

## Efficient Project Management

With our robust project management, we ensure **timely and efficient delivery** of your custom peptides. Our **dedicated project managers** work closely with you to understand your requirements, provide regular updates, and **ensure that all milestones are met.**

## The Benefits of Choosing Eurogentec

### One-Stop Provider

We offer a **5-in-1 solution**, including **GMP manufacturing**, Analytical Method **Validation**, **Stability** Programs, **CMC** assistance for IND filing, and **GLP study** packages.

### Small to Large Scale

Our end-to-end approach, provides a **single point of contact** for process development, manufacturing validation, and quality control—streamlining your workflow.

### Fast Lead-Times

**Quick delivery** of peptide API for clinical use, emergency IND use, API precursors and excipients, and other critical raw materials.

### Quality Management

Our comprehensive **QMS policy** ensures **top-quality products** from raw materials to manufacturing, QA, QC, shipping, and beyond.

### Transparent Communication

We provide **fast, responsive**, and **transparent communication**, with access to key team members. Client audits are always welcome.

### Made in the USA

Our state-of-the-art facility in **Silicon Valley** houses **all departments**, including manufacturing, QA, QC testing, EH&S, and materials management.



 Eurogentec

 eurogentec.com



# GMP Peptides

**One-Stop Service Provider**

We specialize in **GMP peptide manufacturing**, leveraging decades of expertise, a rigorous quality management system, and state-of-the-art facilities to deliver **complex-modified peptides** and **peptidomimetics** tailored to your unique requirements.



SCAN THIS CODE AND

**Discover more about our services**

THE-EN-FLY-GMP-PEPTIDES-A4-OCT2024-V1



**30 years**  
of experience



**mg to kg**  
net quantity



**Fast**  
lead-times



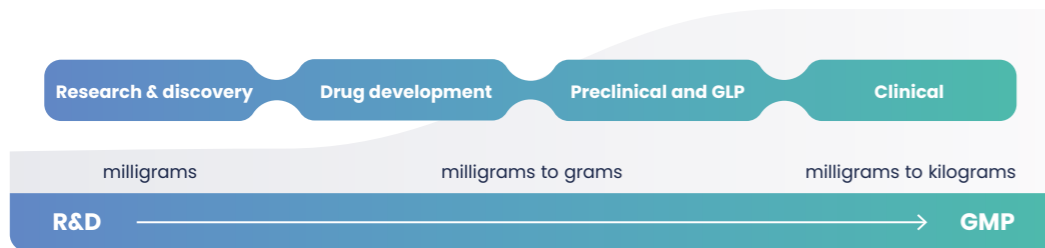
Made  
in the **USA**

Get the right quantity and quality, at the right time.



to  
**98%**  
purity by HPLC

Each peptide we produce undergoes rigorous testing to ensure it meets your pre-established acceptance criteria.



## Comprehensive GMP Peptide Manufacturing Services

Our fully customizable GMP peptide service is designed to meet the diverse requirements of our clients, delivering high-quality peptides at the scale that suits your specific needs.

We support each stage of your GMP peptide drug development, from early R&D to clinical, ensuring consistent quality and compliance throughout your project's lifecycle.

## From milligrams to kilograms

We produce peptides in quantities ranging from milligrams to kilograms, in our facilities equipped to handle both small-scale production and large-scale manufacturing.

We offer both Gross and Net quantities of peptides, packaged in the amount per vial as requested.

Our flexible production capabilities ensure you receive the right peptide quantity and quality according to your specifications.

### IN-VIVO & GLP STUDIES

We give you access to *in vivo* and *in vitro* preclinical studies prior to progress to clinical trials.

### PEPTIDE MODIFICATION EXPERTS

Our expert team brings decades of experience in peptide synthesis and purification of highly modified peptides including:

#### Labeled

- Fluorescent Dyes
- FRET
- TR-FRET
- Heavy isotope

#### Structural

- Lactam ring cyclic
- Disulfide-bridge
- Stapled
- Thioether-bridge
- Thiolactone cyclization

#### Conjugated

- Drug-Peptide (PDC)
- Chelation
- Peptide-Oligo Conjugates
- Carrier Proteins

#### Specialized

- Unusual Amino Acids
- Lipopeptides
- Phosphorylated
- Glycosylated
- Peptidomimetics



## State-of-the-Art Manufacturing Facilities

Our cutting-edge 44,000-square-foot facility in California's Silicon Valley is designed to support high-quality GMP peptide production with state-of-the-art equipment and ISO 7 cleanrooms.

We adhere to strict GMP guidelines, including secure airlock entry, rigorous environmental monitoring, and thorough segregation of processing areas.

## Solid Quality Assurance

Our comprehensive Quality Management System (QMS) adheres to 21 CFR 210 and 211 as well as ICH Q7 guidelines, applicable for products intended use.

Key elements of our QMS include:

- Regulatory Compliance:** We follow GMP standards and undergo regular audits to ensure quality.
- Raw Material Control:** We verify the quality of starting materials and qualify suppliers.
- In-Process Control:** Robust controls during production ensure our products meet specifications.
- Quality Control:** We use qualified analytical methods to verify peptide quality.
- Stability Testing:** Stability samples are maintained and monitored in validated storage chambers, upon request.
- Documented Process:** Our GMP manufacturing process is meticulously documented and archived for five years, with redacted records available for transparency.