

PCR & qPCR REAGENTS AND SERVICES



GRT-GEN-0509-V1

EUROGENTEC, YOUR UNIVERSAL PCR & qPCR SUPPLIER
ALL YOU NEED IS HERE, SIMPLY ASK!

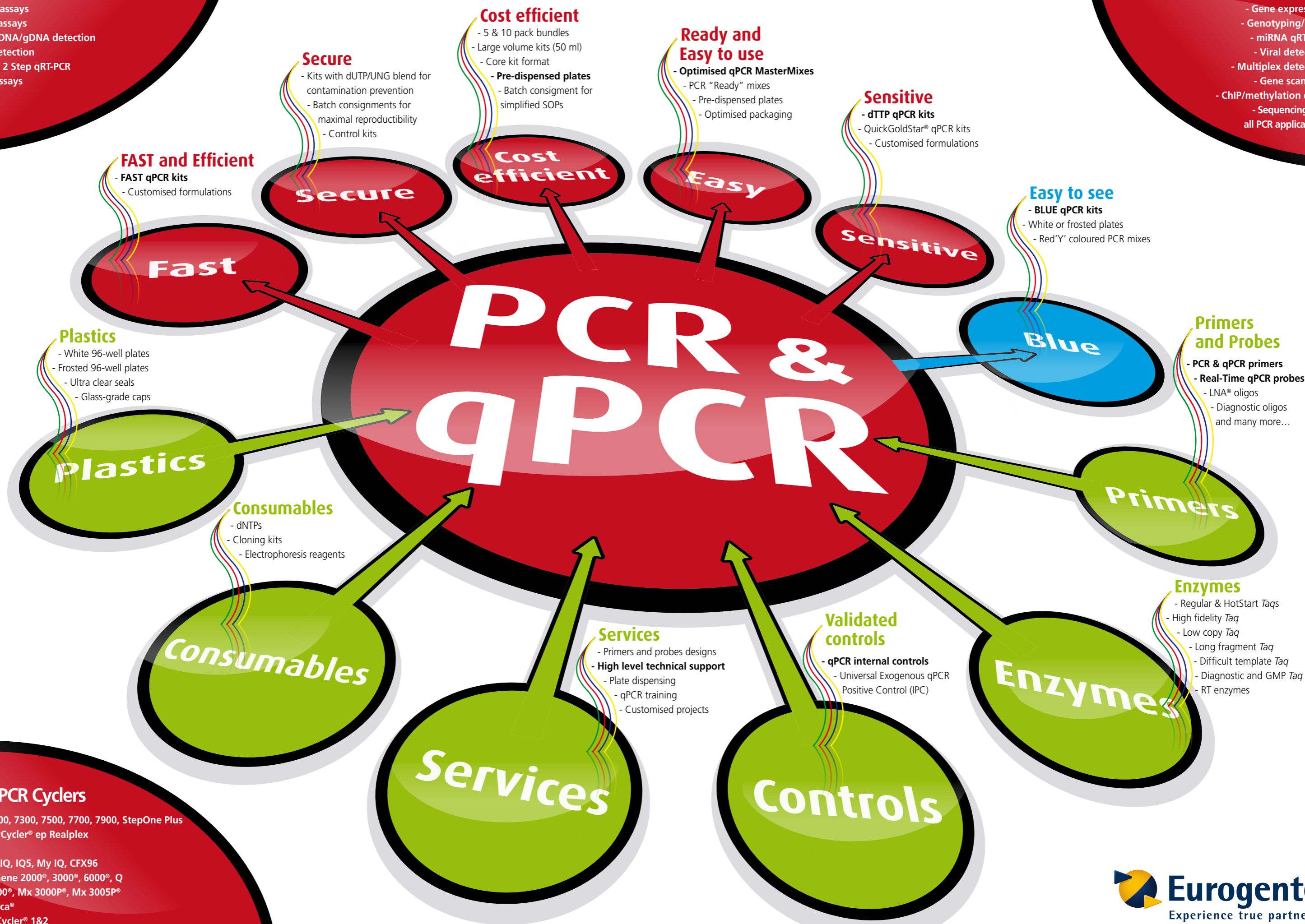
Eurogentec S.A.
LIEGE SCIENCE PARK
Rue du Bois St. Jean, 5
4102 Seraing, Belgium
Tel: +32 (0)4 372 74 00
Fax: +32 (0)4 372 75 00
info@eurogentec.com
www.eurogentec.com

All Techniques

- SYBR® assays
- Probe assays
- DNA/cDNA/gDNA detection
- RNA detection
- One or 2 Step qRT-PCR
- HRM assays

All Applications

- Gene expression
- Genotyping/SNPs
- miRNA qRT-PCR
- Viral detection
- Multiplex detection
- Gene scanning
- ChIP/methylation qPCR
- Sequencing and all PCR applications



All qPCR Cyclers

- ABI 7000, 7300, 7500, 7700, 7900, StepOne Plus
- MasterCycler® ep Realplex
- LC 480
- Icyler IQ, IQ5, My IQ, CFX96
- RotorGene 2000®, 3000®, 6000®, Q
- Mx 4000®, Mx 3000P®, Mx 3005P®
- Quantica®
- SmartCycler® 1&2
- Opticon® 1&2, Chromo 4, MiniOpticon