

Therapeutic Oligonucleotides

Research & GMP grade Manufacturing

We support scientists worldwide in their quest to develop and produce **therapeutic oligonucleotides** either if they are in **early discovery**, preclinical studies or **clinical trials**.



40 years
of experience



10 M oligos
manufactured



250+
Modifications



End-to-end
assistance

Our expertise in therapeutic oligonucleotide

We offer comprehensive and customized solutions for the **development, manufacturing, and testing** of oligonucleotides. Our services cover every stage from **research and discovery** to oligo **screening libraries** and the production of **larger quantities** needed for **preclinical** and **early clinical** studies.

With us, get your oligo in the quality and quantity adapted to your needs.

Diverse chemistries & modifications

We specialize in designing and synthesizing a wide range of oligonucleotides, including **DNA, RNA, ASOs, siRNA, miRNA mimics, aptamers** and **CpG oligos**.

Our expertise covers a broad range of chemistries (**2'O-Me-RNA, 2'F-RNA, etc.**) and modifications (**Fluorescent dyes, GalNAc, PEG, lipids** etc.) to enhance stability, binding affinity, specificity, and delivery of your therapeutic oligonucleotides.

Peptide Oligonucleotide Conjugates

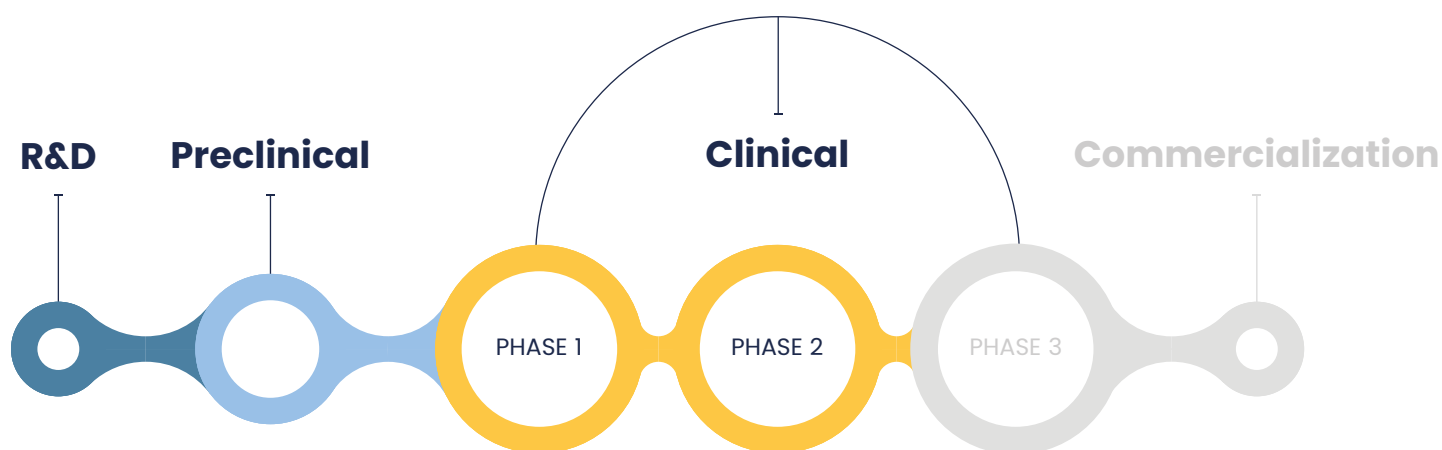
Leverage our years of expertise in peptides and oligonucleotides for custom production of innovative peptide-oligonucleotide conjugates.

From micrograms to hundreds of grams

We produce oligonucleotides in the quantities and quality **that best meet your application needs**. Whether for research, preclinical studies, or clinical phases, you can be confident that you will receive exactly what your application deserves.

IN-VIVO & GLP STUDIES

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





High quality manufacturing process

We ensure the highest quality and efficacy for every oligonucleotide we deliver. Our **documented** and tested manufacturing process guarantees **reliability** at each step.

For clinical oligonucleotides, we adhere to a validation approach that **complies with ICH** guidelines, facilitating compliance with FDA, EMA, and other regulatory requirements.

State-of-the-art facilities & equipment

We have **multiple rooms**, from unclassified laboratories to **ISO7** and **ISO8** cleanrooms equipped with the latest technologies, to produce, purify and control the quality of the oligonucleotides we manufacture.

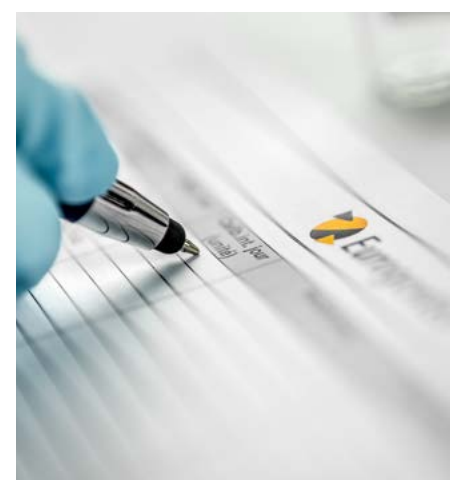
QC analytical methods

QC analytical methods are applied **on raw materials, during the production and on final products.**

They include but are not limited to :

- Appearance, pH, Osmolarity,
- Purity and quantification of the main impurities including profiling by LC-MS,
- Quantification of Bioburden and Endotoxins,
- Identity by Mass Spectrometry (ESI), Sequencing, Extinction coefficient determination,
- Residual solvents, Residual heavy metals, Water content, Sodium content,
- Phosphorothioate - Phosphate content

Available
STABILITY TESTING



Comprehensive Project Oversight and Coordination

Our experienced project managers maintain clear communication, act proactively to **anticipate** potential challenges and implement **strategic solutions** for smooth and efficient progress from initial conception to final delivery.

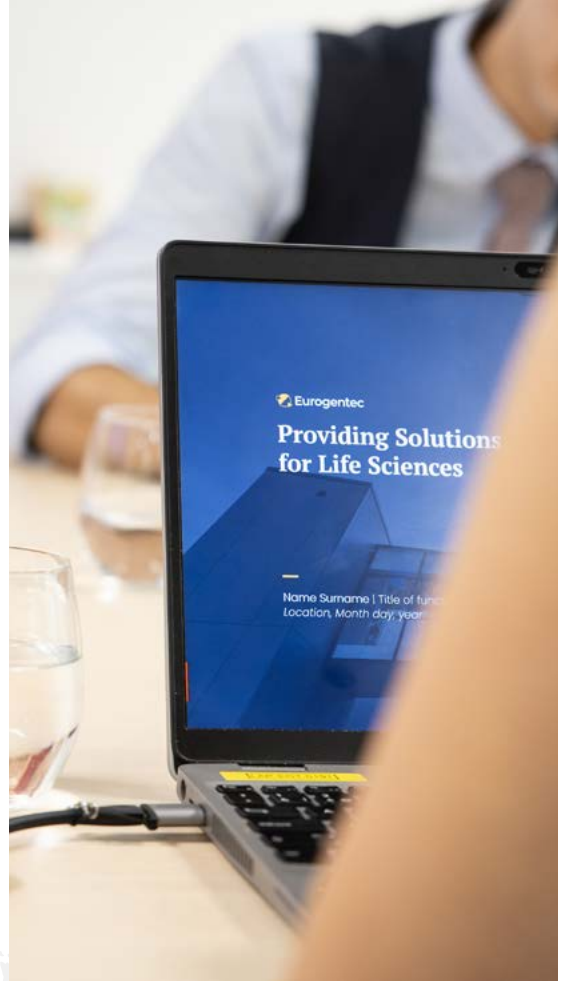
Support in the writing and review of CMC

Our dedicated team of **scientists** and **regulatory experts** ensures that every aspect of your **CMC information submission** meets regulatory standards.

We provide comprehensive assistance for your **IND** (Investigational New Drug) file preparation and deliver **documentation** tailored to your specific needs.



← SCAN THIS CODE AND
DISCOVER MORE ABOUT OUR SERVICES



California

**Kaneka Eurogentec
North America**

Fremont

Belgium

**Kaneka Eurogentec
Headquarters**

Liege