

Dispensing Services

Real-Time qPCR & qRT-PCR assays



PCR & RT-PCR assays



Peptide & protein based assays



Primers & probes blends



High-Throughput screening assays



Simply add your template

- Ready-to-Use
- Cost effective
- Easy-to-Use
- Better productivity
- Faster results

- ✓ Virtually no reaction set-up time
- ✓ Minimal labour required, no more «hidden» costs
- ✓ Minimal expertise required
- ✓ Increase number of data points
- ✓ Publish earlier

Turn homebrew into diagnostic

- Secure
- High consistent quality
- Reproducible
- In line with ISO 15189

- ✓ No contamination risk!
- ✓ Fully controlled equipment, processes and reagents
- ✓ Absolute accuracy & reproducibility guaranteed by our ISO 9001/ ISO 13485 manufacturing

Focus on your data

- Precise
- Flexible
- Cost effective

- ✓ Work confidently with small volumes
- ✓ No concessions with project design, e.g. complex plate set-up, format and support
- ✓ No waste of costly reagents, no investment needed

EUROPEAN OFFICES

BELGIUM

EUROGENTEC HEADQUARTERS
TEL.: +32 4 372 74 00
FAX: +32 4 372 75 00
info@eurogentec.com

EUROGENTEC BIOLOGICS

TEL.: +32 4 372 74 00
FAX: +32 4 365 16 04
info.biologics@eurogentec.com

UNITED KINGDOM

EUROGENTEC LTD.
TEL.: + 44 1794 511 411
FAX: + 44 1794 522 417
info.uk@eurogentec.com

FRANCE

EUROGENTEC FRANCE SASU
TEL.: +33 2 41 73 33 73
FAX: +33 2 41 73 10 26
info@eurogentec.com

GERMANY

EUROGENTEC DEUTSCHLAND GMBH
TEL.: +49 221 258 94 55
FAX: +49 221 258 94 54
info@eurogentec.com

THE NETHERLANDS

EUROGENTEC NEDERLAND B.V.
TEL.: +31 43 352 06 98
FAX: +31 43 354 19 65
info@eurogentec.com

SWITZERLAND

EUROGENTEC S.A.,
SUCCURSALE DE GENÈVE
FRANÇAIS:
TEL.: +32 4 372 74 00
FAX: +32 4 372 75 00
DEUTSCH:
TEL.: +49 221 258 94 55
FAX: +49 221 258 94 54
info@eurogentec.com

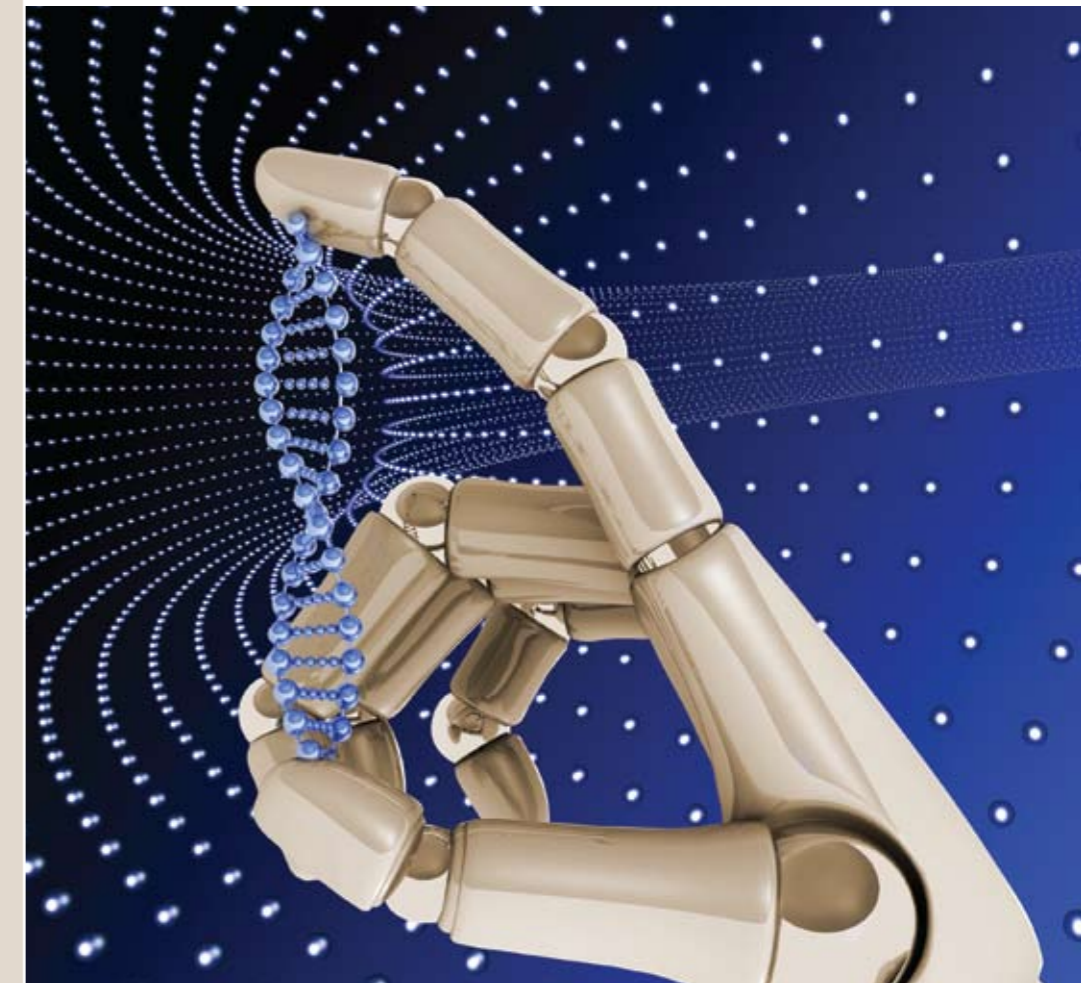
CENTRAL EUROPE

REPRESENTED BY
EUROGENTEC HEADQUARTERS
TEL.: +32 4 372 74 00
FAX: +32 4 372 75 00
info@eurogentec.com

www.eurogentec.com

To start discussing your project and get a quotation adapted to your needs, please contact us at dispensing@eurogentec.com

HIGH-THROUGHPUT DISPENSING SERVICES



MORE DATA, LESS WORK
MORE PRECISION, NO RISK

High-Throughput Dispensing Services

In its constant search for simple, affordable and innovative solutions for the Life Science community, Eurogentec is introducing its dispensing services that bring all advantages of automation within everyone's reach.

Ease your workflow

Automate your routine assays to perfection, with full design and format flexibility. Combine all required reagents, in any type of formulation, support and packaging, to produce your Ready & Easy-to-Use solution, minimizing user input, time and skill level, as well as hidden costs!

Our dispensing services are fully scalable and can be adapted to any requirement, for all types of use ■

Set new quality standards

Reliability, accuracy and risk management are essential parameters in all PCR and qPCR-based applications, especially in this new era of high-throughput projects.

The combination of our state-of-the-art liquid handling platform, with our complete range of high quality reagents gives you the possibility to automate your routine assays with more precision and safety.

Eurogentec can provide you with all needed reagents, from research to IVD/GMP grade, including oligonucleotides, probes, enzymes, MasterMixes and plastics, for single projects up to large OEM productions.

Enforcement of our ISO 9001 and ISO 13485 certifications is a key guarantee of the quality consistency of our products and processes. All production is performed with a perfect control of all contamination risks, with uniform dosage and ultra low wastage ■

Bypass human limits

Make no concession with your project design. Automate your project, whether small or large, beyond human handling limits with utmost reliability, accuracy and no risk of contamination, even at the lowest volumes. Thanks to the flexibility of our integrated production facility, combine all types of reagents and supports, from single-tubes to strips and 96 or 384-well plates, in every possible format. Complex plate structures can easily be automated with uniform dosage and no wasting of costly reagents.

Contrary to current High-Throughput nanoliter-based technologies, our services offer you the possibility to benefit from all advantages of automation without the need of costly investments, assay parameter changes and controversial pre-amplification steps ■

From first contact to mature projects, we consider each request with utmost priority. A dedicated project manager and technical specialists ensure perfect co-ordination between our customer and the production platform to efficiently convert their project into ready and easy-to-use solutions fulfilling all their specific requirements.

↘ QC set-up

A QC Release Method is developed in collaboration with the customer to ensure full compliance with their quality and functionality requirements. We provide a large selection of standard and customized QC methods, in addition to retention samples and complete product documentation traceability from Research to GMP grade.

↘ NDA

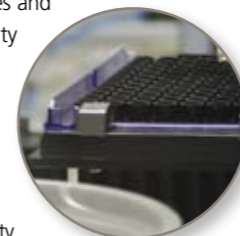
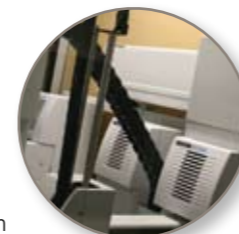
All projects are handled with complete confidentiality. A Non-Disclosure Agreement is executed during the initial stages of communication.

↘ Experienced technical support

From first contact to running projects, we provide our customers with a dedicated, highly experienced technical support. Eurogentec has over 20 years of experience in key-related fields, from nucleic acid design and synthesis, PCR and Real-Time PCR reagents manufacturing, to peptides and custom antibody production.

↘ Quotation

When quality standards and a supply agreement including all manufacturing specifications are defined, a quotation is provided with specific reference, product details, guaranteed price and lead-time information. The quotation includes the production of a first sample batch as well as optional services and QCs.



Project development

Early-stage projects are welcome, from small-scale production to large OEM manufacturing. A feasibility study is performed in collaboration with the customers, to define the reagents and manufacturing specifications that will guarantee the success of their project ■

Project transfer

Mature projects can easily be transferred and implemented into production, thanks to the outstanding flexibility of our manufacturing platform, in terms of supports, reagents, formats and processes ■

Reagents validation

Before the sampling phase, we assist our customers in validating their assays with our reagents, with the support of experienced technical specialists. Eurogentec molecular biology reagents include a wide range of products, many of those can be customized to fulfill all specific requirements. > Best-of-class Real-Time qPCR reagents validated on all cyclers, for all applications.

- > PCR enzymes and mixes adapted to all template types and application requirements, ranging from research to GMP-grade.
- > Synthetic peptides, recombinant proteins, catalogue and custom antibodies.
- > Primer & probes, including hundreds of chemical modifications, from research to IVD/GMP grade, covering early phases of development to clinical validation and routine diagnostic use ■

Reagents synthesis

All reagents undergo rigorous Quality Control testing to guarantee excellent quality and consistency of the finished goods. We offer a large selection of standard and customized QC methods, any of which can be incorporated into the customized QC Release Method:

- > Final mass determination
- > MALDI-TOF or ESI Mass Spectrometry
- > HPLC RP & ion exchange
- > Enzymatic tests e.g. DNA-enzyme conjugates
- > Probe/dye activity: fluorescence, signal to noise ratio, melting profile for MB
- > Functional QCs e.g. amplification assays ■

Sample batch production

When reagents are validated and a quotation is accepted, a sample batch is produced according to defined QA/QC and manufacturing specifications. A full QC, as defined with the client, is performed to ensure quality and functionality of the dispensed reagents. QC-approved samples along with all QC data, including individual reagents QCs, are sent to the customer ■

Sample batch validation

The customer receives the released sample batch, including all QC results, for internal validation. A dedicated project manager and/or technical specialist is available when assistance or advice is needed ■

Master batch production & release

When sample batch is validated, a large batch is produced upon customer request, according to the quotation details and manufacturing specifications. A full QC, as defined in the QC Release Method, is performed to ensure full compliance with QA and product specifications. Our products are provided with a certificate of analysis, comprising a detailed document with all QC results for the research reagents or a detailed document released after review of the batch record and signed by a QC authorized person for the IVD reagents ■