

Peptide Manufacturing

For Discovery to Clinical Phase



Custom Peptide Synthesis

Leverage our vast expertise in liquid and solid phase peptide synthesis, gained from over **30 years** of experience serving **industry, academia, and governmental** agencies. We produce simple **unmodified** to highly **modified** and **hydrophobic peptides** from our manufacturing facility based in the U.S.



LEARN MORE ON

**Custom
Peptides**



Quality

Expert manufacturing by highly experienced chemists under our QMS system.



Fast Delivery

Our processes and facility, are streamlined for efficiency to ensure on-time delivery.



Purity

Up to $\geq 95\%$ purity for Research grade and up to $\geq 98\%$ purity for GMP grade custom peptides.



Specifications

- 1mg-to-hundreds of gram quantities
- Up to 60 amino acids in length
- Crude to $\geq 98\%$
- Purity tested by HPLC analysis
- Identity tested by Mass Spec analysis



Modifications

We can incorporate single or multiple modifications on your peptide, at the sequence N-terminus, C-terminus, and/or internally.

- UAADs & Analogs
- Dye-Labeling and Conjugations
- Cyclic & MAP peptides



Check out our collection of off-the-shelf Catalog Peptides

We maintain one of the world's largest inventories of catalog peptides, with new peptides added regularly. For your convenience, our collection of 1000+ peptides are organized across 11 research areas and subcategories.

Specifications

Bulk quantities available upon request

Purity

>95% by HPLC unless otherwise specified

Identity

Tested by Mass Spec analysis

Quantity

Provided as gross weight as per industry standard, unless otherwise specified



LEARN MORE ON

**Catalog
Peptides**

Research areas



Adhesion &
Extracellular Matrix



Cancer & Apoptosis



Cell permeable &
Penetrating Peptides



Cell Signaling



Cardiovascular



Diabetes



Epigenetics



Immunology
& Infectious Disease



Neuroscience



Peptides Hormones



Peptide & Protein
Analysis

Kaneka Eurogentec SA

5 Rue Bois Saint-Jean, 4102 Seraing, Belgium
T +32 (0) 4 372 74 00

Learn more on www.eurogentec.com

