



SIMPLIFY ROUTINE qPCR

We prepare.

You test.

Custom qPCR assay preparation service

Eurogentec is your one-stop shop for ready-to-use qPCR premix solutions. As a trusted reagent supplier, we also offer services to simplify your routine qPCR testing.

You define your in-house assay content. We prepare it in cleanroom conditions.

Stop mixing. Start testing.

Routine testing demands speed and reliability. However, pipetting multiple oligonucleotides and dispensing qPCR premixes is repetitive and error-prone. **De-risk your workflow** by outsourcing assay preparation at scale in **ready-to-use** format.

Benefit from our expertise in qPCR reagents to build your custom assay

We have been manufacturing high-quality oligonucleotides and state-of-the-art Master Mixes for more than 40 years. To **simplify your routine testing** with **in-house assays**, we premix them according to your specifications.

Whether you require **primers and probes premixed**, or a **full ready-to-use solution** including the master mix, we adapt to your needs.

Custom primers and probes

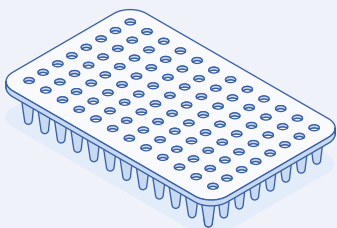
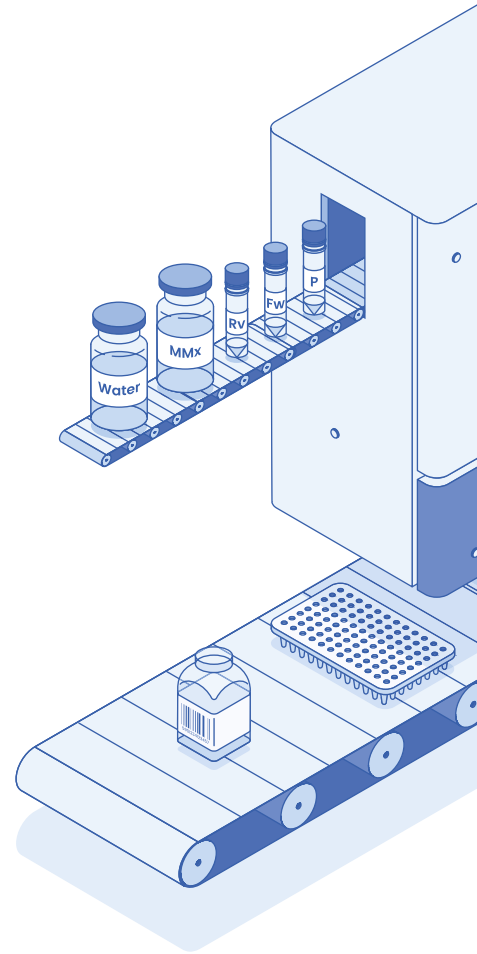
We manufacture custom oligonucleotides with tailored chemistries (DNA, LNA®) and modifications, including double-dye probes (e.g., TaqMan®, FRET), dyes and quenchers to support any assay design.

Enzymes and Master Mixes

We recommend Takyon® Master Mix with validated shelf life and tailored formulations. Combined with EuroScript reverse transcriptases, it also supports robust one-step RT-qPCR applications.

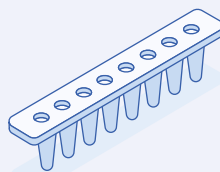
Plastics

We supply your premix in high-quality qPCR plastics and the format adapted to your workflow.



Plates

High-throughput routine testing



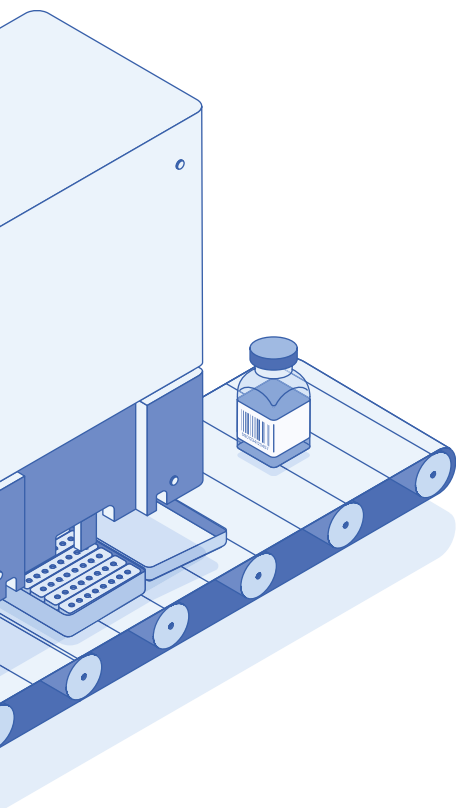
Strips

Modular or multi-assay testing



Tubes | Bottles

Bulk of premix fitting your liquid handler



Why rely on us for your assay preparation?

Save time

No more cumbersome pipetting and rework.
Go for an easy-to-use format.

Simplify your paperwork

1 lot | 1 QC | 1 doc.
Ease your lot and stock management.

Reduce risks

Cleanroom and automated assembly minimize contamination and pipetting error risks.

Keep your design

Stay in control of your assay content, with full design and format flexibility.

Secure your incoming QC

Premixes are tested for innocuity (clean NTC) and qPCR compatibility (positive IPC).

Control your budget

Reduce hidden costs, benefit from economy of scale & avoid costly investments.

Your assay is in good hands.

A dedicated project manager provides expert guidance and attentive support.

- Reviews designs and specifications
- Coordinates the project from manufacturing to delivery
- Ensures regular follow up and responsive answers
- Provides technical expertise when needed

For any type of testing



Human

Molecular assay for clinical diagnostics



Vet

Detect infectious or genetic diseases in animals



Food-feed

Screen pathogens, GMOs, or allergens



Environment

Monitor water, air, or soil contaminants.



Biopharma QC

Control raw materials or in-process samples

The quality expected for your assay

Cleanroom prep & dispensing

We premix your qPCR reagents in ISO-classified cleanrooms, minimizing contamination risks and variability. Our automated liquid handlers deliver precise, reproducible volumes to increase consistency.



Quality controls

We release each premix only after qPCR Quality Control. We verify that there is no contamination and confirm correct amplification with internal positive control. If needed, we can also test your own positive control. You receive your assay with fully documented QC results.



Want to maintain your LDTs under IVDR?

Stay compliant with confidence.

We manufacture your custom in-house IVD assays with documentation, traceability and quality assurance needed to support your IVDR exemption.



Scan this code and
Start your project with us!