpuratus</i>	LGLLLRHLRHHSNLLANI	2090.5	1 mg
62842	Calreticulin (CRT) Binding Peptide 1	GQPMYQQPMY	1171.4	1 mg
62843	Calreticulin (CRT) Binding Peptide 1, Biotin-labeled	Biotin-GQPMYQQPMY	1397.7	1 mg
62622	Hyaluronan Inhibitor	GAHWQFNALTVR	1399.6	1 mg
61664	L1CD Cell Adhesion Molecule (1144-1163)	KRSKGGKYSVKDKEDTQVDS	2255.5	1 mg
61660	L1FLCD (1173-1185), L1 Cell Adhesion Molecule, Cytoplasmic Domain	FGEYRSLESDNEE	1574.6	1 mg
61661	L1ARSLCD, [Ala ¹¹⁷⁶]-L1FLCD (1173-1185), L1 Cell Adhesion Molecule, Cytoplasmic Domain	FGEARSLESDNEE	1482.5	1 mg
61663	L1CD (1163-1176), L1 FL Cell Adhesion Molecule, Cytoplasmic Domain	SEARPMKDETFGEY	1659.8	1 mg
61662	L1 δ SLECD, L1 Cell Adhesion Molecule, Cytoplasmic Domain Mutant	FGEYSDNEE	1089.0	1 mg
62114	pALA, Polyalanine Peptide	AAADAAAAL	743.8	1 mg
62501	S1P, Sphingosine 1-Phosphate Peptide	VSTSIPEVKALRSSVSDYGNVDIIVRHNYTG-KLNIGAEEKDHGIK	5009.6	0.5 mg

Cell Permeable / Drug Delivery Peptides

61376	Antennapedia Leader Peptide (CT)	KKWKMRRNQFWVKVQRG	2275.8	1 mg
-------	----------------------------------	-------------------	--------	------

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61255	Buforin	TRSSRAGLQFPVGRVHRLLRK	2434.9	1 mg
62565	Chimeric Rabies Virus Glycoprotein Fragment (RVG-9R)	YTIWMPENPRPGTPCDIFTNSRG-KRASNGGGGRRRRRRRRR	4843.5	1 mg
62005	[Cys ⁵⁸]-105Y, Cell Penetrating Peptide, α1-antitrypsin (358-374)	CSIPPEVKFNKPFVYLI	1994.4	1 mg
62397	Lipid Membrane Translocating Peptide	KKAAAVLLPVLLAAP	1474.9	1 mg
62398	Lipid Membrane Translocating Peptide, D-isomer	kkaaavllpvllaap	1474.9	1 mg
62646	Mastoparan 7	INLKALAALAKALL-NH ₂	1421.8	1 mg
62647	Mastoparan X	INWKGIAAMAKKLL-NH ₂	1556	1 mg
61257	Nuclear Localization Signal Peptide	PKKKRKV	883.2	1 mg
63849	Pep-1-Cysteamine	Ac-KETWWETWWTEWSQPKKKKRKV-cyste-amine	2950.3	1 mg
62566	Rabies Virus Glycoprotein (RVG)	YTIWMPENPRPGTPCDIFTNSRGKRASNG	3266.7	1 mg
62564	Rabies Virus Matrix Protein Fragment (RV-MAT)	MNLLRKIVKNRREDEDTQKSSPASAPLDDG	3255.7	1 mg
61254	SynB1, Cell Penetration Peptide	RGGRLSYSRRRFSTSTGRA	2171.4	1 mg
62056	TAT (47-57)-Lys(TAMRA)	YGRKKRRQRRRK(TAMRA)	2100.5	1 mg
62064	TAT (48-57)	GRKKRRQRRR	1396.7	1 mg
62063	Cys-TAT (48-57)	CGRKKRRQRRR	1499.8	1 mg
62042	TAT-NR2Bct	YGRKKRRQRRRKLSSIESDV	2518.9	1 mg
62237	TAT-NSF81scr Fusion Polypeptide, scrambled	YGRKKRRQRRR-GGG-QDGCKYFAT-DETIMKLSIAI	3960.6	1 mg
62240	TAT-NSF222 Fusion Peptide	YGRKKRRQRRR-GGG-LDKEFNSIFRRRAFAS-RVFPPE	4239.9	1 mg
62210	TAT-NSF222scr Fusion Polypeptide, scrambled	YGRKKRRQRRR-GGG-ENSFRFLADIFPAK-AFPVRF	4214.9	1 mg
62209	TAT-NSF700scr Fusion Peptide, scrambled	YGRKKRRQRRR-GGG-IPPVYFSRLDLNLV-VLLLAQL	4109.9	1 mg
62238	TAT-NSF700 Fusion Peptide	YGRKKRRQRRR-GGG-LLDYVPIGPRFSNLV-LQALLV	4167	1 mg
62066	Transdermal Peptide	ACSSSPSKHCG	1063.2	1 mg
61256	Transportan	GWTLNSAGYLLGKINLKALAALAKKIL	2841.5	1 mg
62004	105Y, Cell Penetrating Peptide, α1-Antitrypsin (359-374)	SIPPEVKFNKPFVYLI	1891.3	1 mg

Chaperone Peptides

62872	α-A-Crystallin (70-88)	KFVIFLDVKHFSPEDLTVK	2262.7	1 mg
63724	Hsc70-binding Peptide II	NIVRKKK	885.1	1 mg

Chemotaxis Peptides

62009	Ac-PGP, Neutrophil Chemoattractant Tripeptide	Ac-PGP	311.3	1 mg
62719	Formyl-MLF Peptide	Formyl-MLF	437.6	1 mg
62401	IL-8 Inhibitor	Ac-RRWWCR-NH ₂	1003.2	1 mg
62519	MMK-1	LESIFRSLLFRVM	1611	1 mg
62520	MMK-1, amide	LESIFRSLLFRVM-NH ₂	1610	1 mg
62863	Ropalidian Chemotactic Peptide	IVPFLGPLLGLLT-NH ₂	1351.7	1 mg
62864	Vespid Chemotactic Peptide A	FLPMIAKLLGGLL-NH ₂	1384.8	1 mg
62861	Vespid Chemotactic Peptide L	FLPIIAKLVSGLL-NH ₂	1382.8	1 mg
62865	Vespid Chemotactic Peptide M	FLPIIGKLLSGLL-NH ₂	1382.8	1 mg
62869	Vespid Chemotactic Peptide T	FLPILGKILGGLL-NH ₂	1352.8	1 mg
62862	Vespid Chemotactic Peptide X	FLPIIAKLLGGLL-NH ₂	1366.8	1 mg
62638	WRW4, Formyl Peptide Receptor-Like 1 (FPRL1) Antagonist	WRWVVVV	1105.3	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
ClearPoint™-Heavy Isotope Labeled Peptides				
63737	ClearPoint™ β-Amyloid (1-40), ¹³ C, ¹⁵ N-labeled at Arg & Lys	DAEF-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVV [R*=R(U- ¹³ C ₆ , U- ¹⁵ N ₄) & K*=K(U- ¹³ C ₆ , U- ¹⁵ N ₂)]	4355.9	50 µg
63740	ClearPoint™ β-Amyloid (1-40), U- ¹³ C-labeled at Leu ¹⁷	DAEFRHDSGYEVHHQK-L*-VFFAEDVG-SNKGAIIGLMVGGVV [L*=L(U- ¹³ C ₆)]	4335.9	50 µg
63738	ClearPoint™ β-Amyloid (1-42), ¹³ C, ¹⁵ N-labeled at Arg & Lys	DAEF-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVVIA [R*=R(U- ¹³ C ₆ , U- ¹⁵ N ₄) & K*=K(U- ¹³ C ₆ , U- ¹⁵ N ₂)]	4540.1	50 µg
63739	ClearPoint™ β-Amyloid (4-42), ¹³ C, ¹⁵ N-labeled at Arg & Lys	F-R*-HDSGYEVHHQ-K*-LVFFAEDVGSN-K*-GAIIGLMVGGVVIA [R*=R(U- ¹³ C ₆ , U- ¹⁵ N ₄) & K*=K(U- ¹³ C ₆ , U- ¹⁵ N ₂)]	4224.8	50 µg
63741	ClearPoint™ β-Amyloid (4-42), U- ¹³ C, ¹⁵ N-labeled at Val ¹²	FRHDSGYE-V*-HHQKLVFFAEDVG-SNKGAIIGLMVGGVVIA [V*=V(U- ¹³ C ₅ , U- ¹⁵ N)]	4204.8	50 µg
63742	ClearPoint™ β-Amyloid (8-42), U- ¹³ C, ¹⁵ N-labeled at Val ¹²	SGYE-V*-HHQKLVFFAEDVGSNKGAIIGLMVGGVVIA [V*=V(U- ¹³ C ₅ , U- ¹⁵ N)]	3649.2	50 µg
61182	ClearPoint™ Angiotensin I, human, ¹³ C and ¹⁵ N-labeled	DRVY-I*-HPFHL [I*=I(U- ¹³ C ₆ , ¹⁵ N)]	1303.5	0.1 mg
61220	ClearPoint™ BSA (347-359), Isotopic labeled, Mass Spec Standard	DAF-L*-GSF-L*-YEYSR [L* = L(U- ¹³ C ₆ , ¹⁵ N)]	1581.7	0.1 mg
61219	ClearPoint™ C-Reactive Protein-Like Peptide, Isotopic labeled, Mass Spec Standard	GYSI-F*-SYATK [F* = F(U- ¹³ C ₉ , ¹⁵ N)]	1146.3	0.1 mg
Cytochromes and Related Peptides				
62751	Cytochrome b (150-172), Large Subunit, L123	SYLNFARKRIKNIPEGGLYLAVTL	2624.1	1 mg
62752	Cytochrome b (175-194), Small Subunit	AGGPPGGPQVNPIPVTDDEVV	1900.1	1 mg
62750	Cytochrome b (550-569), Large Subunit, Li Peptide	ISNSESGRPGVHFIFNKENF	2279.5	1 mg
62588	Cytochrome P450 4A (431-445), rat	DPSRFAPDSPRHSHS	1692.8	1 mg
Dengue Viral Peptides				
61366	2A/2B Dengue Protease Substrate	Ac-RTSKKR-pNA	937.1	1 mg
61365	2B/3, Dengue Protease Substrate	Ac-EVKKQR-pNA	949.1	1 mg
61364	3/4A, Dengue Protease Substrate	Ac-FAAGRK-pNA	810.9	1 mg
62032	Dengue 1 Fragment	VGIGNRDFV	976.1	1 mg
62033	Dengue 2 Fragment	IGISNRDFV	1020.2	1 mg
Dynorphins				
62682	Deltorphin A	YmFHLMD-NH ₂	955.2	1 mg
62683	Deltorphin B	YaFEVVG-NH ₂	782.9	1 mg
62688	[Dmt ¹]-Deltorphin B	(Dmt)-aFEVVG-NH ₂	810.9	1 mg
62681	Deltorphin C	YaFDVVG-NH ₂	768.9	1 mg
62680	[dAla ² , Cys ⁴]-Deltorphin C	YaFCVVG-NH ₂	756.9	1 mg
62684	Dermorphin	YaFGYPS	803.9	1 mg
62687	Dynorphin B (1-9)	YGGFLRRQF	1143.3	1 mg
62686	Dynorphin B (1-13)	YGGFLRRQFKVVT	1570.9	1 mg
62788	Dynorphin B-29, Leumorphin	YGGFLRRQFKVVTRSQEDPNAYYEELFDV	3527.9	1 mg
Endorphins				
62678	α-Neo-Endorphin	YGGFLRKYPK	1228.5	1 mg
62669	Neoendorphin (166-174)	YGGFLRKYP	1100.3	1 mg
Endothelins				
62946-05	Big Endothelin-1 (1-38), human	CSCSSLMDKECVYFCHLDIIWVNTPEHVVPY-GLGSPRS (Disulfide bridge: 1-15 and 3-11)	4283	0.5 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62946-1	Big Endothelin-1 (1-38), human	CSCSSSLMDKECVYFCHLDIIWVNTPEHVVPY-GLGSPRS (Disulfide bridge: 1-15 and 3-11)	4283	1 mg
62357	Endothelin Antagonist, BQ123, Cyclo(-wdPvL-)	Cyclo(-wdPvL-)	610.7	1 mg
62356	Endothelin-2, mouse	CSCSNWLDKECVYFCHLDIIW (Disulfide bridge: 1-15 and 3-11)	2574	0.5 mg

Enkephalins

62665	[Met ⁶]-Enkephalin-Arg-Phe	YGGFMRF	877	1 mg
62666	Proenkephalin A Precursor (197-208)	SPQLEDEAKEELQ	1386.5	1 mg
62723	Proenkephalin A Precursor (219-229)	VGRPEWWMDYQ	1466.7	1 mg
62728	QF-ERP7, FRET	Abz-GGFLRRV-EDDnp	1132.4	0.5 mg

Enzyme Substrates and Inhibitors

61366	2A/2B Dengue Protease Substrate	Ac-RTSKKR-pNA	937.1	1 mg
61365	2B/3, Dengue Protease Substrate	Ac-EVKKQR-pNA	949.1	1 mg
61364	3/4A, Dengue Protease Substrate	Ac-FAAGRK-pNA	810.9	1 mg
62324	2 CDC25C, CHK1 and CHK2 Substrate, phosphorylated	KKKVSRSGLYR-pS-PSMPENLNRRP-NH ₂	2780.2	1 mg
61862	Activated Protein C (390-404), human	YGVYTKVSRYLWIIH	1900.2	1 mg
62984	APCE, Antiplasmin-Cleaving Enzyme (1-15)	IVLRPSRVHNSEENT	1750.9	1 mg
61318	C1-Inhibitor (Abz/Dnp)	Abz-SVARTLLV-K(Dnp)-NH ₂	1270.5	1 mg
63860	CaMKK Peptide Substrate	AKPKGNKDYHLQTCGSLAYRRR	2666.1	1 mg
61859	Cathepsin S Substrate	Ac-KQKLR-AMC	871.1	1 mg
61898	Dok-4, Downstream of Tyrosine Kinase-4 (130-145)	GVQCEQTDTRFNVFLLP	1866.1	1 mg
61893	Dok-4, Downstream of Tyrosine Kinase-4 (263-275)	LPRSAYWHHITGS	1524.7	1 mg
61895	Dok-5, Downstream of Tyrosine Kinase-5 (130-145)	GVEREQSERFNVYLMP	1954.2	1 mg
61892	Dok-5, Downstream of Tyrosine Kinase-5 (263-275)	LPRSAYWQHITRQ	1655.9	1 mg
61894	Dok-6, Downstream of Tyrosine Kinase-6 (130-145)	GVQREQNERFNVYLMP	1980.2	1 mg
61882	Dok-6, Downstream of Tyrosine Kinase-6 (263-275)	LPRSAYWHHITRQ	1664.9	1 mg
62974	Enterokinase Substrate	GDDDDK-βNA	788.8	1 mg
63854	gp91 ds-tat Peptide 2	RKKRRQRRRCSTRIRRQL-NH ₂	2453.0	1 mg
63855	gp91 ds-tat Peptide 2, sgp91 ds-tat Peptide 2, scrambled	RKKRRQRRRCSTRIRRQL-NH ₂	2453.0	1 mg
61370	HJ Inhibitor Peptide 2	KWWCRW	964.2	1 mg
62990	Matriptase I Substrate II	Boc-AGPR-AMC	656.7	1 mg
61668	NubO (68-75), Ndufa4 (68-75)	VNVDSYSL	937.1	1 mg
63726	O-linked GlcNAc transferase (OGT) Substrate	KKKYPPGGSTPVSSANMM	1783.1	1 mg
63736-01	Peptide S4 (A FRET peptide for O-linked Glycosylation Assay)	FITC-Aha-STPVSRANM-K(QXL™ 520)-R-K(QXL™ 520); Aha=LC=6-Aminohexanoyl	2860.1	0.1 mg
63736-1	Peptide S4 (A FRET peptide for O-linked Glycosylation Assay)	FITC-Aha-STPVSRANM-K(QXL™ 520)-R-K(QXL™ 520); Aha=LC=6-Aminohexanoyl	2860.1	1 mg
61632	proFIX18, pro-Factor IX (-18 - -1)	TVFLDHENANKILNRPKR	2165.5	1 mg
61633	proFIX28, pro-Factor IX (-18 - +10)	TVFLDHENANKILNRPKRYNSGKLEEFV	332.8	1 mg
62113	Proprotein Convertase Substrate, PC4 Substrate	Ac-KTKQLR-AMC	972.1	1 mg
61634	proPT18, pro-Prothrombin (-18 - -1)	HVFLAPQQARSLLRVRR	2175.6	1 mg
61635	proPT28, pro-Prothrombin (-18 - +10)	HVFLAPQQARSLLRVRRANTFLEEVRK	3363.9	1 mg
62051	Transglutaminase Substrate, Biotin-labeled	Biotin-TVQQEL	943.1	1 mg

Exendins

63897	Exendin (4-39)	GTFTSDLSKQMEEEAVRLFIEWLKNGGPSS-GAPPPS-NH ₂	3863.3	1 mg
-------	----------------	---	--------	------

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
63898	Exendin (5-39)	TFTSDLSKQMEEEEAVRLFIEWLKNGGPSS-GAPPPS-NH ₂	3806.3	1 mg
63902	Exendin (7-39)	TSDLSKQMEEEEAVRLFIEWLKNGGPSS-GAPPPS-NH ₂	3558	1 mg
63903	Exendin (10-39)	LSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH ₂	3254.7	1 mg
62072	Exendin 4 Fragment, Trp cage	NLYIQWLKDGPPSSGRPPPS	2169.4	1 mg
63714	[Cys(HiLyte Fluor™ 647 C2 maleimide)]-Exendin-4	C(HiLyteFluor 647 C2 maleimide)-HGEGTFTS-DLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH ₂	5486.3	50 ug
63900	[Des-His ¹ , Glu ⁸]-Exendin-4	GEGTPTSELSKQMEEEEAVRLPIEWLKNGG-PSSGAPPPS-NH ₂	3963.4	1 mg
63899	Fluorescein-Trp ²⁵ -Exendin-4 (FLEX)	HGEGTFTSDLSKQMEEEEAVRLFIE-(Trp-S-FAM)-LKNGGPSSGAPPPS-NH ₂	4606.4	0.5 mg
63901	Glicentin-Related Peptide (21-50), human	RSLQDTEEKSRFSASQADPLSDPDQMNE-NH ₂	3383.5	1 mg
61867-05	Helodermin	HSDAIFTQQYSKLLAKLALQKYLASILGSRT-SPPP	3844.5	0.5 mg

Fibrinogen and Related Peptides

62508	Elastin-Like Octapeptide	GVGVPGVGVPGVGVPGVG	1459.7	1 mg
62128	γ-Fibrinogen (377-395)	YSMKETTMMKIIPFNRLSIG	2229.7	1 mg
62129	γ-Fibrinogen (377-395), scrambled	KMMISYTFPIERTGLISNK	2229.7	1 mg
62704	Heparin-Binding Peptide V	WQPPRARI	1023.2	1 mg
62349	Integrin Binding Peptide	Ac-GCGYGRGDSPG-NH ₂	1066.1	1 mg
62338	TP508, Thrombin-derived Peptide	AGYKPDEGKRGDACEGDSGGPFV	2312.5	1 mg

Fibronectin Fragments

62352	Cyclo (-RADfE), RGD negative control	Cyclo(-RADfE-)	618.7	1 mg
62351	Cyclo (-RADfK-), RGD negative control	Cyclo(-RADfK-)	617.7	1 mg
62333-01	Cyclo[-RGDyK(HiLyte Fluor™ 750)]	Cyclo[-RGDy-K(HiLyte Fluor™ 750)]	1630.5	0.1 mg
62333-1	Cyclo[-RGDyK(HiLyte Fluor™ 750)]	Cyclo[-RGDy-K(HiLyte Fluor™ 750)]	1630.5	1 mg
62581	Fibronectin (1906-1924)	YEKPGSPPREVPRPRPGV	2117.5	1 mg
62578	Fibronectin (1946-1960), FN-C/H2	KNNQKSEPLIGRKKT	1741	1 mg
62580	Fibronectin (1951-1960), FN-C/H2-COOH	SEPLIGRKKT	1128.3	1 mg
62579	Fibronectin (1954-1960)	LIGRKKT	815	1 mg
61384	Fibrinopeptide B, Bovine	QFPTYDEGQDDRPKVGLGAR	2364.5	1 mg
62350	GGGGRGDS	GGGGRGDS	661.6	1 mg
62525	GRGDK, FAM-labeled	GRGDK(5-FAM)	889.9	1 mg
62347	GRGDS, Biotin-LC-labeled	Biotin-LC-GRGDS	830	1 mg
62705	Heparin-Binding Peptide V, scrambled	RPQIPWAR	1023.2	1 mg
62527	[Glu ³]-RGES, Control for RGD Peptides	RGES	447.5	1 mg
62982	THS, Fibronectin Type II Ligand	THSHQWRHHQFPAPT	1867	1 mg
62700	V14 Peptide	VVGSPSAQDEASPL	1356.5	1 mg
62980	WHW, Fibronectin Type II Ligand	WHWRHRIPLQLAAGR	1897.2	1 mg
62654	YRGDS Fibronectin Fragment	YRGDS	596.6	1 mg

FRET Peptides

62231	Bacterial Sortase Substrate I, FRET	DABCYL-LPETG-EDANS	1015.2	1 mg
61318	C1-Inhibitor (Abz/Dnp)	Abz-SVARTLLV-K(Dnp)-NH ₂	1270.5	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61793	Cathepsin D and E Substrate	Mca-GKPILFFRLK(Dnp)-r-NH ₂	1756.1	1 mg
62368	Cathepsin K Substrate	Abz-HPGGPQ-EDDnp	920	0.1 mg
61859	Cathepsin S Substrate	Ac-KQKLR-AMC	871.1	1 mg
61314	Cls Substrate C2, Second Complement Component (5-FAM/QXL™520), amide	5-FAM-SLGRKIQI-K(QXL™ 520)-NH ₂	2019.2	0.5 mg
61315	Cls Substrate C2, Second Complement Component (Abz/Dnp), amide	Abz-SLGRKIQI-K(Dnp)-NH ₂	1326.5	1 mg
61317	Cls Substrate C2, Second Complement Component (DABCYL/EDANS)	DABCYL-SLGRKIQI-EDANS	1413.7	1 mg
61316	Cls Substrate C4, Fourth Complement Component (Abz/Dnp), amide	Abz-GLQRALEI-Lys(Dnp)-NH ₂	1311.5	1 mg
63798	Endothelin Converting Enzyme, ECE-1 FRET Substrate	Mca-RPPGFSAF-K(Dnp)	1388.5	0.5 mg
62015	Hemoglobin, 2837a	DABCYL-GABA-ALERMFLSFP-EDANS	1795.2	1 mg
62014	Hemoglobin, 3037a, Malaria FRET Substrate II	DABCYL-GABA-ERMFLSFP-EDANS	1610.9	1 mg
62050	Hemoglobin Fragment, Plasmeprin II (PM II) FRET, Substrate, 2837b	EDANS-CO-CH ₂ -CH ₂ -CO-ALERMFLSFP-Dap(DABCYL)	1896.2	1 mg
62166	Interferon α A (101-110), FRET Peptide	GVGVTETPLME(EDANS)GVGVTETPLM-K(DABCYL)K	2874.4	1 mg
63736-01	Peptide S4 (A FRET peptide for O-linked Glycosylation Assay)	FITC-Aha-STPVSRANM-K(QXL™ 520)-R-K(QXL™ 520); Aha=LC=6-Aminohexanoyl	2860.1	0.1 mg
63736-1	Peptide S4 (A FRET peptide for O-linked Glycosylation Assay)	FITC-Aha-STPVSRANM-K(QXL™ 520)-R-K(QXL™ 520); Aha=LC=6-Aminohexanoyl	2860.1	1 mg
61872-01	Peptide Substrate for Renin 520 Assay kit		2000 - 2200	0.1 mg
61872	Peptide Substrate for Renin 520 Assay kit		2000 - 2200	1 mg
62334-01	Rat Renin FRET Substrate (5-FAM/QXL™520)		2000 - 2200	0.1 mg
62334	Rat Renin FRET Substrate (5-FAM/QXL™520)		2000 - 2200	1 mg
62022	Renin Substrate 1, FRET	R-E(EDANS)-IHPFHLVIHT-K(DABCYL)-R	2282.7	1 mg
61618	SARS-CoV PLP2 Optimal Substrate (813-823)	FRLKGGAPIKGV	1139.0	1 mg
61619	SARS-CoV PLP2 Optimal Substrate (813-823), FRET	Abz-FRLKGGAPIKGV-EDDnp	1569.9	1 mg
63732-01	Tat FRET Peptide for RNA Binding Assays	FITC-AAARKKRRQRRRAAAC(TAMRA-Mal)-NH ₂		0.1 mg
63732-05	Tat FRET Peptide for RNA Binding Assays	FITC-AAARKKRRQRRRAAAC(TAMRA-Mal)-NH ₂		0.5 mg

Galanins

61845-05	Galanin Like Peptide (GALP), human	APAHRGRGGWTLNSAGYLLGPVLHLPQMG-DQDGKRETALEILDWLKIDGLPYSHPPQPS	6500.4	0.5 mg
61846	Galanin (1-13) / Galanin Like Peptide (GALP) (9-21), human	GWTLNSAGYLLGP	1348.5	1 mg
61848	Galanin Like Peptide (GALP) (41-60), human, porcine, rat	ILDWLKIDGLPYSRSPRMT	2332.8	1 mg
61850-05	Galanin Like Peptide (GALP), N-terminal fragment, porcine	APVHRGRGGWTLNSAGYLLGPVLHPPSRAE	3166.6	1 mg
61849-05	Galanin Like Peptide (GALP), porcine	APVHRGRGGWTLNSAGYLLGPVLHPPSRAE-GGGKGTALGILDWLKIDGLPYPSQLAS	6204.1	0.5 mg
61847-05	Galanin Like Peptide (GALP), rat	APAHRGRGGWTLNSAGYLLGPVLHLLSSKAN-QGRKTDSEILDWLKIDGLPYSRSPRMT	6502.4	0.5 mg

Gastrins

61745	Gastrin derived peptide	GPWLEEEEEAYGWMDFK-NH ₂	2115.3	1 mg
61746	Gastrin derived peptide, FAM-labeled	5-FAM-GPWLEEEEEAYGWMDFK-NH ₂	2473.6	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62095	Gastrin-releasing Peptide 10 / Neuromedin C (GRP10)	GNHWAVGHLM-NH ₂	1120.3	1 mg
62096	Gastrin-releasing Peptide (1-17)	VPLPAGGGTVLTKMYPR	1757.1	1 mg
62558	Leptin (22-56), human	VPIQKVQDDTKTLIKTIVTRINDISHTQS-VSSKQK	3950.6	1 mg
62521	Leptin (57-74)	VTGLDFIPGLHPILTLSK	1921.3	1 mg
62853	Leptin (93-105), human	NVIQISNDLENLR	1527.7	1 mg
62559	Leptin (116-130), mouse	SCSLPQTSGLQKPES	1561.7	1 mg

Ghrelin and Obestatin

61714	Biotin-Ghrelin, bovine	Biotin-GS-S(n-octanoylated)-FLSPE-HQKLQRKEAKKPSGRLKPR	3442.8	1 mg
61905	C-terminal Pro-ghrelin Isoform Peptide, mouse	RRQLTSNHGQA	1267.4	1 mg
62358	Ghrelin (1-18)	GS-S(n-octanoyl)-FLSPEHQRVQQRKES	2227.5	1 mg
62931	[Ala ¹ , D-Trp ^{2,4} , Leu ⁶]-Ghrelin Receptor Agonist	AwFwLL	835	1 mg
62930	[D-Trp ^{1,3} , Leu ⁵]-Ghrelin Core-Ligand	wFwLL	763.9	1 mg
61434-1	Obestatin, human	FNAPFDVGIKLSGVYQQHSQAL-NH ₂	2547.9	1 mg
61434-5	Obestatin, human	FNAPFDVGIKLSGVYQQHSQAL-NH ₂	2547.9	5 mg
61435-1	Obestatin, rat, mouse	FNAPFDVGIKLSGAQYQQHGRAL-NH ₂	2517.9	1 mg
61435-5	Obestatin, rat, mouse	FNAPFDVGIKLSGAQYQQHGRAL-NH ₂	2517.9	5 mg
62041	Obestatin (11-23), human	LSGVYQQHSQAL	1458.6	1 mg
62040	Obestatin (11-23), rat, mouse	LSGAQYQQHGRAL	1428.6	1 mg

Glucagon-Like peptides

62070	GLP-2, Glucagon-like Peptide-2 (1-33), human	HADGSFSDEMNTILDNLAARDFINWLIQTKITD	3766.2	1 mg
62067	GLP-2, Glucagon-like Peptide-2 (1-33), rat	HADGSFSDEMNTILDNLATRDFINWLIQTKITD	3796.2	1 mg
62068	GLP-2, Glucagon-like Peptide-2 (1-34), human	HADGSFSDEMNTILDNLAARDFINWLIQTKITDR	3922.4	1 mg
62069	GLP-2, Glucagon-like Peptide-2 (1-34), rat	HADGSFSDEMNTILDNLATRDFINWLIQTKITDR	3952.4	1 mg

Growth Factor Fragments

61790	EGFR (688-701)-Lys	LVEPLTPSGEAPNQ-K	1579.8	1 mg
61788	EGFR (688-701)-Lys(FAM)	LVEPLTPSGEAPNQ-K(5-FAM)	1939.1	1 mg
61789	EGFR (688-701)-Lys(TAMRA), amide	LVEPLTPSGEAPNQ-K(5-TAMRA)-NH ₂	1993.3	1 mg
62989	Gelsolin (135-142)	KSGLKYKK	951.2	1 mg
62418	IGFBP6, Insulin Growth Factor Binding Protein 6, mouse	GPSEETTK	847.9	1 mg
62417	IGFBP6, Insulin Growth Factor Binding Protein 6, human	APAVAEENPK	1025.1	1 mg
62518	Link N, Cartilage Proteoglycan A Link Component	DHLSDNYTLDDHRAIH	1922	1 mg
62192	PDGF-BB, Platelet-Derived Growth Factor-BB, P62541	CSRNLIDC (Disulfide bridge:1-8)	921.1	1 mg

Guanylin

61643	Guanylin, rat	PNTCEICAYAAGTGC (Disulfide bridges: 4-12 and 7-15)	1515.8	1 mg
61617	Uroguanylin, human	NDDCELCVNVACTGCL (Disulfide bridges: 4-12 and 7-15)	1667.9	1 mg
61645	Uroguanylin, rat	TDECELCINVACTGC (Disulfide bridges: 4-12 and 7-15)	1569.8	1 mg

Heat Shock Protein Fragments

61900	HSP, Heat Shock Protein (3-13)	KTIAYDEEARR	1351.5	1 mg
62594	HSP-B8, Heat Shock Protein-B8 (10-29)	CHYPSRLRRDPFRDPSLSSR	2445.8	1 mg
62593	HSP-B8, Heat Shock Protein-B8 (181-194)	SFNNELPQDSQEVV	1607.7	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62592	HSP-B8, Heat Shock Protein-B8 (184-196), human	NELPQDSQEVTCT	1463.6	1 mg
62035	HSP-65, Heat Shock Protein-65 (458-474), human	NEDQKIGIEIHKRTLKI	2011.4	1 mg
61685	P38, HSP, <i>M. Leprae</i> Heat Shock Protein (411-425)	AGGGVTLTQAAPALD	1353.5	1 mg
61687	P38, HSP, <i>M. Leprae</i> Heat Shock Protein (417-429)	LLQAAPALDKLKL	1393.7	1 mg
61688	P38, HSP, <i>M. Leprae</i> Heat Shock Protein (522-534)	PEKTAAPASDPTG	1241.3	1 mg
61686	P61, HSP, <i>M. Leprae</i> Heat Shock Protein (343-355)	RVAQIRTEIENS	1530.7	1 mg
62748	Shepherdin (79-87)	KHSSGCAFL	949.1	1 mg

Hepatitis C Virus (HCV) Related Peptides

61863	HCV Core Protein (1-20)	MSTNPKPQRKTKRNTNRRPQ	2438.8	1 mg
61710	HCV, NS4A-Like Peptide	KKGSVVIVGRIVLSGKPAIIPKK	2388.0	1 mg
61864	HCV-1 e2 Protein (484-499)	PYCWHYPPKPCGIVPA	1828.2	1 mg
61865	HCV-1 e2 Protein (554-569)	WMNSTGFTKVCGAPPC	1699.0	1 mg

Histone Peptides

61855	Histone Deacetylase, HDAC Substrate	Ac-RGK(Ac)-AMC	600.7	1 mg
61761	Histone H1-derived Peptide	GGGPATPKKAKKL	1252.5	1 mg
61762	Histone H1-derived Peptide, FAM-labeled	5-FAM-GGGPATPKKAKKL	1610.8	1 mg
61741	Histone H1-derived Peptide, phosphorylated	GGGPA-pT-PKKAKKL	1332.5	1 mg
61742	Histone H1-derived peptide, phosphorylated, FAM-labeled	5-FAM-GGGPA-pT-PKKAKKL	1690.8	1 mg
61763	Histone H1-derived Peptide, TAMRA-labeled	5-TAMRA-GGGPATPKKAKKL	1665.0	1 mg
63674	Histone H1' (22-42), human, mouse	STDHPKYSDMIVAAIQAEKNR	2374.7	1 mg
63675	[Lys(Biotin) ⁹]-Histone H2A (1-12)-Cys	SGRGKQGG-K(Biotin)-ARAC	1501.8	1 mg
63676	Histone H2A (1-20)	SGRGKQGGKARAKAKTRSSR	2087.4	1 mg
62753	Histone H3 (1-20), N-Terminal	ARTKQTARKSTGGKAPRKQL	2183.6	1 mg
61701	Histone H3 (1-21), N-Terminal	ARTKQTARKSTGGKAPRKQLA	2254.6	1 mg
61702	Histone H3 (1-21)-GGK(Biotin), amide	ARTKQTARKSTGGKAPRKQLAGGK(Biotin)-NH ₂	2723.2	1 mg
61703	Histone H3 (1-25), amide	ARTKQTARKSTGGKAPRKQLATKAA-NH ₂	2625.1	1 mg
61704	Histone H3 (116-136), C116-136	KRVTIMPKDIQLARRIGERA	2508.0	1 mg
61705	Histone H3 (116-136), N15-39	APRKQLATKAARKSAPATGGVKKPH	2570.1	1 mg
63682	[Lys(Ac) ¹⁴]-Histone H3 (1-19), H3K14(Ac)	ARTKQTARKSTGG-K(Ac)-APRKQ	2112.4	1 mg
63680	[Lys(Ac) ⁹]-Histone H3 (1-20), H3K9(Ac)	ARTKQTAR-K(Ac)-STGGKAPRKQL	2225.6	1 mg
63677	[Lys(Me ₂) ⁴]-Histone H3 (1-21), H3K4(Me ₂)	ART-K(Me ₂)-QTARKSTGGKAPRKQLA	2282.6	1 mg
63679	[Lys(Me ₃) ⁹]-Histone H3 (1-24), H3K9(Me ₃)	ARTKQTAR-K(Me ₃)-STGGKAPRKQLATKA	2597	0.5 mg
63681	[Lys(Ac) ¹⁴]-Histone H3 (9-19), H3K14(Ac)	KSTGG-K(Ac)-APRKQ	1199.4	1 mg
62754	Histone H4 (1-7), N-Terminal	SGRGKGG	617.7	1 mg
62498	Histone H4 (1-20), PRMT7 Substrate, M1	SGRGKGGKGLGKGGAKRHRK-NH ₂	1991.3	1 mg
62500	Histone 4-12, p300/CBP Substrate	RGKGGKGLGKGA	1085.3	1 mg
62499	Histone 4-21, p300/CBP Substrate	SGRGKGGKGLGKGGAKRHRKV	2091.5	1 mg
62555	Histone 4 (1-21)-Lys(Biotin)	Ac-SGRGKGGKGLGKGGAKRHRKVGG-K(Biotin)	2602.1	1 mg
63685	[Lys(Ac) ¹²]-Histone H4 (1-20), H4K12(Ac)	SGRGKGGKGLG-K(Ac)-GGAKRHRK	2034.4	1 mg
63687	[Lys(Ac) ¹⁶]-Histone H4 (1-20), H4K16(Ac)	SGRGKGGKGLGKGGK-Ac)-RHRK	2034.4	1 mg
63686	[Lys(Ac) ¹⁶]-Histone H4 (14-21), Mca-labeled, H4K16(Ac)(14-21), pepH4-Ack16	Mca-GA-K(Ac)-RHRKV-NH ₂	1208.4	1 mg
63683	[Lys(Ac) ⁵]-Histone H4 (1-20), H4K5(Ac)	SGRG-K(Ac)-GGKGLGKGGAKRHRK	2034.4	1 mg
63684	[Lys(Ac) ⁸]-Histone H4 (1-20), H4K8(Ac)	SGRGKGG-K(Ac)-GLGKGGAKRHRK	2034.4	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
HIV Related Peptides				
61439	cDDX4, HIV Fusion Molecule Analog	cyclo(DNVSEADDRYIG)	1335.4	1 mg
61599	G9-209-2L3W HIV	ILWQVPFSV	1088.3	1 mg
62435	HIV-1, Envelope Glycoprotein (187-203)	RGPGRAFVTIGKIGNMR-NH ₂	1829.2	1 mg
62436	HIV-1, Envelope Glycoprotein (187-203), Biotin-labeled	Biotin-RGPGRAFVTIGKIGNMR-NH ₂	2055.5	1 mg
61527	HIV-1 p17 Gag (20-28)	RLRPGGKKK	1039.3	1 mg
61526	HIV-1 p17 Gag (77-85), SL9 Epitope	SLYNTVATL	981.1	1 mg
61674	HIV Nef (90-100)	FLKEKGGLEGL	1190.4	1 mg
61672	HIV QK10	QVPLRMTYK	1135.4	1 mg
62800	HIV V3 Loop P18 Peptide	RIQRGPGRAFVTIGK	1656	1 mg
62101	gp41 HIV Fragment, C34	WMEWDREINNYTSLIHSLIEESQN-QQEKNQELL	4248.6	1 mg
62103	gp41 HIV Fragment, C34-LC-Biotin	WMEWDREINNYTSLIHSLIEESQNQQEKNQ ELLK(LC-Biotin)-NH ₂	4715.3	0.5 mg
62100	gp41 HIV Fragment, N36, HR1	SGIVQQQNLLRAIEAQQHLLQLTVW-GIKQLQARIL	4122.8	1 mg
61353	HIV gag p24 Peptide (197-205)	AMQMLKETI	1064.3	1 mg
61673	HIV Nef (68-76)	FPVTPQVPL	997.2	1 mg
62290	[Pro ¹⁵]-Gag Region of HIV-1 Subtype C	HKGRPGNFLQNRPEPTAP	2016.3	1 mg
62694	HIV-1 env (584-594)	RYLRDQQLLGI	1374.6	1 mg
61524	HIV-1 RT, Reverse Transcriptase IV9 Epitope	ILKEPVHGV	991.2	1 mg
61525	HIV-1 RT, Reverse Transcriptase VL9 Epitope	VIYQYMDL	1159.3	1 mg
62389	RANTES (11-22)	CFAYIARPLPRA	1377.7	1 mg
62583	V3 gp120 HIV (AD-11)	SIPIGPGRAFYT	1379.6	1 mg
62585	V3 gp120 HIV (BK-130)	RIHIGPGRALYT	1454.7	1 mg
62582	V3 gp120 HIV (JR-FL)	SIHIGPGRAFYT	1419.6	1 mg
62584	V3 gp120 HIV (WM)	GIHIGPGKAFYT	1361.6	1 mg
62434	VIR-165, α 1-Antitrypsin Modification (353-372)	LEAIPCSIPPCFAFNKPFVF	2240.7	1 mg
63719	VprBP (1493-1507)	DNSDLEDDIILSLNE	1704.8	1 mg
Hydrocarbon-Stapled Peptides				
63733-025	FITC- β -Ala-NYAD-1 (NYAD-2)	FITC-(β -A)-ITFXDLLXYYGP-NH ₂ (with special cyclization to get double bond, X= (S)- α -(2'-pentenyl)alanine)	1911.3	0.25 mg
63735-1	Hydrocarbon-Stapled Helix SAHB _A Peptide	EDIIRNIARHLAXVGDX-NIe-DRSIW (with special cyclization to get double bond, X= (S)-2-(2'-pentenyl)alanine)	2713.1	1 mg
63735-5	Hydrocarbon-Stapled Helix SAHB _A Peptide	EDIIRNIARHLAXVGDX-NIe-DRSIW (with special cyclization to get double bond, X= (S)-2-(2'-pentenyl)alanine)	2713.1	5 mg
63734-1	NYAD-1	ITFXDLLXYYGP-NH ₂ (with special cyclization to get double bond, X= (S)- α -(2'-pentenyl)alanine)	1450.7	1 mg
63734-5	NYAD-1	ITFXDLLXYYGP-NH ₂ (with special cyclization to get double bond, X= (S)- α -(2'-pentenyl)alanine)	1450.7	5 mg
Ion Channel Peptides				
63697	C-Peptide-1, Rat	EVEDPQVPQLELGGGPEAGDLQTLALEVARQ	3259.6	0.5 mg
62696	Kv3, Channel Containing Protein (567-585)	CKESPVIKYMPTAEVRVT	2122.5	1 mg
61952	Microtubule-associated Protein 7, MAP7, (CT), mouse	TQQTAIEVI	889.0	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62378	Optimal NHERF-1 PDZ1-Binding Motif	VQDTRL	730.8	1 mg
62390	Skeletal Dihydropyridine Receptor (671-690)	TSAQKAKAEERKRRKMSRGL	2331.7	1 mg
62061	TRPA1, Transient Receptor Potential A1 (CT), human	CVLNAVKTkTHCSISHPDI	2066.4	1 mg
61951	TRPV4, Transient Receptor Potential Vanilloid 4 (CT), human	CDGHQQGYAPK	1203.3	1 mg
61953	TRPV4, Transient Receptor Potential Vanilloid 4 (NT), human	MDQAKSAE	879.0	1 mg
61954	TRPV4, Transient Receptor Potential Vanilloid 4 (CT), rat	CDGHQQGYAPKWRAEDAPL	2142.3	1 mg
62226	V-ATPase a-subunit	FRSEEMCLSQLFLQ	1731	1 mg

Influenza Viral Peptides

62406	Influenza A M2 Coat Protein (22-46)	SSDPLVVAASIIGILHLILWILDRL-NH ₂	2699.3	1 mg
61666	Influenza A M2 Coat Protein (22-46), acetylated, amide	Ac-SSDPLVVAASIIGILHLILWILDRL-NH ₂	2769.4	1 mg
61670	Influenza A NP (366-374)	ASNEMNDAM	982.1	1 mg
62036	Influenza Gene Product, PB1F2 (62-70)	LSLRNPILV	1024.3	1 mg
62288	Influenza Matrix Protein (61-72), MP (61-72)	GFVFTLTPSER	1352.6	1 mg
61519	Influenza Peptide, NS2 (114-121)	RTFSFQLI	1011.2	1 mg
61636	Influenza Peptide, PA (224-233)	SSLENFRAYV	1185.3	1 mg
61517	Influenza Peptide, PB1 (703-711)	SSYRRPVGI	1034.2	1 mg
61696	Influenza Virus A/Aichi/2/68 HA1 (195-209), X-31	YVQASGRVTVSTRRS	1666.9	1 mg
62423	Influenza Virus Hemagglutinin, HA (326-340)	KQNTLKLATGMRNVP	1671	1 mg
62421	Influenza Virus Nucleoprotein, NP (266-274)	LRGSVAHKS	954.1	1 mg
62422	Influenza Virus Nucleoprotein, NP (277-286)	PACVYGPVA	947.1	1 mg
62420	Influenza Virus Nucleoprotein, NP (311-325)	QVYSLIRPNENPAHK	1767	1 mg

Kinase / Phosphatase Substrates

61749	Abltide	KKGEAIYAAPFA-NH ₂	1264.5	1 mg
61751	Abltide, FAM-labeled	5-FAM-KKGEAIYAAPFA-NH ₂	1622.8	1 mg
61750	Abltide, TAMRA-labeled	5-TAMRA-KKGEAIYAAPFA-NH ₂	1677.0	1 mg
61784	ADR1, Yeast Alcohol Dehydrogenase (ADH2) Transcription Activator (222-234)	LKKLTRRASFSGQ	1491.8	1 mg
61783	ADR1, Yeast Alcohol Dehydrogenase (ADH2) Transcription Activator (222-234), FAM-labeled	5-FAM-LKKLTRRASFSGQ	1850.1	1 mg
61732	Alternate Syntide	PLSRTLVSSSLPGL-NH ₂	1425.7	1 mg
61733	Alternate Syntide, FAM-labeled	5-FAM-PLSRTLVSSSLPGL-NH ₂	1784.0	1 mg
62596	AMARA Peptide	AMARAASAAALARRR	1542.8	1 mg
61747	AMPKtide	LKKLTRRASFSAQ	1505.8	1 mg
61748	AMPKtide, FAM-labeled	5-FAM-LKKLTRRASFSAQ	1864.1	1 mg
62808-1	AMPKtide, ADR1, FAM-labeled	5-FAM-LKKLTRRRPSFSAQ	1890.1	1 mg
62808-5	AMPKtide, ADR1, FAM-labeled	5-FAM-LKKLTRRRPSFSAQ	1890.1	1 mg
62407	Aurora Kinase Substrate, M peptide, Biotin-LC-labeled	Biotin-LC-LRRWSLGLRRWSLGLRRWSL-GLRRWSLG	3833.7	1 mg
61728	BTK-derived Peptide	KKVVALYDYMPMN-NH ₂	1571.0	1 mg
61729	BTK-derived Peptide, FAM-labeled	5-FAM-KKVVALYDYMPMN-NH ₂	1929.3	1 mg
61726	CDc25C-derived Peptide	VSRSGLYRSPSPENLNRP	2316.6	1 mg
61727	CDc25C-derived Peptide, FAM-labeled	5-FAM-VSRSGLYRSPSPENLNRP	2675.0	1 mg
61758	CDK7tide	YSPTSPSYPTSPSYPTSPS	2177.3	1 mg
61760	CDK7tide derived Peptide, FAM-labeled	5-FAM-FLAKSFGSPNRAYKK	2072.3	1 mg
61759	CDK7tide, TAMRA-labeled	5-TAMRA-YSPTSPSYPTSPSYPTSPS	2589.8	1 mg
63367	CDK7/9 tide	YSPTSPSYPTSPSYPTSPSKKKK	2690	1 mg
61738	CDPKS, Syntide analog	PLRRTLSVAA	1083.3	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62807-1	CHK1tide, TAMRA-labeled	5-TAMRA-ALKLVRYPSFVITAK-NH ₂	2117.6	1 mg
62807-5	CHK1tide, TAMRA-labeled	5-TAMRA-ALKLVRYPSFVITAK-NH ₂	2117.6	5 mg
62806-1	CHK1tide, FAM-labeled	5-FAM-ALKLVRYPSFVITAK-NH ₂	2063.4	1 mg
62806-5	CHK1tide, FAM-labeled	5-FAM-ALKLVRYPSFVITAK-NH ₂	2063.4	5 mg
62819	[pSer ³⁴⁵]-Chk1	QGISF-pS-QPTC	1147.2	1 mg
62820	[Ser ³⁴⁵]-Chk1 (Paired)	QGISFSQPTC	1067.2	1 mg
63790	CHK 1/2 Peptide Substrate, CDC25C (206-225), FAM-labeled	5-FAM-KKKVSRSGLYRSPSPENLNRP	3059.5	1 mg
62826	CHK2tide	AMRLERQDSIFYPK-NH ₂	1753.1	1 mg
61767	CK1tide, amide	HAAIGDDDDAYSITS-NH ₂	1549.6	1 mg
61769	CK1tide, TAMRA-labeled	5-TAMRA-HAAIGDDDDAYSITS-NH ₂	1962.1	1 mg
61768	CK1tide, FAM-labeled	5-FAM-HAAIGDDDDAYSITS-NH ₂	1907.9	1 mg
61773	CK2tide, FAM-labeled	5-FAM-RRRADDSDDDDD-NH ₂	1807.7	1 mg
63795	CK1 Peptide Substrate [pSer ⁷], FAM-labeled	5-FAM-KRRRAL-pS-VASLPGL-NH ₂	1961.2	1 mg
61286	[pSer ¹¹⁹]-CREB327 (113-122)	ILSRRP-pS-YRK	1355.5	1 mg
61287	[pSer ¹¹⁹]-CREB327 (113-123)	ILSRRP-pS-YRKI	1468.7	1 mg
61285	CREB327 (113-126)	ILSRRPSYRKILND	1731.0	1 mg
61288	[pSer ¹¹⁹]-CREB327 (113-126), Biotin-labeled	Biotin-ILSRRP-pS-YRKILND	2037.3	1 mg
61770	CREBtide	GEILSRRPSYRK-NH ₂	1460.7	1 mg
61771	CREBtide, FAM-labeled	5-FAM-GEILSRRPSYRK-NH ₂	1819.0	1 mg
62595	[pSer ¹³³ , Leu ¹²³]-CREBtide Peptide Substrate, Biotin-LC-labeled	Biotin-LC-LRREILSRRP-pS-YRK-NH ₂	2248.7	1 mg
63706	Crosstide[GRPRTSSFAEG-NH ₂], FAM-labeled	5-FAM-GRPRTSSFAEG-NH ₂	1521.6	1 mg
61781	DAP-kinase/MLCK, Death-Associated Protein Kinase Substrate / Myosin Light Chain Kinase Substrate	KKRPQRRYSNVF	1578.9	1 mg
61782	DAP-kinase/MLCK, Death-Associated Protein Kinase Substrate / Myosin Light Chain Kinase Substrate, FAM-labeled	5-FAM-KKRPQRRYSNVF	1937.2	1 mg
62034	<i>Dictyostelium</i> Myosin II Heavy Chain (2020-2035), MH-1	RKKFGESEKTKKEFL	1956.3	1 mg
62698	Dyrktide	RRRFRPASPLRGPPK	1791.2	1 mg
62716	Dyrktide, FAM-labeled	5-FAM-RRRFRPASPLRGPPK	2149.5	1 mg
62715	Dyrktide, TAMRA-labeled	5-TAMRA-RRRFRPASPLRGPPK	2203.6	1 mg
61777	Erktide	IPTTPITTTYFFFK	1677.0	1 mg
61778	Erktide, FAM-labeled	IPTTPITTTYFFFK-5-FAM	2036.3	1 mg
62342	ERK 1/2 /MAP Kinase Substrate, FAM-labeled	5-FAM-IPTSPITTTYFFFKKK	2277.6	1 mg
61744	Forkhead-derived Peptide, Woodtide, FAM-labeled	5-FAM-KKISGRLSPIMTEQ-NH ₂	1945.2	1 mg
61743	Forkhead-derived Peptide, Woodtide	KKISGRLSPIMTEQ-NH ₂	1586.9	1 mg
61875	Generic Protein Kinase Substrate	RARTLSFAEPG	1204.4	1 mg
61785	Glycogen Synthase-derived Peptide	KKLNRTLVA	1129.4	1 mg
62630	Glycogen Synthase derived Peptide, Biotin-labeled	Biotin-KKLNRTLVA	1355.7	1 mg
61786	Glycogen Synthase derived Peptide, FAM-labeled	5-FAM-KKLNRTLVA	1487.7	1 mg
61787	Glycogen Synthase derived Peptide, TAMRA-labeled	5-TAMRA-KKLNRTLVA	1541.9	1 mg
61724	Glycogen Synthase-derived Peptide, phosphorylated	KKLNRTL-pS-VA	1209.4	1 mg
61725	Glycogen Synthase-derived Peptide, phosphorylated, TAMRA-labeled	5-TAMRA-KKLNRTL-pS-VA	1621.9	1 mg
61740	Glycogen Synthase Kinase 3 (GSK-3), FAM-labeled	5-FAM-RRAAEELDSRAG-pS-PQL-NH ₂	2193.2	1 mg
62354	Glycogen Synthase Kinase 3 α (GSK-3 α) (16-26), phosphorylated	RARTS-pS-FAEPG	1258.3	1 mg
62384	GSM Peptide	RRRPASVPPSPSLSRHSSHQRR	2551.9	1 mg
61730	IP3R, Inositol 1,4,5-Triphosphate Receptor	GRRESLTSFG-NH ₂	1108.2	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61731	Inositol 1,4,5-Triphosphate Receptor (IP3R), FAM-labeled	5-FAM-GRRESLTSFG-NH ₂	1466.5	1 mg
61764	IRS1-derived Peptide	KKSRGDYMTMQIG-NH ₂	1513.8	1 mg
61765	IRS1-derived Peptide, FAM-labeled, amide	5-FAM-KKSRGDYMTMQIG-NH ₂	1872.1	1 mg
61766	IRS1-derived Peptide, TAMRA-labeled, amide	5-TAMRA-KKSRGDYMTMQIG-NH ₂	1926.3	1 mg
62814	IRS-1 (650-666)	KKRVDPNGYMMMSPSGV	1897.3	1 mg
62813	IRS-1 (718-735)	KKKLPATGDYMNMSVGD	1952.3	1 mg
62811	IRS-1 (933-948)	KKGSEEYMNMDLGPGR	1812.1	1 mg
62546	IRS-1 (979-989), Peptide-983, amide	SRGDYMTMQIG-NH ₂	1257.5	1 mg
62812	IRS-1 (1165-1177)	KKKNGLNIDLDL	1533.8	1 mg
61774	ITAM, Immunoreceptor Tyrosine-Based Activation Motif, IgE (FcεRI) Receptor	pY-TGLSTRNQET-pY-ETL-NH ₂	1934.9	1 mg
61775	ITAM, Immunoreceptor Tyrosine-Based Activation Motif, IgE (FcεRI) Receptor, FAM-labeled	5-FAM-pY-TGLSTRNQET-pY-ETL-NH ₂	2293.2	1 mg
61776	Immunoreceptor Tyrosine-Based Activation Motif (ITAM), IgE (FcεRI) Receptor, TAMRA-labeled	5-TAMRA-pY-TGLSTRNQET-pY-ETL-NH ₂	2347.4	1 mg
62537	Jak3tide	GGEEEEYFELVKKKK	1813	1 mg
62825-1	JJMO-5	FAM-AMRLERQDSIFYPK-NH ₂	2111.4	1 mg
62825-5	JJMO-5	FAM-AMRLERQDSIFYPK-NH ₂	2111.4	5 mg
62017-1	Kinase Substrates Library, Group I, biotinylated, 180 distinct peptide mixtures			1 Set
62017-5	Kinase Substrates Library, Group I, biotinylated, 180 distinct peptide mixtures			1 Set
62335	Kinase Substrates Library, Group II, biotinylated, 18 distinct peptide mixtures			1 Set
62219	Leukocyte Common Antigen-related (LAR) PTP Receptor, LARFN5C (29-48), rat	PGPLQAKPFTILSPFLSSRC	2159.6	1 mg
62157	MAPK-activated Protein Kinase 3 (MK3)	DYDQVKIKDLKTSNNRLLNKRKKQAGSS-SAS	3650.1	1 mg
62156	MAPK-activated Protein Kinase 5 (MK5)	IQDLKVSLKPLHSVNNPILRKRKLLGTPKDSV	3736.5	1 mg
62154	Mitogen-and Stress-activated Kinase1 (MSK1), D-domain Fragment 1	FNKYKREGFCLQNVDKAPLAKRRKMKKTST-STE	3904.6	1 mg
62155	Mitogen-and Stress-activated Kinase2 (MSK2), D-domain Fragment 2	FNRGKREGFFLKSVENAPLAKRRKQKLRSATAS	3792.4	1 mg
62400	Mitogen-activated Protein Kinase Kinase 2 (363-374), MAPK Kinase (363-374), human, monkey	LNQPGTPTRTAV	1254.4	1 mg
62153	Mitogen-activated Protein (MAP) Kinase-activated Protein Kinase 2a (MK2a), (370-400)	IKIKKIEDASNPLLLKRRKKARALEAAALAH	3467.2	1 mg
62249	MNK1, MAPK-interacting Kinase 1 Fragment	ALADGLCSMKLSPCKSRLARRRALAQA-GRGED (S-S bond)	3497.1	1 mg
61734	p34cdc2-derived Peptide	KVEKIGEGTYGVV-NH ₂	1377.6	1 mg
61736	p34cdc2-derived Peptide, FAM-labeled, amide	5-FAM-KVEKIGEGTYGVV-NH ₂	1735.9	1 mg
61735	p34cdc2-derived Peptide, TAMRA-labeled, amide	5-TAMRA-KVEKIGEGTYGVV-NH ₂	1790.1	1 mg
61752	p38tide	IPTTPITTTYFFFK-NH ₂	1676.0	1 mg
61754	p38tide, FAM-labeled	5-FAM-IPTTPITTTYFFFK-NH ₂	2034.3	1 mg
61753	p38tide, TAMRA-labeled	5-TAMRA-IPTTPITTTYFFFK-NH ₂	2088.5	1 mg
61876	p120ctn, human	QDEGQESLEEELDVLVLDDEGGQVSYPSM	3211.4	1 mg
61874	PAKtide	RRRLSFAEPG	1188.4	1 mg
62948	[Ala147]-PKC-δ (137-153)	AMFPTMNRGSIKQAKI	1949.4	1 mg
61792	PKCε Pseudosubstrate-derived Peptide	ERMRRPRKRQGSVRRRV-NH ₂	2066.5	1 mg
61791	PKCε Pseudosubstrate-derived Peptide, TAMRA-labeled	5-TAMRA-ERMRRPRKRQGSVRRRV-NH ₂	2479.0	1 mg
62025	PKC ζ 410	GDTTSTFCGTPN	1200.3	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62020	PKC ζ 560	EPVQLTPDDEDA	1328.4	1 mg
63695	Myristoylated PKC ζ	Myristoyl-RLYRKRIWRSAGR	1928.5	1 mg
63720	[Ala ¹⁴⁷]-PKC δ (137-153), FAM-labeled	5-FAM-AMFPTMNRGSIKQAKI	2307.7	1 mg
61737	Phospho PKCtide	5-FAM-RFARKG-pS-LRQKNV-NH ₂	1997.2	1 mg
63361	Myristoylated PKC ζ, Pseudosubstrate (ZIP)	Myr-SIYRRGARRWRKL	1928.5	1 mg
61780	PLM-derived Peptide	IRRLSTRRR	1213.5	1 mg
61779	PLM derived Peptide, TAMRA-labeled	5-TAMRA-IRRLSTRRR	1626.0	1 mg
62387	Polo-like Kinase (134-140)	RRRSLLE	929.1	1 mg
62353	Protein Kinase Substrate LSFAEPG Motif 2, FAM-labeled	5-FAM-KKLNRTLSSFAEPG-NH ₂	1818	1 mg
62158	Ribosomal Protein S6 Kinase 2 (RSK2), (708-740)	SALNRNQSPVLEPVGRSTLAQRRGIKKITSTAL	3563.1	1 mg
62079	Ribosomal Protein S6 Kinase 3 (RSK3), (709-741)	ALNRTPQAPRLEPVLSSNLAQRRGMKRLT-STRL	3732.4	1 mg
61722	RS Domain-derived Peptide	GRSRSRSRSR	1204.3	1 mg
61723	RS Domain-derived Peptide, FAM-labeled	5-FAM-GRSRSRSRSR	1562.6	1 mg
62699	SGKtide, SGK Substrate	KKRNRRLSVA	1227.5	1 mg
63771	Soymetide-4, MITL	MITL	476.6	1 mg
62810	Src-tide	GEEIYGEFD-NH ₂	1057.1	1 mg
62343	Src Optimal Peptide Substrate, FITC-LC-labeled	FITC-LC-AEEIYGEFEAKKKK	2301.6	1 mg
61755	Synapsin I-derived Peptide	LRRRLSDANF-NH ₂	1246.4	1 mg
61756	Synapsin I-derived Peptide, FAM-labeled	5-FAM-LRRRLSDANF-NH ₂	1604.8	1 mg
61757	Synapsin I-derived Peptide, TAMRA-labeled	5-TAMRA-LRRRLSDANF-NH ₂	1658.9	1 mg
63673	T308tide	KTFCGTPEYLAPEVRR	1967.2	1 mg
61320	Zipper-interacting Protein Kinase (ZIPK), (264-276)	MTIAQSLEHSHWIK	1543.8	1 mg
61319	Zipper-interacting Protein Kinase (ZIPK), (309-327)	SHSSLPPNNSYADFERSK	2183.3	1 mg
63366	ZIptide	KKLNRTLSSFAEPG	1460.7	1 mg

LL-37 / CRAMP Peptides

63710	LL-37, acetylated & amidated	Ac-LLGDFFRKSKEKIGKEFKRIVQRIKDFL-RNLVPRTES-NH ₂	4534.4	1 mg
61302	LL-37, Antimicrobial Peptide, human	LLGDFFRKSKEKIGKEFKRIVQRIKDFL-RNLVPRTES	4493.3	1 mg
63694	LL-37, FAM-labeled with 6-carbon spacer(LC)	5-FAM-LC-LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLVPRTES	4964.9	0.1 mg
63711	LL-37 fragment (8-37), LL-8-37	KSKEKIGKEFKRIVQRIKDFLRNLVPRTES	3644.3	1 mg
63712	LL-37 fragment (18-37), LL-18-37	KRIVQRIKDFLRNLVPRTES	2468.9	1 mg
62208	LL-37, reverse sequence	SETRPVLNRLFDKIRQVIRKFEKGIKEK-SKRFFDGLL	4493.3	0.5 mg
61309	LL-37 pentamide	LLGNFFRKSQKIG-KQFKRIVQRIKNFFRNLVPRTS	4522.4	1 mg
63708	LL-37, scrambled	GLKLRFEFSKIKGEFLKTPEVRFDRDIKLDN-RISVQR	4493.3	1 mg
63693	Biotin-LC-LL-37; LL-37, biotin labeled with 6-carbon spacer(LC)	Biotin-LC-LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLVPRTES	4832.8	0.5 mg
63692	Cys-LC-LL-37; LL-37 with Cys and 6-carbon spacer(LC) at the N-Terminal	C-LC-LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLVPRTES	4709.7	1 mg
61313	RL-37, Rhesus Monkey Antimicrobial Protein	RLGNFFRKYKEKIGGGLKKVGGQIKDFL-GNLVPRTS	4100.9	0.5 mg
61305	mCRAMP, mouse	GLLRKGGEKIGEKLK-KIGQKIKNFFQKLVQPPEQ	3878.7	1 mg
63709-05	mCRAMP, mouse, biotin labeled with 6-carbon spacer(LC)	Biotin-LC- GLLRKGGEKIGEKLK-KIGQKIKNFFQKLVQPPEQ	4218.2	0.5 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
63709-1	mCRAMP, mouse, biotin labeled with 6-carbon spacer(LC)	Biotin-LC- GLLRKGGEKIGEKLK-KIGQKIKNFFQKLVPQPEQ	4218.2	1 mg
61306	rCRAMP	GLVRKGGEKFGEKLRKIGQKIKEFFQKLA-LEIEQ	3946.7	1 mg

Lymphocytic Choroemeningitis Virus (LCMV) Peptides

61296	LCMV (33-41), gp33	KAVYNFATM	1044.3	1 mg
62538	LCMV (70-77), gp70	GVYQFKSV	927.1	1 mg
62541	LCMV (166-173), gp166	ITIQYNLT	965.1	1 mg
62542	LCMV (221-228), gp221	SQTSYQYL	989.1	1 mg
62539	LCMV (276-286), gp276	SGVENPGGYCL	1095.2	1 mg
62540	LCMV (365-372), gp365	MGVPYCNV	946.1	1 mg
62543	LCMV (2062-2069), L2062	RSIDFERV	1021.2	1 mg
61669	LCMV Glycoprotein 33, (LCMV gp1), gp 33 (33-41)	KAVYNFATC	1016.2	1 mg
61699	LCMV Nucleoprotein Fragment (LCMV-NP 205)	YTVKYPNL	997.2	1 mg
61700	LCMV Nucleoprotein Fragment (LCMV-NP 396 H-2Db)	FQPQNGQFI	1078.2	1 mg

MAGE Peptides

61469	MAGE-1 (161-169)	EADPTGHSY	976.0	1 mg
61356	MAGE-1 (230-238)	STAPPAHGV	1020.2	1 mg
61490	MAGE-3 (112-120)	KVAELVHFL	1055.3	1 mg
62702	MAGE-3 (114-127)	AELVHFLLLKYRAR	1729.1	1 mg
62701	MAGE-3 (121-134)	LLKYRAREPVTKAE	1674	1 mg
61354	MAGE-3 (161-169)	EVDPIGHLY	1042.2	1 mg
61355	MAGE-3 (271-279)	FLWGPRALV	1058.3	1 mg
62709	MAGE-A1 (96-104)	SLFRAVITK	1034.3	1 mg
61488	MAGE-A1 (237-245)	KLLTQDLVQ	1057.3	1 mg
62365	MAGE-A1 Antigen (278-286), human	KVLEYVIKV	1090.4	1 mg
61489	MAGE-A2 (112-120)	KMVELVHFL	1115.4	1 mg
61503	MAGE-A2 (157-166)	YLQLVFGIEV	1180.4	1 mg
61504	[Ile ¹⁶¹]-MAGE-A2 (157-166)	YLQLIFGIEV	1194.4	1 mg
61498	MAGE-A3 (167-176)	MEVDPIGHLY	1173.4	1 mg
61499	MAGE-A3 (195-203)	IMPKAGLLI	955.3	1 mg
62364	MAGE-A4 Antigen (230-239), human	GVYDGREHTV	1132.2	1 mg
62708	MAGE-A10 (183-191)	MLLVFGIDV	1006.3	1 mg
61457	MAGE-C2 (336-344)	ALKDVEERV	1058.2	1 mg

Matrix Metalloproteinase Substrates and Related Peptides

62027	390 MMP FRET Substrate V	Mca-PLGLEEA-Dap(Dnp)-NH ₂	1193.2	1 mg
62028	390 MMP FRET Substrate VI	Mca-PLA-C(OMeBn)-WAR-Dap(Dnp)-NH ₂	1403.5	1 mg
62052	390 MMP FRET Substrate VII	Mca-PLGLAR-Dap(Dnp)	1094.2	1 mg
63728-01	FRETS-VWF73 (Fluorescence-Quenching Substrate for ADAMTS-13)	DRE-Dap(Nma)-APNLVYMTG-Dap(Dnp)-PASDEIKRLPGDIQVVPIGVGPAN-VQELERIGWPNAPILIQDFETLPREAPDLVLQR	8314.3	0.1 mg
63728-05	FRETS-VWF73 (Fluorescence-Quenching Substrate for ADAMTS-13)	DRE-Dap(Nma)-APNLVYMTG-Dap(Dnp)-PASDEIKRLPGDIQVVPIGVGPAN-VQELERIGWPNAPILIQDFETLPREAPDLVLQR	8314.3	0.5 mg
62054	Gelatinase Inhibitor, CTT	CTTHWGFTLC (Disulfide bridge:1-10)	1166.4	1 mg
62053	Gelatinase Inhibitor Negative Control	STTHWGFTLS	1136.2	1 mg

CAT#	Product Name	Sequence	MW	Size
Melan-A and Mucin Related Peptides				
61476	Melan-A/MART-1 (24-34)	AEEAAGIGILT	1044.2	1 mg
61497	Melanosomal Antigen II	DAEKCDKTDEY	1316.4	1 mg
61357	MUC1 (130-138)	STAPPAHGV	835.9	1 mg
61451	MUC1, tandem repeat fragment	PDTRPAPGSTAPPAHGVTS	1887.0	1 mg
62648	MUC1 TR	PAPGSTAPPAHGVTSAPDTR	1887	1 mg
61337	MUC2	PTTTPISSTTVPTPTPTGTQT	2201.4	1 mg
61338	MUC3	HSTPSFTSSITTETTS	1784.9	1 mg
61339	MUC4	TSSASTGHATPLPVD	1541.6	1 mg
61334	MUC5AC 3	GT-T*-PSPVPTTSTTSAP (T*= GalNAc-modified residue)	1703.9	1 mg
61329	MUC5AC Analog 1	GTTTSPVPTTSTTSAP	1501.6	1 mg
61330	[Ala ⁶ , Val ¹⁵] MUC5AC Analog 2	GTTTSAVPTTSTTSVP	1503.6	1 mg
61331	[Ile ¹² , Val ¹⁵] MUC5AC Analog 3	GTTTSPVPTTSTTSVP	1541.7	1 mg
61335	MUC5AC, Analog B	TTSTTSAPTTS	1054.1	1 mg
61333	MUC5AC-13	GTTTSPVPTTST-T*-SAP (T*= GalNAc-modified residue)	1704.6	1 mg
61332	MUC5AC-3/13	GT-T*-PSPVPTTST-T*-SAP (T*= GalNAc-modified residue)	1907.6	1 mg
61340	MUC6a	SFQTTTTYPTPSHPQTTL	2105.3	1 mg
61341	MUC6b	TSLSHTSSTHHPEVPT	1855.0	1 mg
61342	MUC7	PTPSATTPAPPSSSAPPETAA	2037.2	1 mg
Melanocyte Stimulating Hormones and Analogs				
61856-5	Melanotropin-Potentiating Factor, MPF	KKGE	460.5	5 mg
62367	RAB38/NY-MEL-1 (50-58)	VLHWDPETV	1095.2	1 mg
61906	VA-β-MSH, Lipotropin-7, Proopiomelanocortin-derived	VAAEKKDEGYPYRMEHFRWGSPPKD	2831.2	1 mg
MHC Related Peptides				
62552	2Wp, Class II MHC Molecule	AWGALANWAVDS	1260.4	1 mg
62522	Adeno-Associated Virus Epitope 2 (AAV2)	VPQYGYLTL	1053.2	1 mg
62388	Apolipoprotein B-100 (3136-3155), human	KTTKQSFDSLVSQAQYKKNKH	2379.8	1 mg
62756	BDC2.5 mimotope 1040-31	YVRPLWVRME	1348.6	1 mg
62755	BDC2.5 mimotope 1040-51	RVLPLWVRME	1298.6	1 mg
61624	CLIP, Class II-Associated Invariant Chain Peptide (85-99)	KPVSQMRMATPLLMR	1759.3	1 mg
62567	CLIP (85-100), Biotin-LC-labeled	Biotin-LC-KPVSKMRMATPLLMQA-NH ₂	2140.8	1 mg
62438	CLIP (85-101), CD74 Antigen (85-101)	KPVSQMRMATPLLMRPM	1987.6	1 mg
62569	CLIP, Class II-Associated Invariant Chain Peptide (86-100)	PVSKMRMATPLLMQA	1674.1	1 mg
62568	CLIP, Class II-Associated Invariant Chain Peptide (88-99)	SKMRMATPLLMQ	1406.8	1 mg
61667	CPα1, F-actin Capping Protein Subunit 1 (92-99)	ISFKFDHL	1006.2	1 mg
62689	Cytomegalovirus (CMV) CTL Epitope (pp65)	QYDPVAALF	1023.2	1 mg
61046	Dby HY Peptide	NAGFNSNRANSSRSS	1568.6	1 mg
62195	dsDNA Mimotope Peptide - on 8-branch MAPS	DWEYSVWLSN-on 8 branch MAPS	11229.5	1 mg
62371	DQ (65-79), α Chain HLA II fragment	NIAVLKHNLNIVIKR	1745.2	1 mg
62828	DWEYS-Peptide --on 8 branch MAPS	DWEYS-on 8 branch MAPS	6432	1 mg
61621	Eα, Class II MHC, Eα Chain Peptide (52-68)	ASFEAQGALANIAVDKA	1675.9	1 mg
62690	EBV 2A.32	TYGPFVMSL	1014.2	1 mg
62218	EYK(EYA) ₃ EYK	EYKEYAEYAEYAEYK	1949.1	1 mg
63718	α-9 gliadin Peptide	QVLQQSTYQLLQELCCQHLW	2461.9	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61899	α -gliadin (57-73)	QLQPFQPELPYPQPQS	1994.2	1 mg
62757	Glutamic Acid Decarboxylase (GAD65) (524-543), p524-543	SRLSKVAPVIKARMMEYGTT	2238.7	1 mg
62402	Immunodominant Mouse Minor Histocompatibility Antigen (MiHA)	AAPDNRETF	1020.1	1 mg
62297	IRBP (1-20), human	GPTHLFQPSLVLDMAKVLLD	2194.6	1 mg
62553	3Kp2, Class II MHC Molecule	ASFEAQKAKANKAVD	1578.8	1 mg
62362	Myelopeptide-1 (MP-1)	FLGFPT	680.8	1 mg
62363	Myelopeptide-2 (MP-2)	LVVYPW	776	1 mg
62655	NY-ESO-1 (87-111)	LLEFYLAMPFATPMEAEELARRSLAQ	2869.4	1 mg
62656	NY-ESO-1 (158-180)	LLMWITQCFLPVFLAQPPSGQRR	2702.3	1 mg
61622	p3K, [Lys ^{58,60,63}]-E α (52-68), Class II MHC E α Chain Peptide	ASFEAQKAKANKAVDKA	1777.0	1 mg
61623	p3K truncated, [Lys ^{58,60,63}], Class II MHC E α Chain Peptide	FEAKGAKANKAVD	1348.5	1 mg
61059	QL9 Peptide	QLSPFPFDL	1063.2	1 mg
62749	Retinal S-Antigen (141-160)	GFLGELTSSEVATEVPFRLM	2183.5	1 mg
61625	RNase (90-105)	SKYPNCAYKTTQANKH	1854.1	1 mg
61043	Smcy HY Peptide (738-746)	KCSRNRQYL	1167.4	1 mg
62746	SmD1 (83-119)	VEPKVKSKKREAV- AGRGRGRGRGRGRGRGRGGPRR	4011.6	1 mg
61058	TRP-2 (180-188)	SVYDFVWL	1175.4	1 mg

Microbial Peptides

62845	PvMSP-1 Peptide 4 (78-97)	NFVGKFLLEQIPGHTDLLH	2178.5	1 mg
62846	PvMSP-1 Peptide 6 (118-137)	FNQLMHVINFHYDLLRANVH	2481.9	1 mg
62847	PvMSP-1 Peptide 8 (158-177)	LDMLKKVVVLGLWKPLDNIKD	2338.9	1 mg
62841	PvMSP-1 peptide 19 (378-397)	LEYLREKAKMAGTLIIPES	2325.8	1 mg
62844	PvMSP-1 Peptide 53 (1058-1077)	SKDQIKLTSLNKLERRQN	2427.8	1 mg
61351	SP1	GFRLTNFGYFEPG	1504.7	1 mg

Myelin Oligodendrocyte Glycoprotein (MOG) / Myelin Basic Protein (MBP) Fragments

62730	MBP (1-20)	ASQKRPSQRSKYLATASTMD	2226.5	1 mg
62734	MBP (14-33)	KYLATASTMDHARHGFLPRH	2309.7	1 mg
62082	MBP (64-80)	ARTTHYGSLPQKSQRSQ	1945.1	1 mg
62081	MBP (69-85)	YGSLPQKSQRSQDENPV	1933.1	1 mg
62085	MBP (69-88)	YGSLPQKSQRSQDENPVVHF	2316.5	1 mg
62733	MBP (79-87)	DENPVVHFF	1103.2	1 mg
62731	MBP (84-97)	VVHFFKNIVTPRTP	1655	1 mg
62739	MBP (84-105)	VVHFFKNIVTPRTPPPSQGKGR	2462.9	1 mg
62729	MBP (92-111)	VTPRTPPPSQGKGRGLSLSR	2091.4	1 mg
62742	MBP (111-129)	LSRFSWGAEGQRPGFGYGG	2029.2	1 mg
62598	MBP (131-155)	ASDYKSAHKGLKGVDAQGTLSKIFK	2650	1 mg
62591	MBP (146-170), Myelin Basic Protein (146-170)	AQGTLKIFKLGGRDSRSGSPMARR	2677.1	1 mg
62369	MOG (8-21)	PGYPIRALVGDEAE	1486.7	1 mg
61605	MOG (35-51)	MEVGWYRSPFSRVVHLY	2126.5	1 mg
61603	MOG (35-52)	MEVGWYRSPFSRVVHLYR	2282.7	1 mg
61604	MOG (35-53)	MEVGWYRSPFSRVVHLYRN	2396.8	1 mg
62737	MOG (37-54), mouse, rat	VGWYRSPFSRVVHLYRNG	2193.5	1 mg
61613	MOG (38-53)	GWYRSPFSRVVHLYRN	2037.3	1 mg
61606	MOG (38-55)	GWYRSPFSRVVHLYRNGK	2222.6	1 mg
61626	MOG (40-54)	YRSPFSRVVHLYRNG	1851.1	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61607	MOG (40-55)	YRSPFSRVVHLYRNGK	1979.3	1 mg
61608	MOG (41-54)	RSPFSRVVHLYRNG	1687.9	1 mg
61609	MOG (42-54)	SPFSRVVHLYRNG	1531.7	1 mg
61610	MOG (43-54)	PFSRVVHLYRNG	1444.7	1 mg
61620	MOG (44-54)	FSRVVHLYRNG	1347.6	1 mg
61611	MOG (45-54)	SRVVHLYRNG	1200.4	1 mg
61612	MOG (46-54)	RVVHLYRNG	1113.3	1 mg
62735	MOG (91-108), rat	SDEGGYTCFFRDHSYQEE	2170.2	1 mg
62738	MOG (91-114), rat	SDEGGYTCFFRDHSYQEEAAVELK	2782	1 mg
62732	MOG (92-106)	DEGGYTCFFRDHSYQ	1824.9	1 mg
62736	MOG (101-108), rat	RDHSYQEE	1063.1	1 mg
62370	MOG (101-120), human, mouse	RDHSYQEEAAMELKVEDPFY	2457.7	1 mg
62083	Peripheral Myelin Protein (P2), (57-81)	FKNTEISFKLQGFEETTADNRKTK	2815.1	1 mg

Myosin Peptides

62126	α -Myosin (614-643)	Ac-SLKLMATLFSTYASADTGDSGKGGKGGK-KKG	3047.5	1 mg
62127	α -Myosin (614-643), Modified	Ac-SLKLLSNLFANYAGADTAPVDKGGK-KAKKG	3105.6	1 mg
62124	α -Myosin (735-748)	GQFIDSGKAGAEKL	1420.6	1 mg
62125	α -Myosin (947-960)	DECSELKKDIDDLE	1651.8	1 mg
62554	Myosin H Chain Fragment, mouse	Ac-RSLKLMATLFSTYASADR	2073.4	1 mg

Neuropeptides

62797	Achatin-1	GfAD	408.4	1 mg
62793	Achatina Cardio-Excitatory Peptide-1 (ACEP-1)	SGQSWRPQGRF-NH ₂	1304.4	1 mg
61858	Activity-Dependent Neurotrophic Factor (1-14), ADNF (1-14)	VLGGGSALLRSIPA	1310.6	1 mg
62094	Adrenal Medulla Chromaffin Vesicles Derived Peptide, (ARP)	ARPVKEP	795.9	1 mg
61531	Alytesin	pE-GRLGTQWAVGHLM-NH ₂	1535.8	1 mg
62795	Aplysia Pedal Peptide	PLDSVYGTHGMSGFA	1538.7	1 mg
61647	Brain Neuropeptide I	AGEGLSSPFWSLAAPQRF-NH ₂	1920.2	1 mg
61648	Brain Neuropeptide II	FLFQPPRF-NH ₂	1050.3	1 mg
62956	Calcitonin Gene Related Peptide, CGRP (8-37), rat	VTHRLAGLLSRSGGVKDNFVPTNVGSEAF-NH ₂	3127.6	1 mg
62433	Catestatin, Chromogranin A (352-372), human	SSMKLSFRARAYGFRGPGPQ	2213.6	1 mg
62513	Catestatin, human	SSMKLSFRARAYGFRGPGPL	2198.6	1 mg
62707	CART Protein Precursor (82-86)	IPIYE	633.8	1 mg
62661	Corticotropin-Lipotropin Intermediary Peptide, CLIP (103-120)	AEEETAGDGRPEPSPRE-NH ₂	1882.9	1 mg
62173	Claudin-like Protein 24, (CLP24), (155-169), rat	TFYRIGPYTNLSWSC	1808.1	1 mg
62172	Claudin-like Protein 24, (CLP24), (187-201), human	IWNILHKREDCMAPR	1882.3	1 mg
62171	Claudin-like Protein 24, (CLP24), (212-226), rat	RFRRGLDNDYVESPC	1827.0	1 mg
62801	Ecdysis-Triggering Hormone 1 (ETH1), <i>Drosophila</i>	DDSSPGFFLKITKNVPRL-NH ₂	2033.4	1 mg
62803	Ecdysis-Triggering Hormone AP (ETH-AP) (1-12), <i>Bombyx mori</i>	NYDSGNHFDIPK	1406.5	1 mg
62802	Ecdysis-Triggering Hormone-AP (ETH-AP) Fragment, <i>Manduca sexta</i>	NYDSENRFDIPK	1497.6	1 mg
62562	Erythropoietin, Human (hEPO) Fragment	MEVQQQAVEVWQGLALLSEAVLR	2527	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62548	Glutamate Receptor Endocytosis Inhibitor, Control Peptide (GluR23A)	AKEGANVAG	815.9	1 mg
62547	Glutamate Receptor Endocytosis Inhibitor, GluR23Y	YKEGYNVYG	1092.2	1 mg
62671	Gonadoliberin I (24-33)	QHWSYGLRPG	1200.3	1 mg
61907	Hym-33H, <i>Hydra</i> Nerve Cell Differentiation Inhibitor	AALPW	556.7	1 mg
61908	Hym-54, <i>Hydra</i> Signal Muscle Contraction Molecule	GPMTGLW-NH ₂	759.9	1 mg
61909	Hym-301, <i>Hydra</i> Axial Patterning Peptide	KPPRRCYLNGYCSP-NH ₂ (Disulfide bridge: 6-12)	1651.0	1 mg
61868-05	Hypocretin-2 (69-97), human	RPGPPGLQGRLQRLQLQANGNHAAGILTMG	2994.5	1 mg
62799	Leucokinin 3	DQGFNSWG-NH ₂	908.9	1 mg
62772	Leucokinin I, <i>Leucophaea maderae</i>	DPAFNSWG-NH ₂	890.9	1 mg
62767	Leucokinin I, Myokinin I, <i>Anopheles</i>	DTPRYVSKQKFHSWG-NH ₂	1835.1	1 mg
62771	Leucokinin II, <i>Leucophaea maderae</i>	DPGFSSWG-NH ₂	850.9	1 mg
62768	Leucokinin II, Myokinin II, <i>Anopheles</i>	NPFHSWG-NH ₂	842.9	1 mg
62769	Leucokinin III, Myokinin III, <i>Anopheles</i>	NTAQVFYPWG-NH ₂	1181.3	1 mg
62770	Leucokinin, <i>Drosophila</i>	NSVLGKKQRFHSWG-NH ₂	1742	1 mg
61291	Leuprolide acetate, truncated	Y-dL-LRP-NHEt	687.8	1 mg
62773	Locustakinin I	AFSSWG-NH ₂	652.7	1 mg
62796	Myomodulin	PMSMLRL-NH ₂	846.1	1 mg
61436	Neuromedin S, human	ILQRGSGTAAVDFTKKDHTATWGRPFFLFR- PRN-NH ₂	3791.4	1 mg
62821	Neuropeptide-FF	FLFQPQRF-NH ₂	1081.3	1 mg
61653	Neuropeptide NPW-23, human	WYKHVASPRYHTVGRAAGLLMGL	2584.1	1 mg
61642	Neuropeptide NPW-30, human	WYKHVASPRYHTVGRAAGLLMGLRRSPYLW	3543.2	1 mg
61852	Neuropeptide W-23 (NPW23), porcine	WYKHTASPRYHTVGRAAGLLMGL	2566.0	1 mg
61853	Neuropeptide W-23 (NPW23), rat	WYKHVASPRYHTVGRASGLLMGL	2600.1	1 mg
61870-05	Neuropeptide W-30 (NPW30), rat	WYKHVASPRYHTVGRASGLLMGLRRSPYLW	3559.2	0.5 mg
62599	Neuropeptide S, NPS (1-10), rat	SFRNGVGSGV	979.1	1 mg
62608	Neuropeptide S, hNPS (1-12), human	SFRNGVGTGMKK	1281.5	1 mg
62618	Neuropeptide S, NPS, chicken	SFRNGVGSGIKKTSFRRAKS	2183.5	1 mg
62617	Neuropeptide S, NPS, chimpanzee, dog	SFRNGVGTGMKKSFRRAKS	2215.6	1 mg
62016	NPSF, Neuropeptide SF (1-37)	SLNFEELKDWGPKNVIKMSTPAVNKMPHS- FANLPLRF-NH ₂	4257.0	1 mg
62006	NPVF, Neuropeptide VF (1-37)	VPNLPQRF-NH ₂	969.2	1 mg
61936	Ac-[Leu ^{28,31}]-Neuropeptide Y (24-36)	Ac-LRHLYNLLTRQRY-NH ₂	1787.1	1 mg
62677	Neurotensin	QLYENKPRRPYIL	1690	1 mg
62928	[Trp ⁵]-Oryzatesin (5-9), rice	WPLPR	667.8	1 mg
62774	Pevpyrokinin-1 (PPK-1, Pev-PK 1)	DFAFSPRL-NH ₂	951.1	1 mg
62775	Pevpyrokinin-2 (PPK-2, Pev-PK 2)	ADFAFSPRL-NH ₂	1022.2	1 mg
62675	Pituitary Adenylate Cyclase-Activating Polypeptide Precursor (111-128)	GMGENLAAAVIDDRAPLT	1772	1 mg
62804	Pre-Ecdysis-Triggering Hormone (PETH)	SFIKPNVPRV-NH ₂	1269.5	1 mg
62419	Rhodopsin (338-348), human, mouse, rat, dog, bovine, pig	KTETSQVAPA	1031.1	1 mg
62662	Secretogranin I Precursor (585-594)	SFAKAPHLDL	1098.3	1 mg
62706	Secretogranin-5, C-Terminal Peptide (198-210), rat	SVPHFSEEEKEPE	1543.6	1 mg
62673	Secretoneurin	TNEIVEEQYTPQSLATLESVFQELGKLTGP- SNQ	3652	1 mg
62787	Δ-Sleep-Inducing Peptide (DSIP)	WAGGDASGE	848.8	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62798	Small Cardioactive Peptide B (SCPB)	MNYLAFPRM-NH ₂	1141.4	1 mg
62674	SMR1-Related Undecapeptide	VRGPRRQHNPR	1372.6	1 mg
62823	Tyr-K-MIF-1	YPKG-NH ₂	462.6	1 mg
62822	Tyr-MIF-1	YPLG-NH ₂	447.5	1 mg
62703	VEGF Protein Precursor (491-507)	PPEPVPPPRAAPATHV	1730	1 mg
62672	WE-14	WSRMDQLAKELTAE	1677.9	1 mg

NF-κB/Transcription Factors Related Peptides

62116	EBIP37, Biotin-LC, Nuclear Receptor-Interacting LXXLL Peptide	Biotin-LC-TGGGVSLLLHLLNTEQGES	2264.6	1 mg
61445	FACTP140 (838-845), Chromatin-Specific Transcription Elongation Factor (838-845)	DEVELIHF	1001.1	1 mg
61671	IκBα, Nuclear Factor κB (20-39), human	LKKERLLDDRHSGLDSMKD	2371.7	1 mg
61444	SOAT1/ACAT1 (163 - 171)	DEGRLVLEF	1077.2	1 mg

Nuclear Export Signal Peptides

61885	NES, Nuclear Export Signal, Adenoviral E1A Protein	VMLAVQEGIDL	1187.4	1 mg
61886	NES, Nuclear Export Signal, Dok1 (348-359)	LLKAKLTDPKED	1370.6	1 mg
61890	NES, Nuclear Export Signal, HIV Rev	LQLPPLERLTLTD	1407.7	1 mg
61888	NES, Nuclear Export Signal, MAPKK	ALQKKLEEELELD	1428.7	1 mg
61884	NES, Nuclear Export Signal, MDM2	SLSFDESLALC	1184.3	1 mg
61889	NES, Nuclear Export Signal, Nmd3p (491-500)	INIDELLDEL	1186.3	1 mg
61883	NES, Nuclear Export Signal, p53	FRELNEALELKD	1476.7	1 mg
61878	NES, Nuclear Export Signal, p120 Catenin Fragment	CSLEEEELDVLVLDDEGG	1835.0	1 mg
61887	NES, Nuclear Export Signal, PKI	ELALKLAGLDIN	1269.5	1 mg
61897	NES, Nuclear Export Signal, Topoisomerase II α (1017-1028)	DILRDFELRLK	1564.9	1 mg

Opioid Related Peptides

62653	Lactoferrin C, Lactoferrin (673-679)	KYLG PQY	868	1 mg
62790	LVV-Hemorphin-6, Leu-Valorphin-Arg	LVVYPWTQR	1161.4	1 mg
62789	Valorphin	VVYPWTQ	892	1 mg
62611	Nociceptin (1-7), Orphanin FQ, or OFQ/N (1-7)	FGGFTGA	655.7	1 mg
62610	Nociceptin (1-11), Orphanin FQ, or OFQ/N (1-11)	FGGFTGARKSA	1098.2	1 mg
61950-5	Nociceptin (1-13), amide	FGGFTGARKSARK-NH ₂	1381.6	5 mg
62609	Orphanin FQ2, (OFQ2, NOCII)	FSEFMRQYLVLSMQSSQ	2081.4	1 mg
62612	Prepronociceptin (169-176), human	TLHQNGNV	882	1 mg

OVA Peptides

61659	β-catenin Peptide, βCATP	RTYTYEKL	1073.2	1 mg
62194	[Glu ²⁵⁷]-OVA (257-264), E1	EIINFEKL	1005.2	1 mg
62573-1	MSL8, [Met ²⁵⁶]-OVA (257-264)	MSIINFEKL	1094.4	1 mg
62573-5	MSL8, [Met ²⁵⁶]-OVA (257-264)	MSIINFEKL	1094.4	5 mg
62577	OVA (327-333)	VHAAHAE	733.8	1 mg
62571-1	OVA (324-339), amide	SQAVHAAHAEINEAGR-NH ₂	1773	1 mg
62571-5	OVA (324-339), amide	SQAVHAAHAEINEAGR-NH ₂	1773	5 mg
62572-1	OVA (257-264), amide	SIINFEKL-NH ₂	962.2	1 mg
62572-5	OVA (257-264), amide	SIINFEKL-NH ₂	962.2	5 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
Oxytocins, Vasopressin and Related Peptides				
62097	3,4 Dehydro-Pro-Arg-Vasotocin	CYIQNC-(3,4-Dehydro-P)-RG-NH ₂ (Disulfide bridge: 1-6)	1048.3	1 mg
62007	Aquaporin 2 (AQP2), (241-271)	EPDTDWEEREVRRRQSVLHSPQSLPRG-SKA	3676.0	1 mg
62535	Serum Albumin (202-226)	ADDKETCFEEGKLVAAASQAALGL	2565.9	1 mg
62664	Vasopressin-Neurophysin 2-Copeptin Precursor (151-168)	VQLAGTQESVDSAKPRVY	1948.2	1 mg
p53 Peptides				
61374	p53 (12-20)	PPLSQETFS	1005.1	1 mg
61373	p53 (12-26)	PPLSQETFSDLWKLL	1774.1	1 mg
61375	p53 (17-26)	ETFSDLWKLL	1251.4	1 mg
62386	p53 (17-26), FITC-LC-labeled	FITC-LC-ETFSDLWKLL-NH ₂	1753.1	1 mg
62122	p53 Mutant Form (361-371), Pab 421	KKGQSTSRHKK-NH ₂	1283.5	1 mg
62121	[Lys(Ac) ³⁷¹]-p53 Mutant Form (361-377), TP53 Q9NP68	KKGQSTSRHK-K(Ac)-LMFKTEG	2133.5	1 mg
62120	p53 Mutant Form (361-377), TP53 Q9NP68 (361-377)	KKGQSTSRHKKLMFKTEG	2091.5	1 mg
62529	p53 Tumor Suppressor (361-393), human	GSLRAHSSHLKSKKGQSTSRHKKLMFKTEG-PDSD	3653.1	1 mg
62530	p53 Tumor Suppressor (361-393), Biotin-LC, human	Biotin-LC-GSLRAHSSHLKSKKGQSTSRHKKLM-FKTEGPDSD-NH ₂	3992.6	0.5 mg
Pancreatic Polypeptides				
62516	Apolipoprotein J (215-222)	RPHFFFPK	1075.3	1 mg
62517	Apolipoprotein L (306-316)	VNEPSILEMSR	1274.5	1 mg
62405	Insulin α-chain (85-99)	KRGIVEQCCTSICSL	1640	1 mg
63773	Insulin β Chain (15-23)	LYLVCGERG	1009.2	1 mg
61532	Insulin B (9-23)	SHLVEALYLVCGERG	1645.9	1 mg
62645	Insulin B (13-23)	EALYLVCGERG	1209.4	1 mg
62431	Preptin, Insulin Like Growth Factor 2 (67-100), mouse	DVSTSQAVLPDDFPRYPVVGKFFQYDT-WRQSAGRL	3948.4	1 mg
62432	Preptin, Insulin Like Growth Factor 2 (67-100), rat	DVSTSQAVLPDDFPRYPVVGKFFKFDT-WRQSAGRL	3932.4	1 mg
62430	Preptin, Insulin Like Growth Factor 2 (96-129), human	DVSTPPTVLPDNFPRYPVVGKFFQYDT-WKQSTQRL	4029.5	1 mg
62545	Insulin Receptor Substrate 1 (IRS-1) (979-989), Peptide-983	SRGDYMTMQIG	1258.5	1 mg
62544	CKK-Insulin Receptor Substrate 1 (IRS-1) (979-989), mouse	CKKSRGDYMTMQIG	1617.9	1 mg
62092	Pancreatic Polypeptide (30-53), human	APLEPVYPGDNATPEQMAQYAADL	2561.8	1 mg
Parathyroid Hormone Related Peptides				
62093	Parathyroid Hormone-related Protein (PTHrP), (107-111)	TRSAW	619.7	1 mg
62286	[pSer ⁸⁸⁴]-Thyroid Hormone Binding Protein (TRBP), Biotin-labeled	Biotin-KDVTLT(pSer)PLLVNLLQSDIS	2362.7	1 mg
Peptide Standards				
61220	ClearPoint™ BSA (347-359), Isotopic labeled, Mass Spec Standard	DAF-L*-GSF-L*-YEYSR [L* = L(U- ¹³ C ₆ , ¹⁵ N)]	1581.7	0.1 mg
61219	ClearPoint™ C-Reactive Protein-Like Peptide, Isotopic labeled, Mass Spec Standard	GYSI-F*-SYATK [F* = F(U- ¹³ C ₉ , ¹⁵ N)]	1146.3	0.1 mg
Phytochelatin Peptides				
61843	Homo-Phytochelatin-2	(γE-C) ₂ -β-A	553.6	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61842	Hydroxymethyl-Phytochelatin 2	(γ E-C) ₂ -S	569.6	1 mg
61844	Iso-Phytochelatin-2	(γ E-C) ₂ -E	611.7	1 mg
61795	Phytochelatin 2, PC2, Biotin-labeled	Biotin-LC-(γ E-C) ₂ -G	879.1	1 mg
61839	Phytochelatin 2, PC2, HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-LC-(γ E-C) ₂ -G	1009.2	0.5 mg
61796	Phytochelatin 3, PC3, Biotin-labeled	Biotin-LC-(γ E-C) ₃ -G	1111.4	1 mg
61840	Phytochelatin 3, PC3, HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-LC-(γ E-C) ₃ -G	1241.5	0.5 mg
61797	Phytochelatin 4, PC4, Biotin-labeled	Biotin-LC-(γ E-C) ₄ -G	1343.6	1 mg
61841	Phytochelatin 4, PC4, HiLyte Fluor™ 488-labeled	HiLyte Fluor™ 488-LC-(γ E-C) ₄ -G	1473.7	0.5 mg

Prion Protein (PrP) Fragments

62112	Prion Protein, (PrP), (117-136), hamster	AAGAVVGGGLGGYMLGSAMSR	1825.2	1 mg
62110	Prion Protein, (PrP), (166-179), hamster	VDQYNNQNNFVHDC	1709.8	1 mg
62111	Prion Protein, (PrP), (166-179), hamster, Biotin-LC-labeled	Biotin-LC-VDQYNNQNNFVHDC	2049.3	1 mg
62109	Prion Protein, (PrP), (200-223), hamster	ETDIKIMERVVEQMCTTQYQKESQ	2918.3	1 mg

Protease Inhibitors

61832-5	Apstatin, Aminopeptidase P Inhibitor	N-[(2S,3R)-3-amino-2-hydroxy-4-phenylbutanoyl]-PPA-NH ₂	459.6	5 mg
61832-25	Apstatin, Aminopeptidase P Inhibitor	N-[(2S,3R)-3-amino-2-hydroxy-4-phenylbutanoyl]-PPA-NH ₂	459.6	25 mg

Protein Structure / Folding Related Peptides

62563	Elution (EY) Peptide	EYMPD	754.8	5 mg
62311	GALA, Pore-Forming Peptide	WEAALAEALAEALAEHLAEALAEALAEALAA	3032.4	1 mg
62532	β -hairpin (BHA), Protein G B1 Domain (41-56)	GEWTYDDATKTFTVTE	1864	1 mg
62533	[Cys40]- β -hairpin (BHA), Protein G B1 Domain (41-56)	CGEWTYDDATKTFTVTE	1967.1	1 mg

Protein Phosphorylation Related Peptides

61284	CREB2 (91-116)	ESEDSQESVDSVTDSQKRREILSRRP	3034.2	1 mg
61809	Erythrocyte G-protein, α -subunit, FAM-labeled	5-FAM-DLLRCRVLTS	1590.8	1 mg
61242	DAM1 (221-241) Peptide	SFVLNPTNIGM-pS-KSSQGHVTK	2312.6	1 mg
62510	Dominant H2-Kd Ligand	SYFPEITHI	1106.3	1 mg
62346	Erythropoietin Receptor Peptide (EpoR), pTyr429, acetylated, amide	Ac-DPPHLK-pY-LYLVVSDSK-NH ₂	1995.2	1 mg
62744	[GalNAc-Ser]-Erythropoietin (Epo) (117-131)	EAISPPDAA-*S-AAPLR (*S=GalNAc-Ser)	1668.6	0.5 mg
61528	HIF-1 α (556-574), Transcription Factor	DLGLEMLAPYIPMDDDFQL	2254.6	1 mg
61529	HIF-2 (66-84), Transcription Factor	SLEAQGIKADRETVAVKPT	2013.3	1 mg
62594	HSP-B8, Heat Shock Protein-B8 (10-29)	CHYPSRLRRDPFRDPSLSSR	2445.8	1 mg
62593	HSP-B8, Heat Shock Protein-B8 (181-194)	SFNNELPQDSQEV	1607.7	1 mg
62166	Interferon α A (101-110), FRET Peptide	GVGVTETPLME(EDANS)GVGVETETPLM-K(DABCYL)K	2874.4	1 mg
62713	[Asp ⁴⁵] p21 (140-147) Phosphomimic Peptide	RKRRQDSM	1076.3	1 mg
61737	PKCtide, Phosphorylated, FAM-labeled	5-FAM-RFARKG-pS-LRQKNV-NH ₂	1997.2	1 mg
62809	Phospholemman, PLM (58-72)	GTFRSSIRRLSTRRR	1849.1	1 mg
62185	δ -PKC Agonist, MRA	MRAAEDPM	920.1	1 mg
62176	ϵ PKC, C2 Domain, ϵ -C2/V1 (4-12)	FNGLLKIKI	1045.3	1 mg
62179	ϵ PKC, C2 Domain, ϵ -C2/V1 (28-35)	AVGPRPQT	824.9	1 mg
62178	ϵ PKC, C2 Domain, ϵ -C2/V1 (36-41)	FLLDPY	766.9	1 mg
62175	ϵ PKC, C2 Domain, ϵ -C2/V1 (49-53)	SRIGQ	559.6	1 mg
62184	ϵ PKC, C2 Domain, ϵ -C2/V1 (62-67)	SPAWHD	711.7	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62183	ε PKC, C2 Domain, ε-C2/V1 (68-73)	EFVTDV	708.8	1 mg
62174	ε PKC, C2 Domain, ε-C2/V1 (74-79)	CNGRKI	689.8	1 mg
62182	ε PKC, C2 Domain, ε-C2/V1 (79-84)	IELAVF	690.8	1 mg
62181	ε PKC, C2 Domain, ε-C2/V1 (112-118)	HFEDWID	961.0	1 mg
62177	ε PKC, C2 Domain, ε-C2/V1 (119-124)	LEPEGK	671.8	1 mg
62180	ε PKC, C2 Domain, ε-C2/V1 (125-131)	VYVIIDL	834.0	1 mg
62188	ε PKC, C2 Domain, ψ-RACK, ε-C2/V1 (82-92)	HDAPIGYD	886.9	1 mg
62295	ε-V1-2, ε-PKC Specific Inhibitor	CEAVSLKPT	947.1	1 mg
62187	ε-PKC Inhibitor, Cys conjugated, ε-V1-2	EAVSLKPTC	947.1	1 mg
62186	ε-PKC Inhibitor, ε-V1-2	EAVSLKPT	844.0	1 mg
62345	[pTyr ⁴²⁹]-Erythropoietin Receptor Peptide (EpoR), (423-438)	DPPHLK-pY-LYLVVSDSK	1954.2	1 mg
62160	R18, 14-3-3/Raf-1 Kinase Binding Inhibitor	PHCVPRDLSWLDLEANMCLP	2309.7	1 mg
62637	S3 Fragment, ADF/cofilin	MASGVMVSDDVVKVFN	1698	1 mg
61740	Glycogen Synthase Kinase 3 (GSK-3), FAM-labeled	5-FAM-RRAAEELDSRAG-pS-PQL-NH ₂	2193.2	1 mg
62287	TRBP, Thyroid Hormone (TR)-Binding Protein, Biotin-labeled	Biotin-KDVTLTSPLLVNLQSDIS	2282.7	1 mg
62286	[pSer ⁸⁸⁴]-TRBP, Thyroid Hormone (TR)-Binding Protein, Biotin-labeled	Biotin-KDVTLT-pS-PLLVNLQSDIS	2362.7	1 mg

Proteolipid Protein (PLP) Fragments

62740	PLP (57-70)	YEYLINVIHAFQYV	1772	1 mg
62741	PLP (178-191)	NTWTTCQSIAPFSK	1583.8	1 mg

Salusin Peptides

61282	Salusin-α	SGALPPAPAAPRPALRAQRAGPAGPGAK-NH ₂	2603.0	1 mg
61283	Salusin-β	AIFIFIRWLLKLGHHGRAPP	2342.9	1 mg

Signal Transduction Peptides

62191	60S Ribosomal Epitope, L3 (162-170), human	NRAVVGVA	884.1	1 mg
63315	Bacterial Trypsin Fragment	TCQGDSGGPMFR	1255.4	1 mg
62524	bhpBR3	CHWDLLVRHWVC (Disulfide bridge: 1-12)	1564.9	1 mg
62910	C3f fragment, Human c3 (1286-1297)	THRIHWESASLL	1449.6	1 mg
62909	C3f, Human c3 (1282-1298)	SSKITHRIHWESASLLR	2021.3	1 mg
63775	Cardiac Troponin I (41-49)	SASRKLQLK	1030.2	1 mg
63753	CaMKI, Calmodulin-Dependent Protein Kinase I (299-320) Binding Domain	AKSKWKQAFNATAVVRHMRKIQ	2599.1	1 mg
62255	Caveolin-1 Scaffolding Domain (82-101)	DGIWKASFTTFTVTKYWFYR	2517.9	1 mg
62019	Caveolin-3 (Cav-3), (55-74)	DGVWRVSYTFTVSKYWCYR	2517.9	1 mg
63316	Chymotrypsin Bovine A Fragment	SCMGDSGGPLVC	1125.3	1 mg
62816	CLA-1	CTSAPKGSVLQEAKL	1531.8	1 mg
62817	CLA-2	CLPDSPSRQPPSPTA	1552.7	1 mg
61446	CUGBP1, RNA CUG Triplet Repeat Binding Protein	TENDIRVMF	1124.3	1 mg
62251	[Cys ¹] Inter-α-Trypsin Inhibitor Heavy Chain 4 (668-681), C-terminal fragment (H4CT)	CLGLPGPPDVPDH	1459.7	1 mg
62250	[Cys ¹] Inter-α-Trypsin Inhibitor Heavy Chain 4, ITIH4 (657-687), C-terminal fragment 1, H4CT	CMNFRPGVLSSRQLGLPGPPDVPDHAAYHPF	3376.9	1 mg
62698	Dyrktide	RRRFRPASPLRGPPK	1791.2	1 mg
62716	Dyrktide, FAM-labeled	5-FAM-RRRFRPASPLRGPPK	2149.5	1 mg
62715	Dyrktide, TAMRA-labeled	5-TAMRA-RRRFRPASPLRGPPK	2203.6	1 mg
63314	Elastase Fragment	GCQGDSGGPLHC	1130.2	1 mg
62234	eNOS (493-512), bovine	TRKKTfKEVANAVKISASLM	2222.7	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62372	Eukaryotic Translation Initiation Factor, 4E b	PGGTRIIYDRKFLMECRNSP	2353.8	1 mg
61739	[pSer ^{8,13}]-Eukaryotic Translation Initiation Factor 2B, eIF2B	RRAAEELD-pS-RAG-pS-PQL	1915.9	1 mg
62123	[pSer ^{8,13}]-Eukaryotic Translation Initiation Factor 2B, eIF2B, FAM-labeled	5-FAM-RRAAEELD-pS-RAG-pS-PQL-NH ₂	2273.2	1 mg
61875	Generic Protein Kinase Substrate	RARTLSFAEPG	1204.4	1 mg
63722	GS2, Glycogen Synthase Peptide-2	YRRAAVPPSPSLSRHSSPHQ-pS-EDEEE	3029.1	1 mg
62048	Human Platelet Membrane Glycoprotein Binding Site, GPIIb (580-590)	LVAGRRPSALS	1126.3	1 mg
62252	Inter- α -trypsin inhibitor heavy chain 4, C-terminal fragment (H4CT)	CLGLPGPPDVPDH	1316.5	1 mg
61298	JAG-1 (188-204), Jagged-1 (188-204), Notch Ligand	CDDYYYGFGCNKFCRPR	2107.4	1 mg
61627	JNK1(180-189)	F-Nle-Nle-pT-P-pY-VVTR	1368.5	1 mg
61628	(pTyr ¹⁸⁵) MAP (177-189) Kinase	DHTGFLTE-pY-VATR	1589.6	1 mg
62001	MKK4 Docking Site	QGKRKALKLNF	1302.6	1 mg
62254	iNOS (507-531), human	RPKRREIPLKVLVKAVLFACMLMRK	2996.9	1 mg
62253	nNOS, Neural Nitric Oxide Synthase (725-747), rat	KRRAIGFKKLAEAVKFSAKLMGQ	2578.2	1 mg
62914	Notch 1 (1735-1752)	VLLFFVGCGLLSRKR	2119.7	1 mg
62534	NR-Box 2, Nuclear Receptor-box 2 Peptide	KHKILHRLQDSS	1574.9	1 mg
62628	P1pal-19	Pal-RCLSSSAVANRSKSRALF-NH ₂	2318.9	1 mg
61903	Pannexin-1 Fragment (4512)	VQQKSSLQSESGN	1391.5	1 mg
61904	Pannexin-1 Fragment (4515)	EKNSRQRLNPS	1441.6	1 mg
61911	Pannexin-1 (Panx1) Mimetic Blocking Peptide	WRQAAFVDSY	1242.4	1 mg
61292	PGC-1 α , Peroxisome Proliferator-activated Receptor Coactivator-1 α (205-216)	RPCSELLKYLTT	1423.7	1 mg
61293	Ac-PGC-1 α , Peroxisome Proliferator-activated Receptor Coactivator-1 α (205-216), amide	Ac-RPCSELLKYLTT-NH ₂	1464.8	1 mg
62025	PKC ζ 410	GDTTSTFCGTPN	1200.3	1 mg
62020	PKC ζ 560	EPVQLTPDDEDA	1328.4	1 mg
62376	Protein A Specific Ligand	GHHKHHGGG	923	1 mg
63755	Protein-Tyrosine Phosphatase-PEST-WT	PLSFTNPLHSDDWH	1665.8	1 mg
62062	Receptor Interacting Protein 140, RIP 140	SFSKNGLLSRLLRQNQDSY	2226.5	1 mg
61995	Receptor Interacting Protein, (RIP), (131-142), L2	TLLASLLQSESS	1248.4	1 mg
61997	Receptor Interacting Protein, (RIP), (498-509), L6	TLLQLLLGHKNE	1378.6	1 mg
61996	Receptor Interacting Protein, (RIP), (817-828), L8	GLLSRLLRQNQD	1412.6	1 mg
62018	RI Anchoring Disruptor, (RIAD), FAM-labeled	LEQYANQLADQIIKEATEK(FAM)-NH ₂	2563.8	1 mg
62000	Signal Transduction/Transcription Factor (STTF) (681-698)	TSLKEKHKILHRLQDSS	2133.5	1 mg
62676	Stathmin (2-20)	ASSDIQVKELEKRASGQAF	2064.3	1 mg
62383	TET 830 modified/T-helper Epitope, FAM-labeled	5-FAM-AQYIKANSKFIGITEL	2154.4	1 mg
61993	Transcriptional Intermediary Factor 2, (TIF2), Box1, (636-649)	KGQTKLLQLLTTKS	1558.9	1 mg
61994	Transcriptional Intermediary Factor 2, (TIF2), (685-697)	EKHILHRLQDSS	1616.9	1 mg
61992	Transcriptional Intermediary Factor 2, (TIF2), Box3, (740-753)	KENALLRYLLDKDD	1705.9	1 mg
62228	V-ATPase 31, C-terminal	CGANANRKFLD	1208.4	1 mg
62227	V-ATPase a3 (OC-116), (820-830)	LSPFTFAATDD	1184.3	1 mg

Somatostatins

62071	Cortistatin-29 (human)	EGAPPQQSARRDRMPCRNFFWKTFSSCK (Disulfide bridge:16-27)	3329.8	1 mg
62663	Somatostatin-28 (1-12)	SANSNPAMAPRE	1244.4	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
Steroid Receptor Coactivators (SRCs)				
62152	SRC-1, Steroid Receptor Coactivator-1 (676-700), Biotin-labeled	Biotin-CPSSHSSLTERHKILHRLLQEGSPS	3026.5	1 mg
62258	SRC-1, Steroid Receptor Coactivator-1 (676-700)-Lys(Biotin), amide	CPSSHSSLTERHKILHRLLQEGSPS-K(Biotin)-NH ₂	3153.7	1 mg
62151	SRC-1, Steroid Receptor Coactivator-1 (676-701)	CPSSHSSLTERHKILHRLLQEGSPSD	2915.3	1 mg
62257	SRC-1, Steroid Receptor Coactivator-1 (690-695)	LHRLQ	779	1 mg
61998	SRC-1, Steroid Receptor Coactivator-1 (1433-1440)	SLLQQLLT	915.1	1 mg
61999	SRC-1, Steroid Receptor Coactivator-1 (1433-1441)	SLLQQLLTE	1044.2	1 mg
Substance P and Analogs				
62824	Ranakinin	KPNPERFYGLM-NH ₂	1350.6	1 mg
61690	Substance P (1-6), amide	RPKPQQ-NH ₂	751.9	1 mg
61691	Substance P (1-7), amide	RPKPQQF-NH ₂	899.1	1 mg
61692	Substance P (5-11), amide	QQFFGLM-NH ₂	869.1	1 mg
61693	Substance P (7-11), amide	FFGLM-NH ₂	612.8	1 mg
61694	Substance P (8-11), amide	FGLM-NH ₂	465.6	1 mg
61695	Substance P (9-11), amide	GLM-NH ₂	318.4	1 mg
62929	[D-Arg ¹ , D-Phe ⁵ , D-Trp ^{7,9} , Leu ¹¹]-Substance P	rPKPqQwFwLL	1517.9	1 mg
Tachykinin and Related Peptides				
62627	[Lys ⁶]-Eledoisin (6-11)	KFIGLM-NH ₂	707	1 mg
61652	Endokinin A/B	GKASQFFGLM-NH ₂	1084.3	1 mg
61638	Endokinin C, human	KKAYQLEHTFQGLL-NH ₂	1675.0	1 mg
61637	Endokinin D, human	VGAYQLEHTFQGLL-NH ₂	1574.8	1 mg
62792	Hylambatin	DPPDPDRFYGMM-NH ₂	1439.6	1 mg
61933	Invertebrate Tachykinin-I, OctTK-I	KPPSSSEFIGLM-NH ₂	1291.5	1 mg
61932	Invertebrate Tachykinin-II, OctTK-II	KPPSSSEFVGLM-NH ₂	1277.5	1 mg
62670	Preprotachykinin A, C-Terminal Flanking Peptide (111-126)	ALNSVAYERSAMQNYE	1846	1 mg
62779	Ranamargarin	DDASDRAKKFYGLM-NH ₂	1615.8	1 mg
62776	Ranatachykinin A	KPSPDRFYGLM-NH ₂	1309.6	1 mg
62777	Ranatachykinin B	YKSDSFYGLM-NH ₂	1209.4	1 mg
62778	Ranatachykinin C	HNPASFIGLM-NH ₂	1085.3	1 mg
61935	Sialokinin-1	NTGDKFYGLM-NH ₂	1144.3	1 mg
61934	Sialokinin-2	DTGDKFYGLM-NH ₂	1145.3	1 mg
61615	γ-TAC4, γ-Tachykinin Precursor (30-61), amide	TEAETWEGAGPSIQLQLQEVTGKASQFF-GLM-NH ₂	3481.9	1 mg
61614	γ-TAC4, γ-Tachykinin Precursor (32-50)	AETWEGAGPSIQLQLQEVK	2084.3	1 mg
62780	Tachykinin OG1	DDTEDLANKFIGLM-NH ₂	1580.8	1 mg
Tag Peptides				
62189	Glu-Glu Epitope Tag	EYMPME	798.9	5 mg
62229	HA 12CA5 Epitope	CYPYDVPDYA	1205.3	1 mg
62190	TETSQVAPA, Rhodopsin Epitope Tag	TETSQVAPA	903.0	5 mg
Thrombin Related Peptides				
62374	Coagulation factor IX Fragment A	TEAETILDNI	1118.2	1 mg
62925	Coagulation Factor V Fragment	SNETALSPDLNQTSPSMSTDR	2251.4	1 mg
62049	GRGDNP, Integrin-Binding Site	GRGDNP	614.6	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62987	Hirudin (54-65)	GDFEEIPEEYLQ	1468.5	1 mg
62986	Hirudin (54-65), sulfated	Ac-GDFEEIPEE-Y(SO3H)-LQ	1590.6	1 mg
62988	Hirullin P18	SDFEEFSLDDIEQ	1573.6	1 mg
62440	α -2 Plasmin Inhibitor, Serpin Peptidase Inhibitor, Clade F (40-46), human	NQEQVSP	800.8	1 mg
62437	Peptide 2A-54, Coagulation Factor IX (11-30), human	TEQKRNVIIRIIPHHNYNAAI	2387.7	1 mg
61530	Protease-Activated Receptor-1, PAR-1-Selective Peptide	TFLLRN	762.9	1 mg
62936	Protease-Activated Receptor-1, PAR-1 Agonist, amide	TFLLRNPNDK-NH ₂	1216.4	1 mg
62937	Protease-Activated Receptor-1, PAR-1 Agonist, amide	TFLLRN-NH ₂	761.9	5 mg
62657	Protease-Activated Receptor-3 (1-6), PAR-3 (1-6), human	TFRGAP-NH ₂	646.8	1 mg
62938	Protease-Activated Receptor-3, PAR-3 Agonist, amide	SFNGGP-NH ₂	576.6	1 mg
63368	Protein C (HPC) Epitope	EDQVDPRLIDGK	1384.5	1 mg
24190-25	Thrombin Receptor Activator for Peptide 6 (TRAP-6)	SFLLRN	748.9	25 mg

TNF Peptides

62098	TRAF - Binding Motif, acetylated, amide	Ac-PVQET-NH ₂	613.7	1 mg
62679	Tumor Necrosis Factor- α Converting Enzyme, TACE (807-823), human	ASFKLQRQNRVDSKETE	2036.2	1 mg
62619	Tumor Necrosis Factor Receptor, TNFR (159-178) Analog	QEKQNTVATAHGFFLRENEG	2276.5	1 mg

Toxins

61646	Agelenin	GGCLPHNRFNCNALSGPRCCSGLKCKELSI-WDSRCL-NH ₂ (Disulfide bridges: 3-19, 10-24, and 18-34)	3818.6	0.1 mg
62506	Bombolitin I	IKITTMLAKLGKVLAVH	1836.4	1 mg
62874	Bombolitin II	SKITDILAKLGKVLAVH	1806.2	1 mg
62875	Bombolitin III	IKIMDILAKLGKVLAVH	1862.4	1 mg
62876	Bombolitin IV	INIKDILAKLVKVLGHV	1873.4	1 mg
62605	Caloxin 1A1	ACPWWSPHACGGG (Disulfide bridge: 2-10)	1326.5	1 mg
62878	Brevinin-2Eg	GIMDTLKNLAKTAGKQALQSLLNHASCKLS-GQC (Disulfide bridge 27-33)	3371	1 mg
62879	Brevinin-2Eh	GIMDTLKNLAKTAG-KGALQSLLNHASCKLSKQC (Disulfide bridge 27-33)	3442.1	1 mg
62604	Caloxin 2A1	VSNSNWSPFPSSGGG	1479.5	1 mg
62606	Caloxin 3A1	WSSTSSVSAPLEFGGGGSAK	1912.1	1 mg
61345	<i>C. difficile</i> Toxin B (8-16)	QLEKMANVR	1088.3	1 mg
62505	α -C-Conotoxin PrXA	TYGIYDAKP-X-FSCAGLRGGCVLP-X-NLR-X-KFKE-NH ₂ (X = 4-trans-hydroxyproline, Disulfide bridge: 13-20)	3541.2	0.5 mg
61346	<i>C. difficile</i> Toxin B (192-207)	AQREENPELIIDDIVK	1882.1	1 mg
61347	<i>C. difficile</i> Toxin B (232-241)	ITQNSGNDVR	1103.2	1 mg
61348	<i>C. difficile</i> Toxin B (248-262)	NGESFNLYEQELVER	1826.9	1 mg
61349	<i>C. difficile</i> Toxin B (392-401)	DSYCSNLIVK	1335.4	1 mg
61350	<i>C. difficile</i> Toxin B (529-536)	AQFEEYKR	1070.2	1 mg
62393	Complement C3a Anaphylatoxin (71-77), porcine	KPLGLAR	754	1 mg
61857	Complement Anaphylatoxin C5a (37-53), human	RAARISLGPRCIKAFTE	1889.3	1 mg
62860	Crabrolin	FLPLILRKIVTAL-NH ₂	1496	1 mg
62856	Granuliberin-R	FGFLPIYRRPAS-NH ₂	1422.7	1 mg
62871	Histamine Releasing Peptide II	FLPLILGKLGKLL-NH ₂	1423.9	1 mg
62866	HP-1 Toxin	LFRLIAKTLGSLM	1462.9	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
62867	HP-2 Toxin	FRLANTLGKIL	1358.7	1 mg
62868	HP-3 Toxin	IFGLLAKTLGNLF	1406.7	1 mg
62849	Kassininin S	FIPVTLALHAIKEKLN-NH ₂	1976.5	1 mg
62855	Mast Cell Degranulating (MCD) Peptide HR-2	FLPLILGKLVKGLL-NH ₂	1523	1 mg
62058	Pertussis Toxin Substrate, Gai3C20, 6-FAM-labeled	6-FAM-VFDAVTDVIKNNLKECGLY	2613.0	1 mg
62859	Pipinin-3	FLPIIASVAAKVFASKIFCAISKKC (Disulfide bridge 18-24)	2583.3	1 mg
61600	pro-ε-TxIX/12 Conotoxin Precursor, ε-TxIX (-12 - -1)	LKRTIRTRLNIR	1539.9	1 mg
62314	SNAP-25 (49-58)	MLDEQGEQLER	1347.5	1 mg
62315	SNAP-25 (181-206)	IMEKADSNKTRIDEANQRATKMLGSG	2865.3	1 mg
61360	Staphylococcal α-Toxin P 1	KTGDLVTYDKENG	1439.6	1 mg
61518	Tetanus Toxin (TT) peptide	QYIKANSKFIGIFE	1657.9	1 mg
62410	Tetanus Toxin (830-844)	QYIKANSKFIGITEL	1725	1 mg
62496	δ-Toxin (1-26), <i>Staphylococcus aureus</i>	MAQDIISTIGDLVKWIIDTVNKFTKK	2978.6	1 mg
62497	δ-Toxin (5-20), <i>Staphylococcus aureus</i>	IISTIGDLVKWIIDTV	1786.2	1 mg
62312	Vesicle-Associated Membrane Protein, VAMP (60-94)	LSELDADRADALQAGASQFETSAAKLKRKYW- WKNLK	4039.6	1 mg
62316	Vesicle-Associated Membrane Protein, VAMP (62-71)	ELDDRADALQ	1145.2	1 mg
62313	Vesicle-Associated Membrane Protein, VAMP (75-94)	SQFETSAAKLKRKYWWKNLK	2513	1 mg

Tyrosinase Peptides

61475	TRP-1, Tyrosinase-Related Protein-1 Fragment	MSLQRQFL	1022.2	1 mg
61474	TRP-2, Tyrosinase-Related Protein-2 Coding Region Fragment	LLPGGRPYR	1028.2	1 mg
62679	Tumor Necrosis Factor-α Converting Enzyme, TACE (807-823), human	ASFKLQRQNRVDSKETE	2036.2	1 mg
61494	Tyrosinase (56-70)	QNILLSNAPLGPQFP	1608.9	1 mg
61491	Tyrosinase (146-156)	SSDYVIPIGTY	1214.3	1 mg
61493	Tyrosinase (192-200)	SEIWRDIDF	1180.3	1 mg
61492	Tyrosinase (206-214), T9206	AFLPWHRLF	1186.4	1 mg
61512	Tyrosinase (240-251)	DAEKCDICTDEY	1404.5	1 mg
62711	[Ser ²⁴⁴] Tyrosinase (240-251)	DAEKSDICTDEY	1388.4	1 mg
61500	Tyrosinase (243-251), core nonamer sequence	KCDICTDEY	1089.2	1 mg
61468	[Asn ³⁷⁰]-Tyrosinase (368-376)	YMNGTMSQV	1030.2	1 mg
61456	[Asp ³⁷⁰]-Tyrosinase (368-376)	YMDGTMSQV	1031.2	1 mg
61495	Tyrosinase (450-462)	SYLQSDPDSFQD	1516.5	1 mg

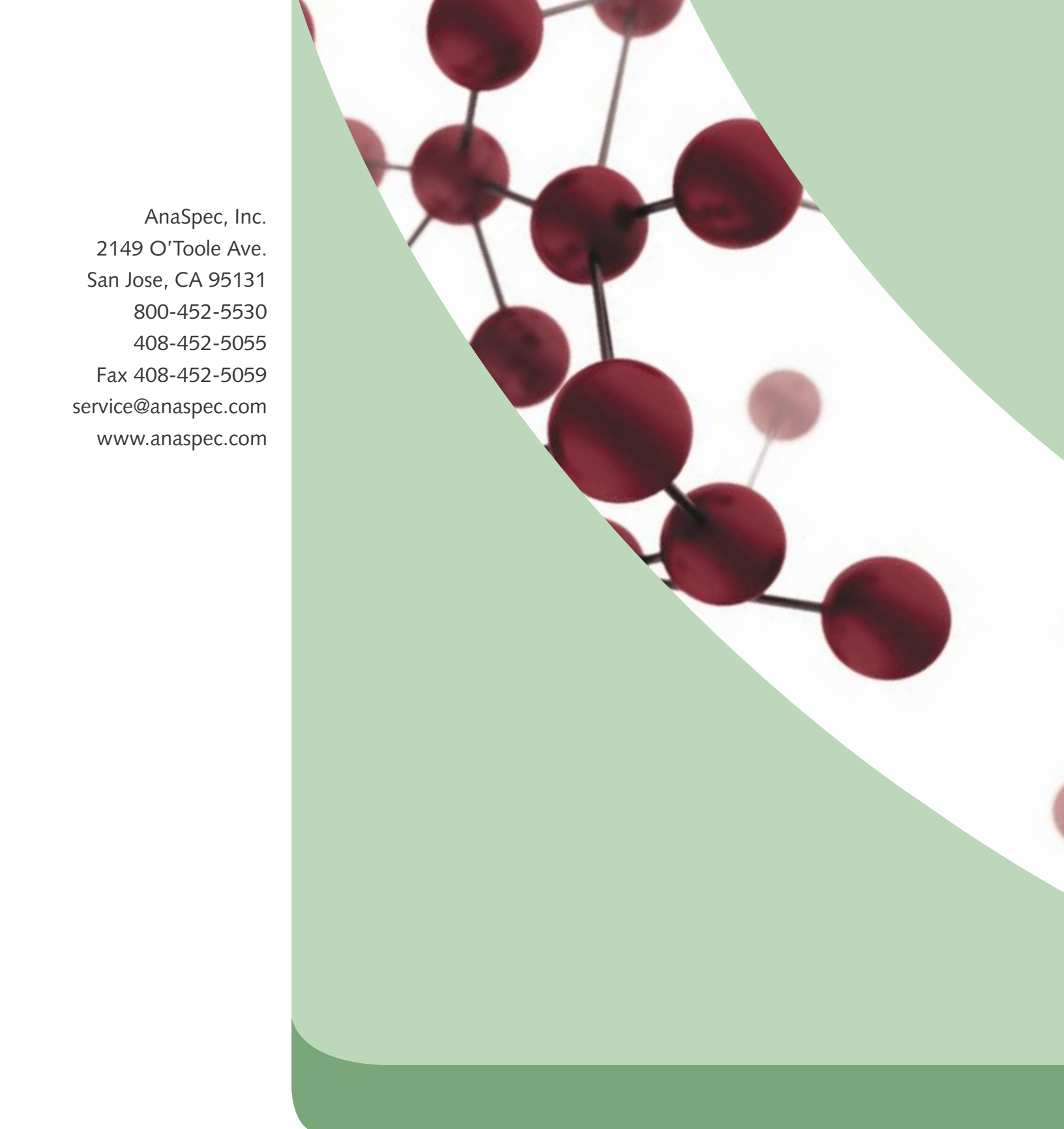
Viral Peptides

62631	Adenovirus Early DNA-Binding Protein, E2A (412-420)	LPKLTFFAL	999.3	1 mg
62782	ADP 715, HIV-1, MN Strain	RKRIHIGPGRAFYTTKN	2015.4	1 mg
62781	Cytomegalovirus Early Protein pp89 Derivative	YPHFMPTNL	1119.3	1 mg
62786	Cytomegalovirus Proteinase Substrate	RGVVNASSRLA	1129.3	1 mg
62032	Dengue 1 Fragment	VGIGNRDFV	976.1	1 mg
62033	Dengue 2 Fragment	IGISNRDFV	1020.2	1 mg
61590	HBcAg HBV, Hepatitis B Core Antigen (18-27)	FLPSDFFPVS	1155.3	1 mg
61594	HBV Core (107-115)	CLTFGRETV	1025.2	1 mg
61597	HBV Polymerase (575-583), HBVpol575	FLLSLGIHL	1012.3	1 mg
61598	HBV Core14, (HBA31)	STLPETTVVRR	1258.5	1 mg
61593	HBV env (183-191)	FLLTRILTI	1089.4	1 mg
61592	HBV env (335-343)	WLSLLVPFV	1073.4	1 mg

GO™ Peptides

CAT#	Product Name	Sequence	MW	Size
61591	HBV pol455	GLSRYVARL	1034.2	1 mg
61595	HBV pol502, HBV polymerase (502-510)	KLHLYSHPI	1107.3	1 mg
61596	HBV pol655, HBV polymerase (655-663)	ALMPYACI	994.3	1 mg
61864	HCV-1 e2 Protein (484-499)	PYCWHYPPKPCGIVPA	1828.2	1 mg
61865	HCV-1 e2 Protein (554-569)	WMNSTGFTKVCGAPPC	1699.0	1 mg
62220	Hepatitis G Virus (HGV), (2268-2276)	CDKCEARQE-NH ₂	1080.2	1 mg
62632	Hexon Protein Fragment	KYSPSNVKI	1035.2	1 mg
62694	HIV-1 env (584-594)	RYLRDQQLLGI	1374.6	1 mg
61708	HSV-gB2 (498-505)	SSIEFARL	922.1	1 mg
61589	Human Papillomavirus 18 (HPV18) E7 protein (86-94)	FQQLFLNTL	1123.3	1 mg
62607	KTL-501, Enterovirus Protein Fragment	PALTAVETGATNPL	1354.5	1 mg
61678	LMP1, Latent Membrane Protein 1 (156-164), IAL, EBV, Epstein-Barr Virus	IALYLQQNW	1148.3	1 mg
61676	LMP1, Latent Membrane Protein 1 (208-224), TDD, EBV, Epstein-Barr Virus	TDDSGHESDSNSNEGRH	1843.7	1 mg
61682	LMP2, Latent Membrane Protein 2 (131-139), PYL, EBV, Epstein-Barr Virus	PYLFWLAAI	1093.3	1 mg
61684	LMP2, Latent Membrane Protein 2 (236-244), RRR, EBV, Epstein-Barr Virus	RRRWRRRLTV	1298.6	1 mg
61681	LMP2, Latent Membrane Protein 2 (340-350), SSC, EBV, Epstein-Barr Virus	SSCSSCPLSKI	1111.3	1 mg
61683	LMP2, Latent Membrane Protein 2 (419-427), EBV, Epstein-Barr Virus	TYGPVFMCL	1030.3	1 mg
62408	Measles Virus Fusion Protein (288-302)	LSEIKGVIVHRLEGV	1649	1 mg
62409	Measles Virus Fusion Protein (421-435)	PVVEVNGVTIQVGSR	1553.8	1 mg
61706	Nuclear Localization Signal (NLS) SV40 Nonviral Gene Delivery Vector	PKKKRKVG	940.2	1 mg
62259	PB2 Polymerase (198-206) Mimotope, Influenza Virus	ISPLMVAYM	1024.3	1 mg
62021	Respiratory Syntical Virus (RSV), (127-135), M2	VYNTVISYI	1071.2	1 mg
62395	Respiratory Syntical Virus (RSV) M2 Protein 6	CHFSHNYFEWPPHAL	1885.1	1 mg
62391	Respiratory Syntical Virus (RSV) M2 Protein 10, bovine	EWPPHALLVR	1217.4	1 mg
62392	Respiratory Syntical Virus (RSV) M2 Protein 15	HNYFEWPPHALLVRQ	1907.2	1 mg
61658	Respiratory Syncytial Virus M Protein (RSV) (187-195)	NAITNAKII	957.1	1 mg
61618	SARS-CoV PLP2 Optimal Substrate (813-823)	FRLKGGAPIKGV	1139.0	1 mg
61619	SARS-CoV PLP2 Optimal Substrate (813-823), FRET	Abz-FRLKGGAPIKGV-EDDnp	1569.9	1 mg
61523	Tax8, HTLV-1, Human T Cell Lymphotropic Virus-1 (12-19)	LFGYPVYV	957.1	1 mg
61522	Tax9, HTLV-1, Human T Cell Lymphotropic Virus-1 (11-19)	LLFGYPVYV	1070.3	1 mg
62031	Viral Matrix Protein (128-135), M1 ¹²⁸	MGLIYNRM	997.3	1 mg
61709	Vesicular Stomatitis Virus (VSV-8) Nucleoprotein (52-59)	RGYGYQGL	913.0	1 mg
62026	Yellow Fever Virus Envelope Protein E1 Fragment	IGITDRDFI	1049.2	1 mg

Check our website for more newly released peptides
www.anaspec.com



AnaSpec, Inc.
2149 O'Toole Ave.
San Jose, CA 95131
800-452-5530
408-452-5055
Fax 408-452-5059
service@anaspec.com
www.anaspec.com



ISO 9001:2000 Company

© 2009 AnaSpec, Inc.