



OEM Services

Custom Molecular Dx Assays

Benefit from our expertise, technologies and facilities to **accelerate the market entry of your nucleic acid-based kit.**

From individual **components** to **ready-to-use kits**

Take advantage of our strong manufacturing background in qPCR with reagents for **research, diagnostics and commercial use.**

We provide and assemble all the necessary components for efficient PCR, qPCR, NGS and other applications.



Up to
40^L
Batch volume



New generation of amplification reagents

Our innovative and robust reagents offer an optimal solution for your amplification assays.



IP-friendly dyes & quenchers

A customizable and attractive offer spanning the full visible spectrum, from blue to far-red-shifted wavelengths.



Assay preparation

With our flexible format and layout approach, we deliver your kit in the quality and design you defined.

Customize your Assay

We offer a **flexible service** to deliver your reagents and assays according to your application and needs.

Reagent Types

Primers, probes, proprietary enzymes, custom buffer, controls, and additives.

Plates, strips & tubes

Get access to a **large choice of plastics** from single tubes up to 96- and 384-well plates directly compatible with your platform.

Quantity

Your reagents and assays are prepared with the utmost accuracy **from small (5 µL) to large (40 liters)** custom batch size.

Labeling

With the labeling option, you can ask for a **laser marking** or a **barcoding** of your plastics.

Packaging

Your reagents can be delivered in a **ready-to-use** format or in **bulk**.



FLUORESCENT DYES | QUENCHERS | qPCR REAGENTS

IP-friendly and IP-free reagents and technologies

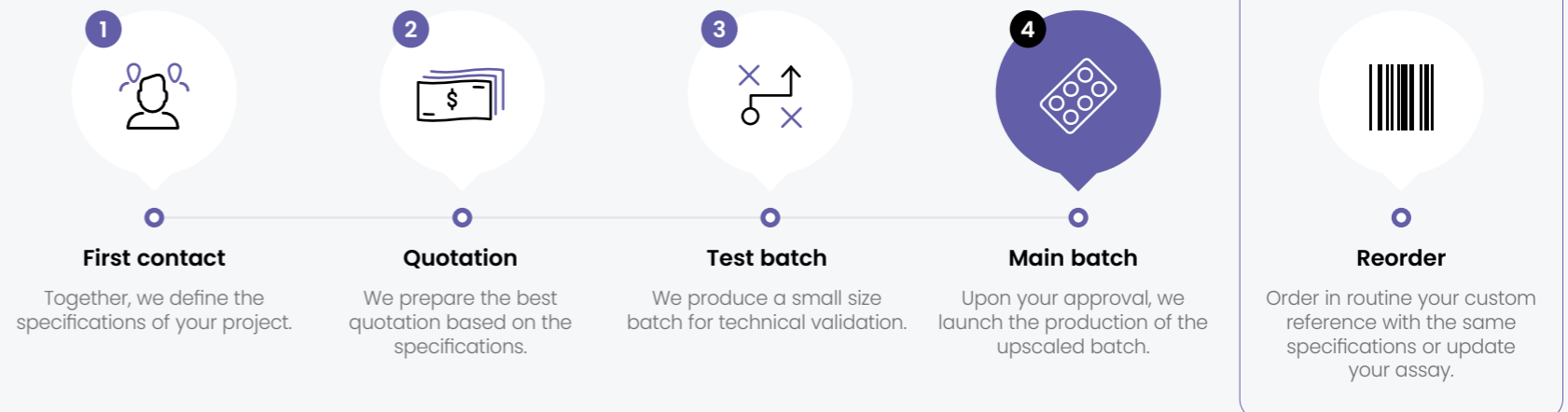
Avoid the burden of royalties reporting

The freedom to use the technologies or tradenames we offer is granted under a simple and convenient single-page license letter.



An easy and flexible manufacturing process

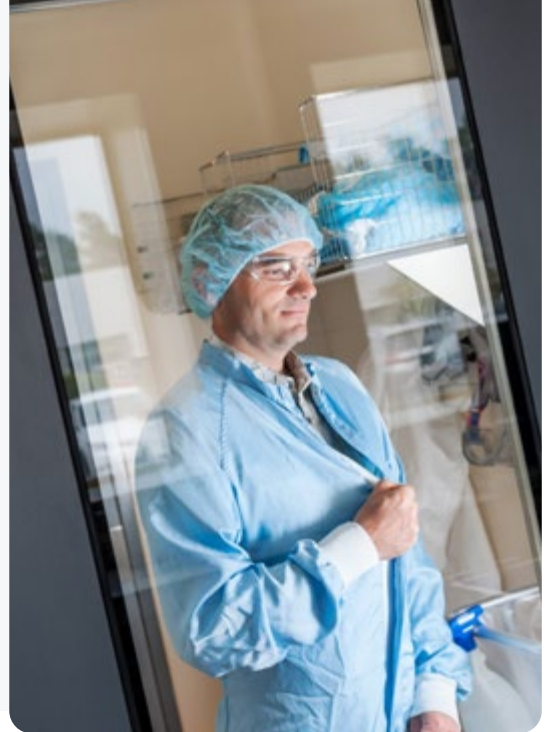
We partner with you from the initial contact to the release of your products. **Your dedicated project manager** supports you all along the process.



Quality all along the process

Our traceability and quality level are in line with IVDR CE-labeling.

- Our Quality Management System (QMS) is **ISO 9001** and **ISO 13485** certified.
- Our **clean rooms** meet ISO8 and ISO7 standards.
- All products are **quality controlled** under specific environments.
- **Tailored QC** available.
- Expertise in **CE-labeled** kits and components.



Your advantages at a glance



Be first on the market

Benefit from our expertise and technology platform to accelerate the release of your assay.



Increase your assay quality

Take advantage of our state-of-the-art equipment dispensing accurate and reproducible aliquots.



Avoid costly investment

Our validated facilities and trained staff are qualified to produce any custom nucleic-acid based diagnostic assay.



Focus on your business

You can keep developing innovative ideas while we take care of the manufacturing of your products.



Access innovative technologies

We commit to continuous investment to overcome any future challenges.



Get support from Experts

Since 2010, we have been refining our expertise and established multiple successful partnerships.



Scan this code and
Discuss your project with us!