

The full value of outsourcing in-house assay preparation

1

Time savings

- Mixing & dispensing handled for you
- Staff time freed for more demanding tasks
- Ready-to-use formats
- Easy-to-use (less training required)
- Reduced re-work
- Reduced batch validation frequency
- Simplified stock management

2

Risk Reduction

- ISO-7/8 cleanroom preparation (minimized contamination)
- Automated dispensing (increased lot-to-lot consistency)
- Professional labeling with UDI-ready traceability
- Stability tested formulations
- qPCR QC release before delivery
- One QC release document per lot (centralized documentation)

3

Cost efficiency

- No capital investment (liquid handler, cleanroom...)
- Economies of scale
- Less reagent waste

4

Full flexibility

- Full control over content, labeling and formulation
- Plates, strips, tubes, or bulk delivery
- Custom layout (for plates and strips)
- Adapt quickly to evolving needs
- Partial deliveries available on demand

5

Technical support

- Support in assay design if needed
- Platform & probes compatibility guidance
- Responsive troubleshooting assistance

Discover more on
www.eurogentec.com