

3

3.5 Catalogue peptides

237 Amyloid Peptides

237 MOG Peptides

237 CEF Control Peptides

237 Histone H3 (1-21), Biotinylated

237 FRET protease substrates



Catalogue Peptides (Go™ Peptides)

**3500 catalogue peptides,
140 categories, ready to be shipped !**

Eurogentec and AnaSpec are pleased to introduce **Go™ Peptides**, the next evolution in catalogue peptides. Our catalogue peptides represent over 17 years of innovative peptide synthesis expertise. This represents one of the world's largest and fastest growing inventories, with new catalogue peptides added monthly. We are able to provide an enormous pool of research ready peptides, in inventory and ready for immediate shipment.

- **The most complete collection of β -Amyloid peptides, kinase/phosphatase substrates, phospho-peptides, FRET peptides and fluorescent imaging peptides.**
- **> 95 % pure, TFA as counter ion.**
- **Guaranteed stability for 1 year from date of purchase if stored at -20 °C as lyophilized powder.**

Our valued customers have chosen Eurogentec/AnaSpec as provider of one of the following products:

Product example	Size	Reference
Beta-Amyloid 1-42	1 mg	AS-20276
MOG (35-55), mouse rat	5 mg	AS-60130-5
Peptide Substrate Renin 520	1 mg	AS-61872
Beta-Amyloid 1-40	1 mg	AS-24236
CEF Control Peptide Pool	16 mg	AS-61036-05
HCV NS3/4A Protease Substrate	1 mg	AS-60798
HCV Protease Substrate (RET-S1)	1 mg	AS-22991
OVA (323-339)	5 mg	AS-27025
Histone H3 (1-21), Biotinylated	1 mg	AS-61702
Beta-Amyloid 1-42	0.5 mg	AS-24224

Amyloid Peptides

β -amyloids are the core components of neuritic plaques seen in Alzheimer's disease brains.

- β -amyloid 1-40, 1-42, 1-28 and dozens of variations
- Labeled peptides (FAM, HiLyte Fluor™, Biotin,...)
- Mutated peptides
- Scrambled and inverted sequences
- APP fragments
- Also mouse and rat β -amyloid
- Other Alzheimer-related peptides

MOG Peptides

- Myelin Oligodendrocyte Glycoprotein (MOG) peptides can induce immune system to damage myelin sheath around axons of spinal cord. MOG peptides are also involved in Experimental Autoimmune Encephalomyelitis (EAE), an animal model for human Multiple Sclerosis.
- MOG 35-55 induced 100 % EAE in mice

CEF Control Peptides

CEF Control Peptide Pool is a group of 32 viral peptides that stimulate gamma-interferon production from CD8+ T cells

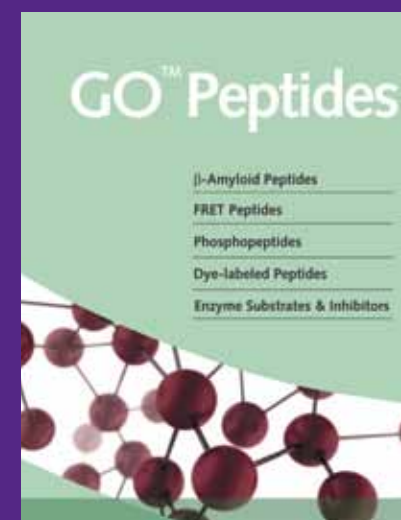
Histone H3 (1-21), Biotinylated

This sequence corresponds to the first 21 amino acids of the NH₂ terminal of histone H3 followed by a GG linker and a biotinylated lysine. This peptide was used to investigate the characteristics and mechanisms of ethanol-induced histone H3 acetylation in rat hepatocytes (Park, P. *et al.* Am. J. Physiol. Gastrointest. Liver Physiol. 289, G1124 (2005)).

FRET protease substrates

In recent years, FRET-based assays have been used extensively in the different proteases, which made the continuous assay of protease activity and HTS of protease inhibitors faster and easier than other analysis methods.

Full list on
Eurogentec website
www.eurogentec.com



See also the related
Neuroscience catalogue
antibodies, page 215