



Therapeutic service overview

Accelerate your drug development process while maintaining the highest GMP standards.

GMP accredited company since 1994

At Eurogentec, we specialize in delivering **multimodality GMP manufacturing** solutions to the **pharmaceutical** and **biotechnology** industries.

Our CDMO expertise is unique, spanning five distinct therapeutic modalities and enabling comprehensive support from early discovery through to commercial supply.

Our advanced facilities and expert team offer a wide range of services for the development and manufacturing of **starting materials** and **Active Pharmaceutical Ingredients (APIs)**, from the discovery phase to preclinical research, clinical trials, and commercialization.



Chemical synthesis

GMP oligonucleotides

We manufacture **small and large scales** oligonucleotides from early-stage development to clinical phases.

We offer a **broad range of chemistries** and **modifications**, including siRNA, antisense, aptamer, and CpG oligonucleotides.

GMP peptides

We offer continuum **from research to clinical trials**.

Our GMP grade peptides can be used as **API drug** substance in Phase 1 Clinical Trials, Emergency IND (EIND) use, API precursors and **starting materials**.

Our **multi-zone facility** houses all depts from manufacturing, QA, QC testing, EH&S, materials management and beyond.

Hundreds of grams

Production scale



Clinical trials



Up to kg production scale



Phase 1 & Clinical trials





Enzymatic synthesis

GMP mRNA

We have combined our expertise in GMP biopharmaceutical manufacturing with our know-how in large scale oligonucleotide synthesis to produce **long** and **stable RNA** molecules.



For human **clinical trials & commercial supply**

50g

Up to **50g mRNA** production scale

Bio manufacturing

GMP plasmid DNA

We manufacture plasmid DNA in bacteria for **vet and human** use. We can deliver **from gram to kg** quantities of plasmid for use as DNA vaccines, non-viral gene therapy or starting material for viral vector and mRNA production.



Up to **Kilo scale**

3

Up to **Phase 3 & Commercial use**



GMP recombinant proteins & conjugates

We produce recombinant proteins for human use in **bacteria and yeasts** with a **highly scalable** approach. We have a strong expertise in **coupling** peptides, proteins, polysaccharides and PEG to proteins.



Scalable process

3

Up to **Phase 3 & Commercial use**



A flexible and transparent process



Our end-to-end solutions ensure consistent quality, safety, and compliance.



Expertise

With 35 years of GMP manufacturing experience, our team of experts offers invaluable guidance on process optimization, and problem-solving to meet your needs.



Quality

We ensure GMP compliance through rigorous quality assurance and a deep understanding of international regulatory guidelines, offering reliable manufacturing solutions.



Comprehensive services

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



Flexibility

Our multimodal manufacturing approach allows scalable production, adapting seamlessly to different stages of drug development.



State-of-the-art facilities

Our facilities are engineered to meet the rigorous requirements of GMP, featuring state-of-the-art technology to ensure precision and consistency.



Transparent communication

Our effective project managers keep you informed at every step, ensuring smooth collaboration and successful project execution from start to finish.



Scan this code and
Discuss your project with us!

Our production sites

America
AnaSpec, Inc. - Fremont CA (USA)

Europe
Kaneka Eurogentec Headquarters
Liege (Belgium)