

**Providing Trusted
and Innovative
Solutions to
the Life Science
Communities**

Eurogentec

We are a **leading supplier** of trusted and innovative reagents, kits, specialty products and custom research/development services. Our 3 Business Units: Research Products, Diagnostic Services and Therapeutic Manufacturing; provide high quality products to support scientists involved in the **Life Science** research, **biotechnology**, **diagnostic** and **pharmaceutical** markets.

Eurogentec is headquartered in the Liege Science Park (Belgium) and has subsidiaries in France, Germany, The Netherlands, the United Kingdom, Switzerland and USA (AnaSpec Inc.) and production sites in Belgium, United Kingdom and USA.

Eurogentec has been a part of Kaneka Corporation since 2010.



kaneka

Kaneka Corporation is an innovation-oriented chemical company. Business activities span a broad spectrum of markets ranging from plastics, EPS resins, chemicals and foodstuffs to pharmaceutical intermediates, medical devices, electrical and electronic materials and synthetic fibers.

Kaneka Corporation is headquartered in Osaka-Japan, with subsidiaries in the United States, Belgium, Singapore, Malaysia, China, Australia, Vietnam, India, South Korea and Taiwan.

Kaneka employs more than 11000 people worldwide.

Research

Provides R&D laboratories with trusted reagents and innovative services for genomics and proteomics.

[Oligonucleotides

[Antibodies

[qPCR and PCR kits and reagents

[Gene analysis

[Protein iDentification

[AnaSpec™ Peptides



Products

> Innovative catalog reagents

- Room temperature stable lyophilized qPCR and Master Mixes
- DNA extraction & purification kits
- Catalog peptides
- Highly sensitive enzymatic assay kits

> Flexible and customized solutions

- From PCR primers to highly modified oligonucleotides (including MGB probes, NGS oligos ...)
- Custom monoclonals, polyclonals & antibody fragments
 - Small to large antibody project management
 - Antibodies against post-translational modifications
- Specialized peptide synthesis
- Lab Developed Test Solutions
- Tailored amplification reagents

> Unique products & services

- Efficient PCR DNA polymerase – Takyon™
- 28 days polyclonal immunisation programmes
- Proprietary dyes & quenchers
- Quantified peptides
- Optimized qPCR reagents for viral RNA detection



FS 638601



info@eurogentec.com



Diagnostic

Guarantees high quality, IVD grade components dedicated for diagnostic applications.

- [IVD oligonucleotides
- [Critical Raw Material peptides
- [Lab developed test solutions
- [Kit Manufacturing for nucleic acid-based assays



Services

- > Broad expertise
 - IVD oligonucleotides, peptides
 - Complex oligo chemistry expertise
 - IP friendly dyes & quenchers
 - State-of-the art quality management system
- > Manufacturing capabilities
 - From small to large scale oligonucleotide production
 - Extended oligo grades (Track™, pre-Diagnostic, Diagnostic)
 - Classified cleanrooms
 - Highly qualified production teams
- > Custom solutions
 - Collaborative project development
 - Research Use Only to GMP components including ASRs
 - Assay dispensing services
 - Custom fill & finish

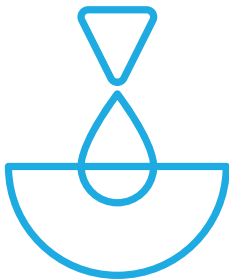


MD 638600 For production & sales of IVD oligonucleotides



FS 638601

 ivd@eurogentec.com



From research to commercialization

Therapeutic

Develops and produces material for clinical trials and commercial products since 1994.

- [Microbial expert
- [Process development
- [GMP plasmid DNA
- [GMP proteins
- [GMP protein conjugates
- [GMP therapeutic oligos
- [GMP mRNA



Manufacturing

- > Broad Product Experience
 - Chemical Synthesis of GMP therapeutic grade oligonucleotides (clinical phases I & II)
 - GMP grade RNA (clinical phases I & II)
 - Sterile injectable products from microbes including :
 - Plasmid DNA
 - Recombinant proteins (antibody fragments, enzymes, cytokines, subunit vaccines, others)
 - Protein conjugated to PEG, peptide, polysaccharide or another protein
 - BSL 2 protein extracts
- > A Comprehensive GMP offering
 - Fermentation, purification & quality control development
 - Tox, clinical & commercial manufacturing
 - ICH stability studies
 - Regulatory affairs
- > Unique Manufacturing Platforms
 - High cell density fed-batch fermentation processes
 - Continuous & batch purification processes
 - Disposables & single-use technologies
 - Large scale manufacturing of Biologics
 - Plasmid DNA manufacturing platform to kg scale



For our GMP BioManufacturing facilities

✉ info.biologics@eurogentec.com



Corporate Social Responsibility



Eurogentec is committed to adhere to and fully adopts Kaneka's code of ethics, fulfilling our corporate social responsibility through the materialisation of our corporate philosophy with the earnest and forward-looking efforts of each employee.

The company continuously strives to improve its processes and practices based on Kaneka's "Zero tolerance" philosophy.

The full version of our CSR policy is available at: www.eurogentec.com

Confidentiality



We believe that being aware of our customers' needs and expectations is essential to build a true partnership. A dedicated Project Manager or Technical Specialist is always available when assistance or advice is needed. Every project is discussed in detail to clearly define timelines and expectations.

We ensure that your project and your data will be handled with complete confidentiality and never shared with third parties.

All data/results related to your project remain your property and a non-disclosure agreement can also be executed if necessary.



Quality

ISO certifications

The Quality Management Systems (QMS) of our Genomics and Proteomics activities are certified according to ISO 9001 standard.

Eurogentec is certified for the "Development, Production and Sales of products and services for Research and Development in Life Science".

AnaSpec, Inc. is certified for the manufacture/ re-packaging and supply of peptides, antibodies, assay kits, synthesis resins, amino acids, dyes, reagents for Life Sciences research and distribution of qPCR reagents.

In addition Eurogentec holds ISO 13485 certification for the contract manufacturing of synthetic oligonucleotides in support of *in vitro* diagnostic and related applications.

Directives for pharmaceuticals

Eurogentec is accredited by the Belgian Federal Agency for Medicines and Health Products for the manufacture of clinical trial and commercial material.

Eurogentec manufactures biopharmaceuticals according to the European Commission's Directive 2003/94/EC, EudraLex Volume 4 regulations as well as the US FDA's 21 CFR Part 210 and 211 GMP manufacturing requirements.

Environment

As a leading provider to the Life Science Community, Eurogentec is actively involved in environmental protection initiatives. We proactively develop new programs to reduce our environmental footprint without compromising the quality and efficiency of our products and services.

In an effort to protect the environment, Eurogentec has taken the following actions:

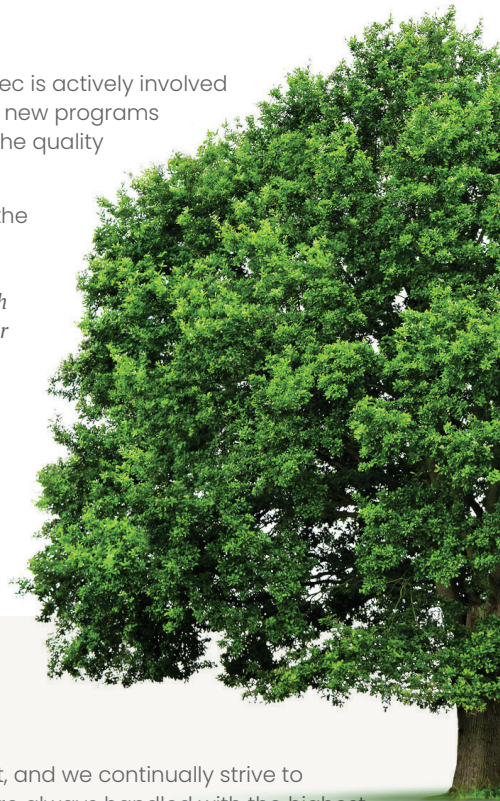
> *New laboratories built following latest energy efficiency standards.*

> *Production of RT stable products that require no cold chain transportation and storage.*

> *Exclusive use of carbon-neutral gas, certified from "Green Origin" (N₂, O₂ and Ar).*

> *Paper product inserts replaced with electronic versions available online or via e-mail.*

> *Use of natural and biodegradable packing peanuts made from cornstarch.*



Animal Welfare

Eurogentec is committed to treating every animal with respect, and we continually strive to enhance animal comfort and reduce animal stress. Animals are always handled with the highest care and respect in compliance with the latest rules/guidelines.

Feedback from customer visits and audits confirm that our animal facility is one of the cleanest and safest animal housings in the industry.

We respect the 3Rs philosophy: reduce the number of animals used; refine space to improve animal comfort; replace animal use by other techniques whenever possible. With this aim, we continue to recommend and suggest alternatives to *in vivo* antibody development and production (e.g. monoclonal antibodies are exclusively produced *in vitro*).



Kaneka Eurogentec S.A. (Headquarters)
LIEGE Science Park, rue Bois Saint-Jean, 5
4102 Seraing - BELGIUM
☎ +32 4 372 74 00
☎ +32 4 372 75 00
✉ info@eurogentec.com
🌐 www.eurogentec.com



Kaneka Eurogentec S.A. - Biologics
Kaneka Eurogentec S.A. - rue Bois Saint-Jean, 14
4102 Seraing - BELGIUM
☎ +32 4 372 74 00
☎ +32 4 372 75 00
✉ info.biologics@eurogentec.com
🌐 www.eurogentec.com



AnaSpec Inc.
34801 Campus Drive - Fremont
CA 94555 - USA
☎ +1-510-791-9560
☎ +1-510-791-9572
☎ **+1-800-452-5530**
✉ service@anaspec.com
🌐 www.anaspec.com



www.eurogentec.com



Follow us on

- www.linkedin.com/company/eurogentec
- www.twitter.com/eurogentec
- www.facebook.com/eurogentec