



Takyon™

Amplification Technologies

qPCR • RT-qPCR • PCR

For Life Science Research

- > qPCR/RT-qPCR Takyon™ kits
- > qPCR control assays
- > qPCR plastics
- > PCR kits and reagents
- > Custom formulations





Highly consistent
quality reagents



Clear detection of
low copy targets



Maximum **reproducibility**
of results



Assay **dispensing services**
for process automation



Custom formulations
for demanding projects



Innovative dry format

RELIABLE TOOLS FOR AMPLIFICATION, DETECTION AND QUANTIFICATION OF NUCLEIC ACIDS

For over a decade Eurogentec has forged a strong reputation for developing trusted tools for qPCR and PCR applications. We bring you solutions, from individually performing reagents and consumables to ready-to-use kits.

qPCR/RT-qPCR/PCR

qPCR
& RT-qPCR

PCR

Custom
services

Primer
& Probes

Plastics

- p5 *From PCR to qPCR*
- p6 **qPCR/RT-qPCR range**
- p6 *Takyon™ Kits for qPCR & RT-qPCR assays*
- p12 *qPCR control assays*
- p14 **PCR range**
- p14 *Ready-to-use PCR Mixes*
- p15 *PCR Reagents*
- p18 **Custom services**
- p18 *Custom formulations*
- p18 *Assay dispensing service*
- p18 *Kit design and manufacturing*
- p19 **Primers & Probes**
- p19 *Primers & probes for qPCR assays*
- p23 **qPCR plastics**

Annexes

- p29 *License statements*
- p29 *Trademarks/labels*



FROM PCR TO qPCR

Since its creation in 1983, the PCR has gained interest in molecular identification, sequencing and genetic engineering and is no more restricted to the amplification of low amount targets. However the PCR technology remains an end-point method which still requires a downstream processing after amplification (e.g.: a running gel) to detect and identify the amplified product.

In contrast Real-time PCR technology relies on the signal intensity of a fluorescent reporter dye which increases proportionally with the sample amount. The reaction progress is then monitored in live in closed tubes, significantly reducing hands-on time and contamination risks.

Real-time PCR is frequently used in research and diagnostic assays to precisely detect or quantify the starting template copy number. Result accuracy is also requested for demanding applications such as gene expression and miRNA analysis, Single Nucleotide Polymorphism (SNP) genotyping... and is mainly linked to the reproducibility of the assay.

Different parameters such as the buffer composition or the efficacy of the polymerase may influence result quality and consistency. To ensure successful qPCR reactions, Eurogentec has developed various kit components including a proprietary Takyon™ fusion enzyme and performing reagents as well as efficient consumables you can rely on durably.

- qPCR/RT-qPCR/PCR
- qPCR & RT-qPCR
- PCR
- Custom services
- Primer & Probes
- Plastics

qPCR/RT-qPCR range

Inspired by the faster than light particle (Tachyon), the Takyon™ qPCR line from Eurogentec has been designed to shorten your qPCR runs. Thanks to a better affinity to DNA, and combined to an optimised reaction buffer, the new Takyon™ fusion polymerase ensures sensitivity and fast delivery of accurate and reproducible results. In this chapter, you will find ready-to-use kits, plastics, controls and reagents to achieve your qPCR and RT-qPCR assays successfully. ■

BENEFITS OF TAKYON™ KITS

Universal capabilities

The Takyon™ kits are compatible with all cyclers and applications including high throughput and on the field testing.

Large dynamic range

The Takyon™ kits give consistent amplification results over a large range of initial target concentrations

Robust formulation

The fusion polymerase is highly efficient and less sensitive to inhibitors. This allows amplification with some difficult templates like biological fluids and viruses.

Small reaction volumes

The high sensitivity of the Takyon™ kits combined to the highly concentrated 5X format give reliable results even in small volume reaction.

Suitable for multiplex

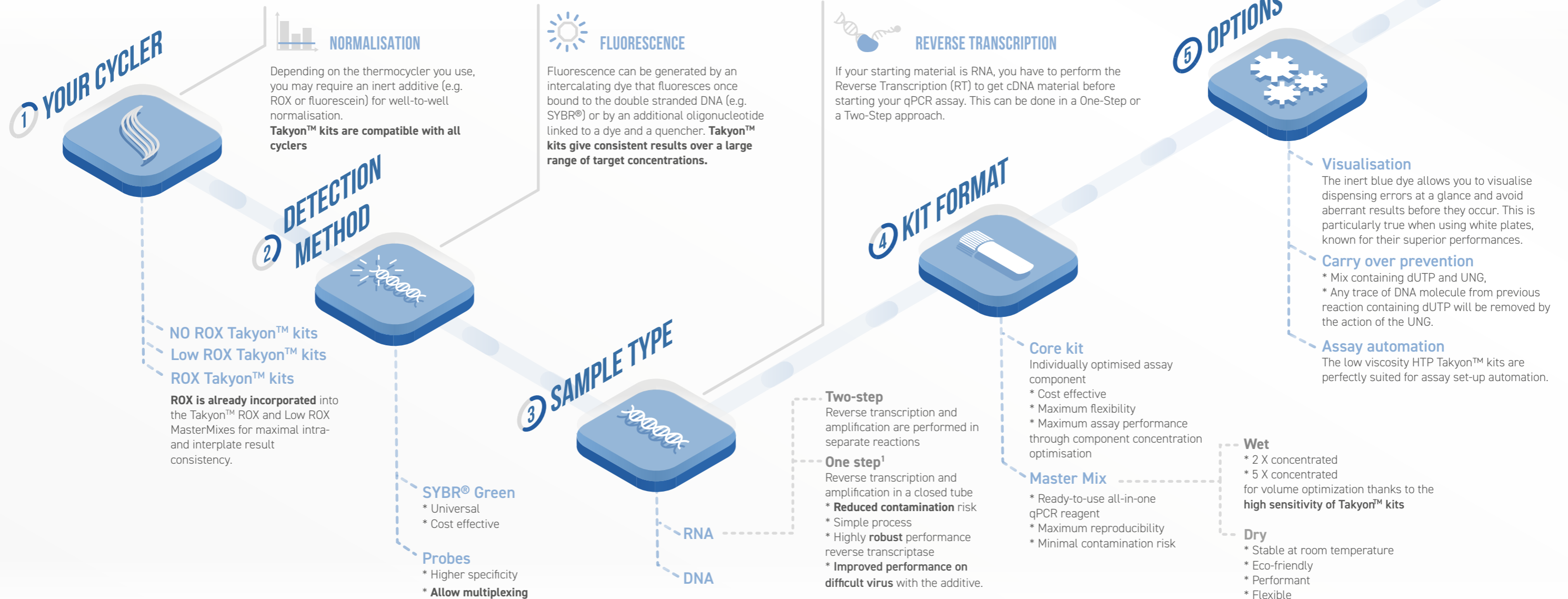
The Takyon™ kits allow the detection of multiple targets in a single reaction tube.

GOOD TO KNOW

To help you in your kit selection, a convenient **online qPCR selector** is available on <http://www.eurogentec.com/qpcr-selector.html>. You can also evaluate a sample of any MasterMix for free and benefit from a 50% discount on your first kit order.

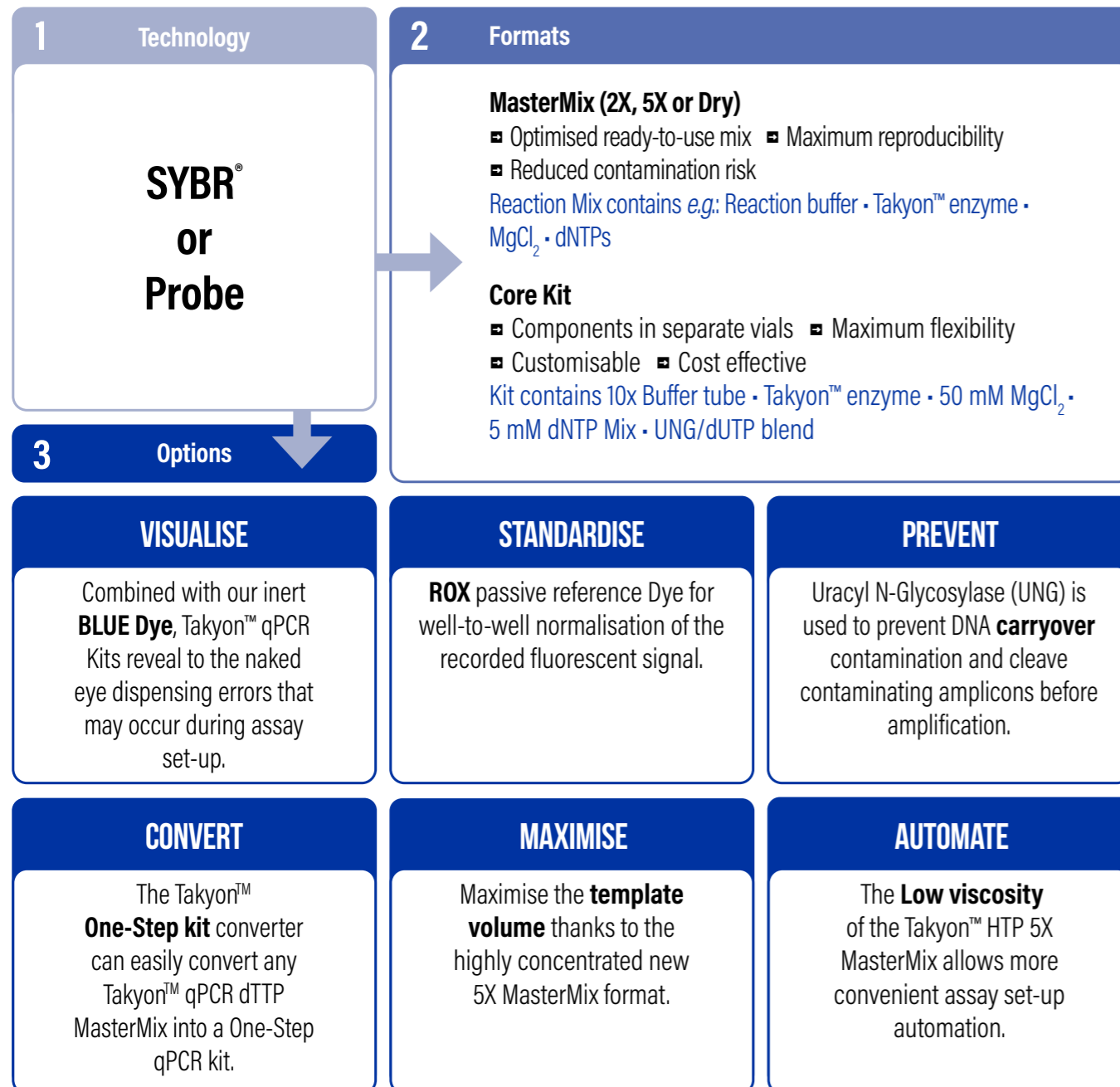


TAKYON™ KITS FOR qPCR & RT-qPCR ASSAYS



1. Any Takyon™ dTTP Master Mix for qPCR assays can easily be converted into a one-step RT-qPCR kit by using the Takyon™ **One-Step kit converter** simply combining the two.

GLOBAL VIEW OF TAKYON KITS



FIND THE PRODUCT MATCHING YOUR THERMOCYCLER

Some thermocyclers use fluorescent dyes to allow well-to-well normalisation. To meet any option, Eurogentec offers kits with passive references (ROX or fluorescein) already incorporated into the mix but also kits without passive reference**.

The lexicon below guides you to the right ordering information in tables.

Cycler Group	Cycler Group	Cycler Group
A	G	P
<ul style="list-style-type: none"> ABI Prism[®] 5700 (p.9) ABI Prism[®] 7000 (p.9) ABI Prism[®] 7300 (p.9) ABI Prism[®] 7500 & FAST 7500 (p.9) ABI Prism[®] 7700 (p.9) ABI Prism[®] 7900 & FAST 7900 (p.9) ABI Step One & Plus (p.9) ABI ViiATM⁷ (p.9) AriaMx (p.10) 	<ul style="list-style-type: none"> GeneAmp[®] SDS 5700 (p.9) iCycler iQ[®] (p.10) iQTM5 (p.10) LC96 & LC480 (p.10) LightCycler[®] Nano System (p.10) Mastercycler[®] ep realplex I & II (p.10) MiniOpticon (p.10) Mx3000P[®] (p.9&10) Mx3005P[®] (p.9&10) Mx4000[®] (p.9&10) My iQ (p.10) 	<ul style="list-style-type: none"> PiKoReal[™] (p.10) Quantica[®] (p.10) QuantStudio[®] (p.9) qTower (p.10) Rotor-Gene[®] 2000/3000/6000/Q (p.10) SmartCycler[®] 1 & 2 (p.10) Swift Spectrum[™] 96 (p.10)
C	L	R
<ul style="list-style-type: none"> CFX384 (p.10) CFX96 (p.10) Chromo 4 (p.10) 		
D	M	S
<ul style="list-style-type: none"> DNA Engine Opticon[®] 1 & 2 (p.10) 		

Legend

- ROX
- Low ROX
- No ROX
- Use Fluorescein with SYBR[®] MMix or No ROX with Probe MMix

RT-qPCR kits

The Takyon™ One-Step MasterMixes contain the new EuroscriptII RT which delivers highly robust performance on classical mRNAs, small RNA and viral RNAs.

The new additive improves significantly the amplification performance on difficult virus types as well as on FFPE samples.

The 5X highly concentrated format allows more flexibility in template volume use especially in the context of small volume reaction.

Product name	1 pack	5 packs	10 packs	# RXNs per single kit ¹	Format	Full carryover (dUTP+UNG)	Corresponding sample
SYBR[®] Assays							
For Two-Step Assays combine any kit mentioned above to our reverse transcriptase core kit RT-RTCK-03							
For One-Step Assays combine any dTTP Takyon™ 2X MasterMix with the Takyon™ One-Step kit converter UF-RTAD-D0101 (for 150 RXN MMix) or UF-RTAD-D0701 (for 750 RXN MMix)							
Probe Assays							
Takyon™ One-Step ROX Probe 5X MasterMix dTTP	UF-RP5X-RT0501	UF-RP5X-RT0505	UF-RP5X-RT0510	1250	MasterMix 5x		UF-RP5X-RT0100
Takyon™ One-Step Low ROX Probe 5X MasterMix dTTP	UF-LP5X-RT0501	UF-LP5X-RT0505	UF-LP5X-RT0510	1250	MasterMix 5x		UF-LP5X-RT0100
Takyon™ One-Step No ROX Probe 5X MasterMix dTTP	UF-NP5X-RT0501	UF-NP5X-RT0505	UF-NP5X-RT0510	1250	MasterMix 5x		UF-LP5X-RT0100

¹ #RXNs is calculated for a final volume of 20 µL.

² Optional: when dUTP/UNG (RT-UTPUNG-020) additive is used. Component included in Core Kits.

³ Combine with fluorescein additive (RT-FLUO-ADD).

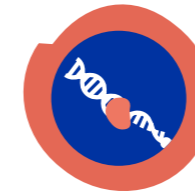
qPCR kits

Product name	1 pack	5 packs	10 packs	# RXNs per single kit ¹	Format	Full carryover (dUTP+UNG)	Corresponding sample
SYBR® Assays							
ROX	Takyon™ ROX SYBR® MasterMix 2X Blue dTTP 75 mL	UF-RSMT-B0701	UF-RSMT-B0705	UF-RSMT-B0710	750	MasterMix 2x	² UF-RSMT-B0100
	Takyon™ ROX SYBR® 2X Core Kit Blue dTTP	UF-RSCT-B0201	UF-RSCT-B0205	UF-RSCT-B0200	1250	Core Kit	²
LOW ROX	Takyon™ Low ROX SYBR® 2X MasterMix blue dTTP	UF-LSMT-B0701	UF-LSMT-B0705	UF-LSMT-B0710	750	MasterMix 2x	²
NO ROX	Takyon™ No ROX SYBR® 2X MasterMix Blue dTTP 75 mL	UF-NSMT-B0701	UF-NSMT-B0705	UF-NSMT-B0710	750	MasterMix 2x	² UF-NSMT-B0100
	Takyon™ No ROX SYBR® Core Kit blue dTTP	UF-NSCT-B0201	UF-NSCT-B0205	UF-NSCT-B0210	1250	Core Kit	²
FLUORESCEN	Takyon™ SYBR® 2X MasterMix blue dTTP with fluorescein	UF-FSMT-B0701	UF-FSMT-B0705	UF-FSMT-B0710	750	MasterMix 2x	²
	Takyon™ No ROX SYBR® Core Kit blue dTTP ³	UF-NSCT-B0201	UF-NSCT-B0205	UF-NSCT-B0210	1250	Core Kit	²
Probe Assays							
ROX	Takyon™ ROX Probe 2X MasterMix UNG 75 mL	UF-RPMU-C0701	UF-RPMU-C0705	UF-RPMU-C0710	750	MasterMix 2x	✓ UF-RPMU-C0100
	Takyon™ ROX Probe 2X MasterMix dTTP Blue 75 mL	UF-RPMT-B0701	UF-RPMT-B0705	UF-RPMT-B0710	750	MasterMix 2x	² UF-RPMT-B0100
	Takyon™ ROX Probe 2X MasterMix dTTP 75 mL	UF-RPMT-C0701	UF-RPMT-C0705	UF-RPMT-C0710	750	MasterMix 2x	² UF-RPMT-C0100
	Takyon™ ROX Probe 5X MasterMix dTTP 5 x 1 mL	UF-RP5X-C0501	UF-RP5X-C0505	UF-RP5X-C0510	1250	MasterMix 5x	² UF-RP5X-C0100
	Takyon™ ROX Probe Core Kit dTTP	UF-RPCT-C0201	UF-RPCT-C0205	UF-RPCT-C0210	1250	Core Kit	²
LOW ROX	Takyon™ Low ROX Probe 2X MasterMix UNG 75 mL	UF-LPMU-C0701	UF-LPMU-C0705	UF-LPMU-C0710	750	MasterMix 2x	✓ UF-LPMU-C0100
	Takyon™ Low ROX Probe 2X MasterMix dTTP Blue 75 mL	UF-LPMT-B0701	UF-LPMT-B0705	UF-LPMT-B0710	750	MasterMix 2x	² UF-LPMT-B0100
	Takyon™ Low ROX Probe 2X MasterMix dTTP 75 mL	UF-LPMT-C0701	UF-LPMT-C0705	UF-LPMT-C0710	750	MasterMix 2X	² UF-LPMT-C0100
	Takyon™ Low ROX Probe 5X MasterMix dTTP 5 x 1 mL	UF-LP5X-C0501	UF-LP5X-C0505	UF-LP5X-C0510	1250	MasterMix 5x	² UF-LP5X-C0100
	Takyon™ No ROX Probe 2X MasterMix UNG 75 mL	UF-NPMU-C0701	UF-NPMU-C0705	UF-NPMU-C0710	750	MasterMix 2x	✓ UF-NPMU-C0100
NO ROX	Takyon™ No ROX Probe 2X MasterMix dTTP BLUE 75 mL	UF-NPMT-B0701	UF-NPMT-B0705	UF-NPMT-B0710	750	MasterMix 2x	² UF-NPMT-B0100
	Takyon™ No ROX Probe 2X MasterMix dTTP 75 mL	UF-NPMT-C0701	UF-NPMT-C0705	UF-NPMT-C0710	750	MasterMix 2x	² UF-NPMT-C0100
	Takyon™ No ROX Probe 5X MasterMix dTTP 5 x 1 mL	UF-NP5X-C0501	UF-NP5X-C0505	UF-NP5X-C0510	1250	MasterMix 5x	² UF-NP5X-C0100
	Takyon™ No ROX Probe Core Kit dTTP	UF-NPCT-C0701	UF-NPCT-C0705	UF-NPCT-C0710	1250	Core Kit	²
	Takyon™ No ROX Probe Core Kit dTTP blue	UF-NPCT-B0701	UF-NPCT-B0705	UF-NPCT-B0710	1250	Core Kit	²

¹ #RXNs is calculated for a final volume of 20 µL.
² Optional: when dUTP/UNG (RT-UTPUNG-020) additive is used. Component included in Core Kits.
³ Combine with fluorescein additive (RT-FLUO-ADD).



Additional options

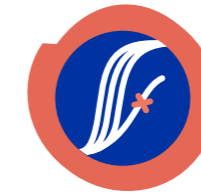


REVERSE TRANSCRIPTION

When working with RNA samples, a reverse transcription is required prior the qPCR reaction. One-step Takyon™ kits containing reverse transcriptase are available for all cyclers. In addition, you can convert any dTTP qPCR Takyon™ kit into a:

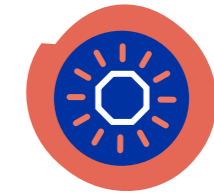
One-step RT-qPCR kit: Use Takyon™ One-Step Kit converter with any dTTP Takyon™ 2X Master Mix to make reverse transcription and qPCR in the same tube.

Two-step RT-qPCR kit: Use the Reverse Transcriptase Core Kit prior the qPCR assay to make a two steps reactions with any qPCR kit.



PREVENT CONTAMINATION

Using a mix containing dUTP and UNG brings security to the assay. Indeed any trace of DNA molecule from previous reaction containing dUTP will be removed by the action of the Uracyl-N-Glycosylase (UNG). Takyon™ 2X Master Mixes for probe assays are available with dUTP/UNG already incorporated. Takyon™ core kits contain dUTP and UNG. Otherwise, dUTP/UNG additive can be used to convert any dTTP kit (e.g. SYBR® kit).



PASSIVE REFERENCE

Passive reference is required for well-to-well normalisation on some thermocyclers. Takyon™ kits with passive reference (ROX or fluorescein) already incorporated into the mix are available. ROX and Fluorescein additives are available as stand-alone reagents.



MISCELLANEOUS

ORDERING INFORMATION

Description	References	Quantity
Takyon™ One-Step Kit converter 20 µl	UF-RTAD-D0101	150 RXNs
	UF-RTAD-D0701	750 RXNs
	UF-RTAD-D0705	5x 750 RXNs
	UF-RTAD-D0710	10x 750 RXNs
Reverse Transcriptase Core Kit 300	RT-RTCK-03	300 RXNs
	05-RTCK-03	5x 300 RXNs
	10-RTCK-03	10x 300 RXNs
	RT-RTCK-05	500 RXNs
Reverse Transcriptase Core Kit 500	05-RTCK-05	5x 500 RXNs
	10-RTCK-05	10x 500 RXNs
	RT-UTPUNG-005	40 µL (150 RXNs)
dUTP/UNG Additive for dTTP kit conversion	RT-UTPUNG-020	200 µL (750 RXNs)
	05-UTPUNG-020	5x 200 µL (5x 750 RXNs)
	10-UTPUNG-020	10x 200 µL (10x 750 RXNs)
dUTP/UNG Additive for 5X MMx kit conversion	RT-UTPUNG-033	1250 rxn

ORDERING INFORMATION

Description	References	Quantity
Fluorescein additive 1 mM (for My iQ / i-Cycler iQ ⁺ / iQ ⁵)	RT-FLUO-ADD	1 mL
ROX Passive Reference	RT-PARE-03	300 µl
	05-PARE-03	5x300 µl
	10-PARE-03	10x300 µl
	RT-PARE-100	10 mL
UNG (Uracil-N-Glycosylase)	RT-0610-03	300 u
	05-0610-03	5x300 u
	10-0610-03	10x300 u
UNG (Uracil-N-Glycosylase)	RT-0610-15	1500 u
	05-0610-15	5x1500 u
	10-0610-15	10x1500 u
Takyon™ Blue Additive	UF-BLUE-ADD	3750 RXN

qPCR CONTROL ASSAYS

qPCR is the method of choice for quantitative gene expression studies. However qPCR assays are prone to inhibition by various substances found in many samples (clinical, soil, plant...).

Internal Positive Controls (IPC) allow you to detect inhibited amplification reactions

due to the carryover of reagents used for the isolation of nucleic acids. Other causes of false-negative results include target nucleic acid degradation, sample processing errors and thermocycler malfunction.

Housekeeping controls allow for gene expression studies. Reliable determination

of the gene expression level across samples requires efficient normalisation of the recorded signals. For that purpose Eurogentec has developed a limited set of optimised and validated housekeeping controls.

Housekeeping internal controls

The accurate and reliable determination of gene expression is commonly achieved by using well characterised internal genes as reference.

Accordingly, Eurogentec has functionally validated each of its controls. Each set has been optimised to a PCR efficiency of 95-100% when combined with a qPCR kit from Eurogentec.

The closer the level of target expression is to the expression of the internal control gene, the more accurate the determination

will be. Therefore the selected control gene should match your target gene expression depending on the cell type or tissue used. The table here below give you an idea of the expression levels of a selection of human housekeeping genes.

ALSO AVAILABLE

You can request your **custom control design** for any other selected gene by sending the gene details to: scientific.support@eurogentec.com

FEATURES

- Validated for human samples
- Sensitive and specific detection
- Full expression range coverage
- Compatible with SYBR® green or probe assays



ORDERING INFORMATION

Human Gene	Relative expression level	cDNA	gDNA	Organism	Amplicon length	#RXN (50µL)	Reference
18S rRNA	Very high	✓	✓	H, Ra, M, Rb	121 bp	250	RT-CKFT-18S RT-CKYD-18S
28S rRNA	Very high	✓	✓	H, M	87 bp	250	RT-CKYD-28S
β-actin	High	✓	✓	H	91 bp	250	RT-CKYD-ACTB
GAPDH	High	✓		H	86 bp	250	RT-CKYD-GAPD
Ubiquitin C	High	✓	✓	H	76 bp	250	RT-CKYD-UBC
β-2-microglobulin	High	✓	✓	H	112 bp	250	RT-CKYD-B2M
Phospholipase A2	Medium	✓		H	93 bp	250	RT-CKYD-YWHAZ
RPL13a	Medium	✓	✓	H	92 bp	250	RT-CKYD-RPL13A
SDHA	Medium	✓		H	118 bp	250	RT-CKYD-SDHA
HMBS	Medium	✓		H	119 bp	250	RT-CKYD-HMBS
HPRT1	Low	✓		H	100 bp	250	RT-CKYD-HPRT1
TBP	Low	✓		H	145 bp	250	RT-CKYD-TBP

Relative expression levels of a selection of human housekeeping genes
H: Human Ra: Rat M: Mouse Rb: Rabbit

Positive Controls

Eurogentec's Positive Controls (PC) are optimised Taqman® assays designed to distinguish true target negatives from false negatives.

The **Sample Processing Control (SPC)** is spiked into samples before extraction to monitor poor extraction yield, PCR

inhibition, incorrect pipetting or cycling parameters.

The **Internal Positive Control (IPC)** is spiked into samples before qPCR assay to confirm the absence of target sequence, inhibition or a reaction set-up/cycling error.



ORDERING INFORMATION

Product	#RXN (50µL)	Thermocycler channel (see page 24)	Reference	
SPC	Sample Processing Control, YY-TAMRA	200	VIC/YY	RT-SPCY-T02
		1000	VIC/YY	RT-SPCY-T10
	Sample Processing Control, YY-BHQ-1*	200	VIC/YY	RT-SPCY-B02
		1000	VIC/YY	RT-SPCY-B12
IPC	Sample Processing Control, HiLyte™647-QXL® 670	200	Cy5	RT-SPCC-Q02
		1000	Cy5	RT-SPCC-Q10
	Universal Exogenous qPCR Positive Control, YY-TAMRA	200	VIC/YY	RT-IPCY-T02
		1000	VIC/YY	RT-IPCY-T10
IPC	Universal Exogenous qPCR Positive Control, YY-BHQ-1*	200	VIC/YY	RT-IPCY-B02
		1000	VIC/YY	RT-IPCY-B10
	Universal Exogenous qPCR Positive Control, HiLyte™647-QXL® 670	200	Cy5	RT-IPCC-Q02
		1000	Cy5	RT-IPCC-Q10

BENEFITS

Specific

Does not interfere with target amplification

Avoid amplification of endogenous genes

Sensitive

Compatible with low copy templates

Available with dark quencher for maximal signal-to-noise ratio

Designed for multiplex assays

Detected in the Yakima Yellow®/VIC®/JOE channel

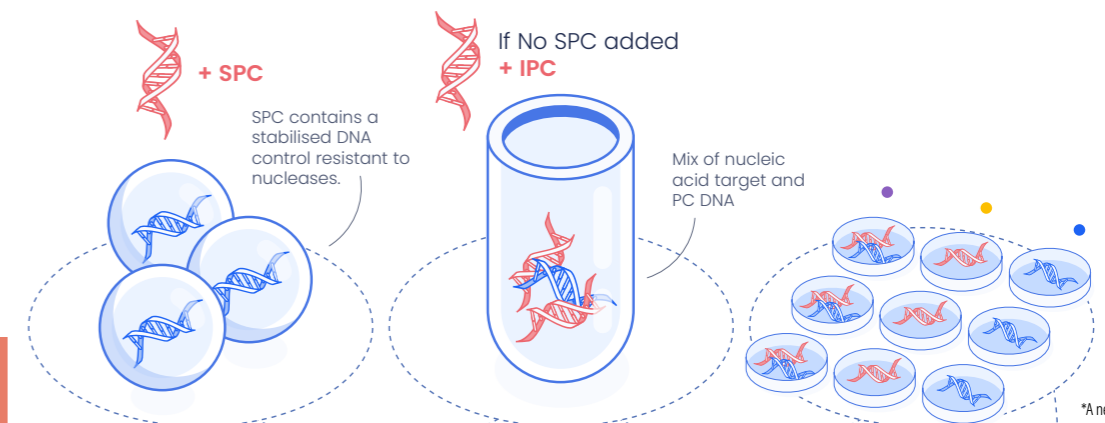
Validated in duplex assays with the target in the FAM channel

Custom labelling is available on request

METHOD

2 controls (SPC or IPC), multiple possibilities

Eurogentec's Positive Controls (PC) are Optimised Taqman® Assays



APPLICATIONS

Distinguish true negative from false negative targets resulting from inhibition.

Monitor loss or degradation of your DNA target during extraction.

Assess the quality of your nucleic acids extracts.

Calibrate interplate runs for data normalisation.

1 Extraction

2 Example of assay set-up

3 Results

Amplification	Results interpretation	
Sample + PC		
+	+	Target Positive
-	+	Target Negative (no target sequence)
-	-	No conclusion on target presence*

A negative PC (IPC or SPC) amplification suggests the presence of inhibitors, a wrong PCR set-up, a defective mix, an incorrect thermal cycling protocol or an extraction failure (in case of SPC use). YY = VIC* equivalent.

● qPCR Assay + Control: Sample + Primers + Probe (FAM) PC DNA + Primers + Probe (YY*)

● No Template Control: No Sample + Primers + Probe (FAM) PC DNA + Primers + Probe (YY*)

● No Positive Control: Sample + Primers + Probe (FAM) No PC DNA + Primers + Probe (YY*)

qPCR/RT-qPCR/PCR

qPCR
& RT-qPCR

PCR

Custom
servicesPrimer
& Probes

Plastics

PCR range

The wide variety of PCR applications requires occasional adjustments to the PCR process. Optimising PCR for your specific application can mean the difference between success and failure. Eurogentec offers you the choice between ready-to-use PCR mixes or separated reagents you can tailor to your need. ■

READY TO USE PCR MIXES

To gain time, prevent contamination risk by multiple pipettings and increase reproducibility of your PCR reactions, we recommend the use of a ready-to-use PCR Mix. **2 x concentrated PCR mixes** are convenient premixed solutions that include in a single tube all common reagents needed for PCR. Simply dilute twice the PCR mix with template, primers and water to start your reaction.

Four PCR mixes are available:

- **GoldStar® Mix** is composed of GoldStar® *Taq* polymerase, dNTPs, MgCl₂ and buffer

- **Red'y'Gold Mix** has the same composition than the GoldStar® Mix except that the loading buffer is **coloured with a red loading dye** allowing the direct loading of the amplicons on a gel.

- **HotGoldStar Mix** is the HotStart version of the GoldStar® Mix, containing a HotGoldStar *Taq* polymerase.

- **Red'y'Star Mix** has a composition similar to the HotGoldStar® Mix with a **buffer coloured with a red loading dye** allowing the direct loading of the amplicons on a gel.

BENEFITS

Easy-to-use

Optimised for a broad range of targets (< 3 Kb) and GC rich sequences

Highest sensitivity and yield

Achieves the highest quality sequencing data

ADDITIONAL INFORMATION

Package Content

5 x 1 mL of PCR Mix comprising of *Taq*® polymerase (1 u/25µl), dNTPs (400 µM), MgCl₂ (3 mM) and buffer

cGMP Mixes

Mixes prepared under ISO 13485 compliant processes (including enzymes of the Diamond *Taq*® family) are available on request at diagnostic. taq@eurogentec.com



PCR REAGENTS

dNTPS

dNTPs are available as convenient **100 mM sets** or in **ready-to-use mixes**. They offer reliable and reproducible performance in any enzymatic reaction, including PCR and reverse transcription.

dNTPs are enzymatically produced from **nucleotide monophosphates** by a process of phosphorylation. This process uses highly specific enzymatic systems, which eliminate impurities and PCR inhibitors such as modified nucleotides and

pyrophosphates. dNTPs are purified with preparative HPLC and are at least **99 % pure**. They are supplied **ready-to-use** as Lithium salts in water at pH 7.0.

Lithium salts have greater resistance to repeated freezing and thawing cycles than sodium salts dNTPs, and remain sterile over the entire shelf-life due to the bacteriostatic activity of lithium towards various microorganisms. ■



BENEFITS

> 99% Pure – free from PCR inhibitors, DNase, RNase and Nickase

Remain sterile over the entire shelf-life

High resistance to repeated freeze-thaw cycles



ORDERING INFORMATION

Description	Quantity	#Ref
dNTP Mix 1 x 20 µmoles	1 x 1 ml	NU-0010-10
dNTP Mix 5 x 20 µmoles	5 x 1 ml	NU-0010-50
dNTP Mix 10 x 20 µmoles	10 x 1 ml	NU-0010-100
dNTP Set 4 x 25 µmoles	4 x 250 µl	NU-0020-50

PACKAGE CONTENT

dNTP mix contains all four nucleotides (dATP, dCTP, dGTP, dTTP) in a premixed solution, ready for immediate use (5 mM each).

dNTP set is delivered as a pack of four vials, each containing 100 mM of one of the four nucleotides (dATP, dCTP, dGTP, dTTP).



Reverse transcriptase

MU-MLV REVERSE TRANSCRIPTASE

Mu-MLV is a reverse transcriptase purified from a recombinant overexpressing clone of *E. coli* that carries the reverse transcriptase gene from Moloney Murine Leukemia Virus. The enzyme synthesises a complementary cDNA strand from a single stranded RNA or DNA template. Mu-MLV lacks the 3'→5' exonuclease activity. It is recommended to use 10 U/μg RNA for optimum reverse transcription.



ORDERING INFORMATION

Description	Quantity (units)	Product Numbers
Mu-MLV reverse transcriptase	40000 U	ME-0125-400

BENEFITS

Lower RNase H activity than AMV (Avian Myeloblastosis Virus) reverse transcriptase
 Suitable for long mRNAs templates (>5kb)

ADDITIONAL INFORMATION

Package Content

Mu-MLV reverse transcriptase (200 U/μl)
 5x cDNA buffer: 250 mM Tris-HCl (pH 8.3), 375 mM KCl, 50 mM DTT, 15 mM MgCl₂

Storage & Shipping conditions

Shipped and Stored at -20°C

Quality

The enzyme is tested for the absence of endonucleases and RNases



Custom services

Each project is different and a unique qPCR solution suiting all requirements does not exist. Eurogentec affords the flexibility to adapt its offer to your specific needs.



Primers & Probes



CUSTOM FORMULATIONS

Because setting up a qPCR or a PCR assay can be a real challenge requiring an optimised Mix composition, Eurogentec offers you the possibility to produce your own Mix.

Combine the following advantages in a single product:

- Maximised sensitivity and specificity of your custom validated Mix
- Absolute reproducibility with the benefits of a unique dedicated batch
- Minimal contamination risks and pipetting errors with ready-to-use Mix
- Consignment and scheduled deliveries
- Various packaging (tube, plate, bottles...)
- Inclusion of external components (enzymes, oligonucleotides...)

Please contact us to develop your own formulation:

custom.qpcr@eurogentec.com



ASSAY DISPENSING SERVICE

In its constant search for simple, affordable and innovative solutions for the Life Science community, Eurogentec has developed a dispensing service that brings all advantages of automation within everyone's reach.

Ease your workflow by automating 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, saving set-up time, minimising user input and skill level, as well as reagent wastage and hidden costs.

The process offers a very high reliability, reproducibility and accuracy. Our dispensing service is in line with ISO15189 quality standards and is fully scalable, taking care of the preparation of any type of routine assays to full kitting solutions.

KIT DESIGN AND MANUFACTURING

A unique offer allowing rapid RUO launch to full kit design and manufacturing, for reduced time-to-market.

We take in charge the preparation of your kits and provide them labelled with your own brand. We offer a flexible production with custom fill & finish release options to fit with your own product lines.

Contact for dispensing & kit design manufacturing: dispensing@eurogentec.com

PRIMERS & PROBES FOR qPCR ASSAYS

Primers

Primers are used to amplify DNA fragment by PCR/qPCR process. Primer sequences need to be chosen to uniquely select for a region of DNA, avoiding the possibility of mishybridisation to a similar unwanted sequence nearby as well as region harbouring highly stable secondary structures.

GOOD TO KNOW

For optimal results in multiplex assays or using SYBR® Green, we recommend RP-cartridge*Gold purification for your primer.

Probes

A probe is a nucleic acid fragment labelled with an internally quenched fluorophore whose fluorescence is restored when hybridised to its target sequence or when physically separated from the quencher due to nuclease activity. Eurogentec offers various probe types (Double Dye, Molecular Beacons, MGB...) with a wide range of fluorophore and quencher pairs covering all the qPCR channels. The use of probes increases the specificity of the assay and allows multiplexing. Please refer to the compatibility table p22. to select the dye adapted to your thermocycler.

MGB PROBES

Eurogentec provides high quality MGB probes perfectly suited for patient management** and available in IVD grade upon request. We provide a complete offer with more than 15 dyes covering all qPCR channels.

PROBES SPECIFICATIONS

Length: From 10 to 50 bases

Synthesis scale: 10 nmol • 40 nmol • 200 nmol • 1000 nmol • 2.5 µmol • 5 µmol • 10 µmol*

Backbone: DNA, LNA®, 2'O-Me RNA and phosphodiester linkage

Modifications: 5': 6-FAM, HEX, Cy3®, TET, Cy5®, AP5, YY... 3': TAMRA, DABCYL, BHQ™, DDQ, MGB-EDQ,...

Purifications: RP-HPLC or Dual HPLC

Quality Control: MALDI-TOF MS and analytical HPLC

Format: Dried

Packaging: 2 mL tube

Documentation: Technical data sheet

Probe Design: Available on request

Shipping: At room temperature



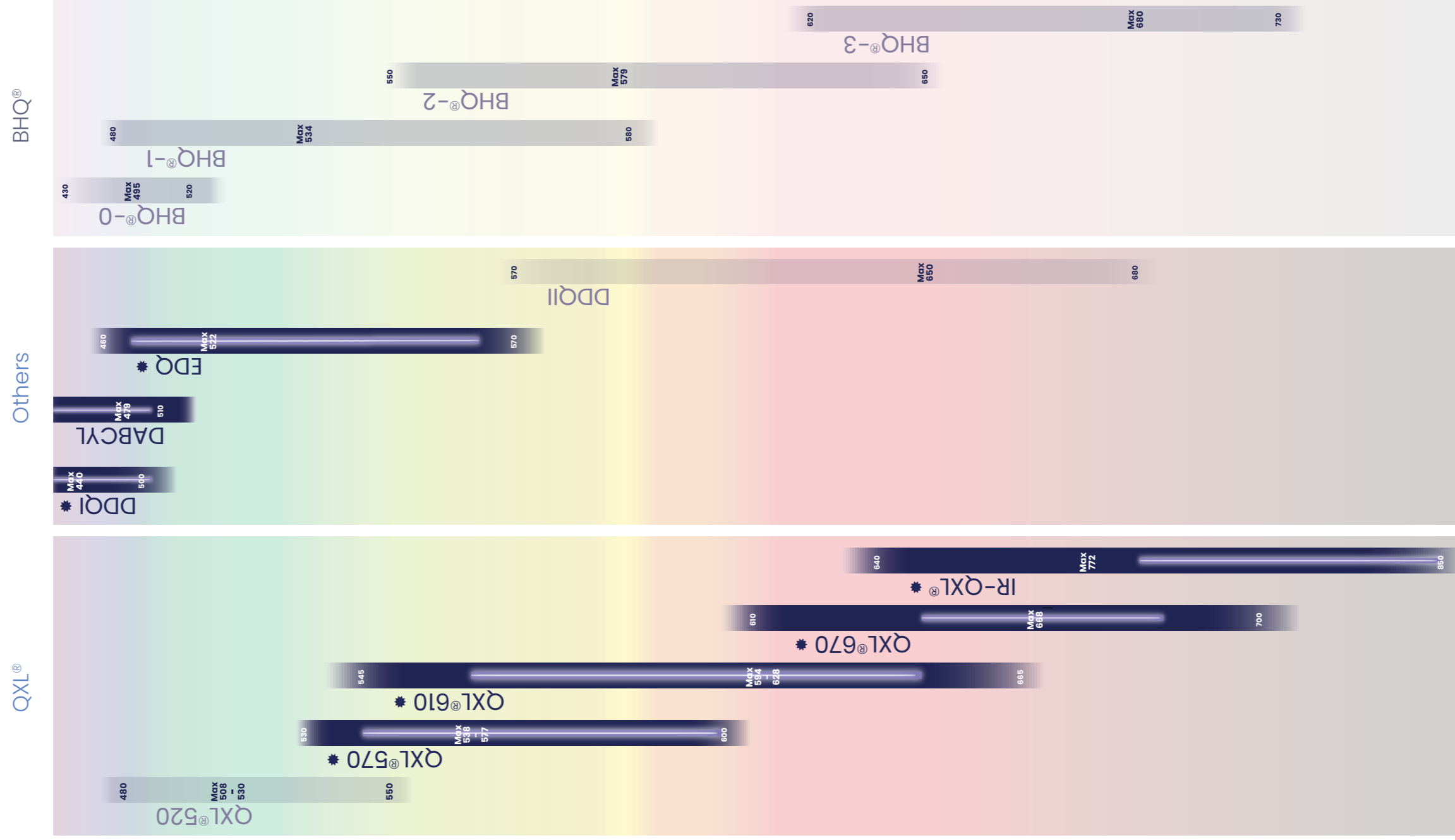
* Double-Dye, Molecular Beacon or MGB Probe. Each primer must be 15-30 DNA bases, unmodified and RP-Cartridge purified. The synthesis scale of the primers must be in accordance with the probe quantity (refer to www.eurogentec.com for details). Offer Valid only for orders placed on the Eurogentec e-commerce platform EOS. General conditions of sale will be applied. ** Restriction of use in the following countries: FR, UK, DE, IT, JP, ES. In these countries MGB probes must only be used for patient management. Use is free of limitation in other countries. End users are covered under Eurogentec's conveyed license for patient management.

FLUORESCENT DYES

Emission color	UV	Fluorescent Dyes	Max Abs in nm	Max Em in nm
550	UV	Crescendo blue	396	410
		Alexa Fluor® 405	401	421
		AMCA-X	353	442
		Alexa Fluor® 350	346	442
		Pacific Blue™	416	451
		Marina blue™	362	459
		Epech Blue	381	460
		ATTO 390	390	478
		ATTO 390	390	478
		ATTO 390	390	478
600	UV	ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
		ATTO 488	488	500
650	UV	Crescendo green	500	527
		Alexa Fluor® 500	500	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
		Rhodamine Green™	505	527
700	UV	Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
		Alexa Fluor® 514	514	530
750	UV	Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
		Alexa Fluor® 546	546	566
800	UV	Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589
		Alexa Fluor® 568	568	589

QUENCHERS

Research & Commercial use



Order

Oligonucleotides can be ordered via the web oligo configurators or on request. For more information, please contact us or visit our website.
www.eurogentec.com

High performance

Access™ Dyes & Quenchers from Eurogentec are a class of high performance proprietary molecules with IP-friendly terms and conditions.

Usage

In bold blue: Dyes & Quenchers for commercial & research use
In black: Dyes & Quenchers for research use only (RUO)
In grey: Not available at Eurogentec

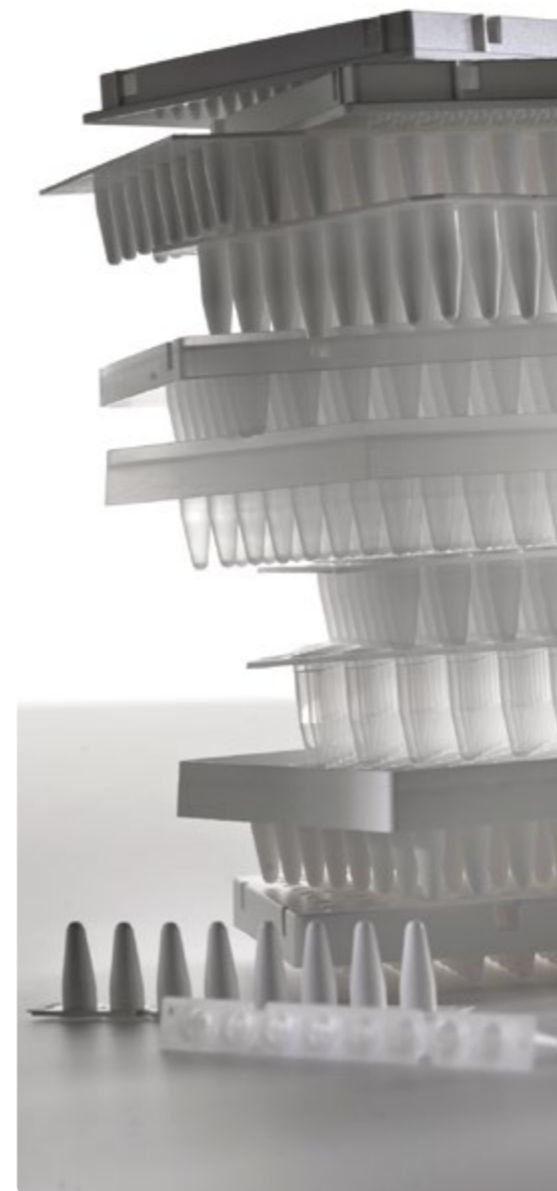
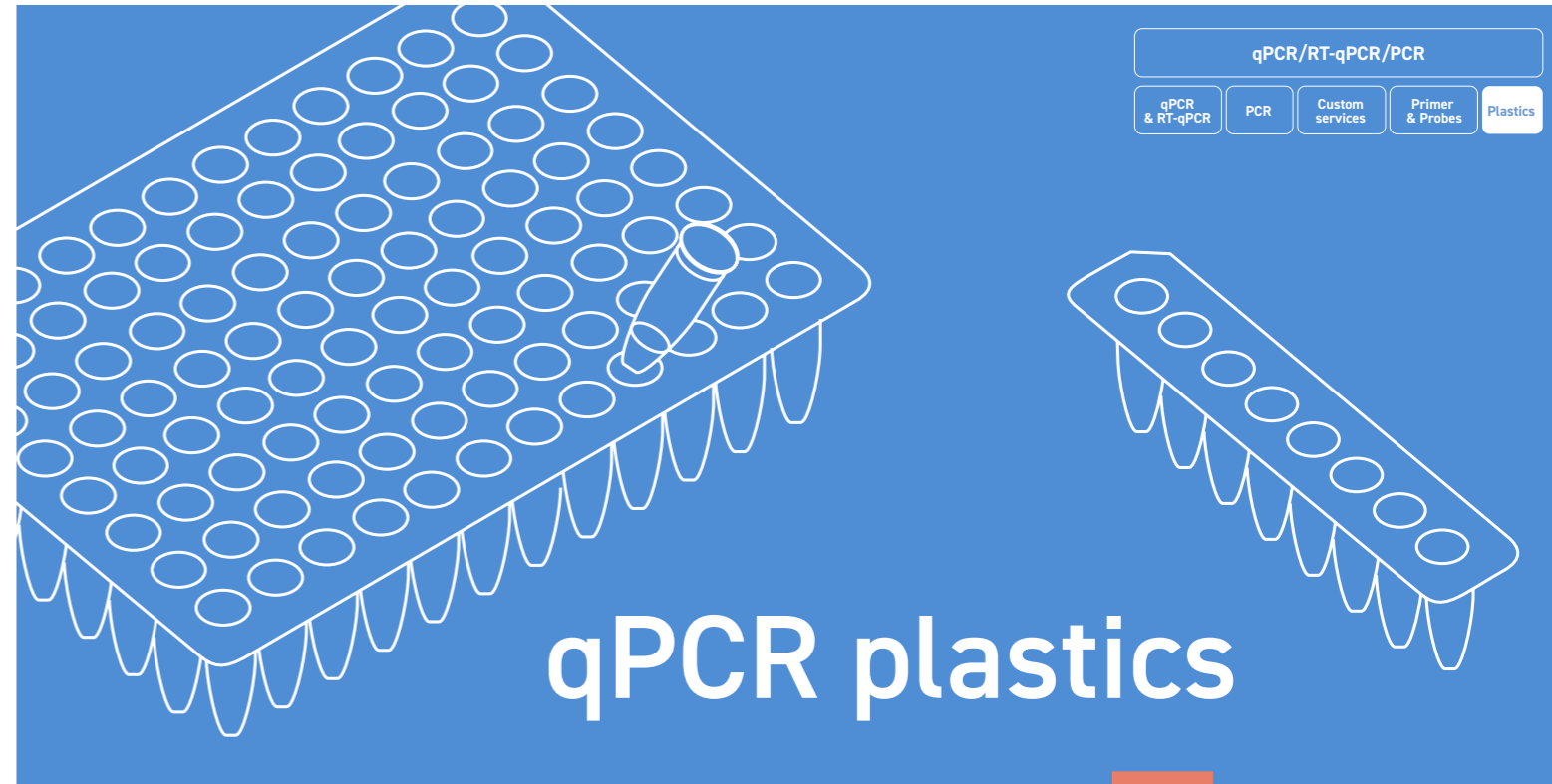
- Dark blue bands represent the indicated quenching range of recommended quenchers
- Grey bands represent the indicated quenching range of alternate quenchers.
- Purple bars represent the optimal recommended quenching range.
- Access™ Dyes & Quenchers

All wavelength values are in nm.
Note: Spectral properties of a dye or a quencher may vary when coupled to another molecule

* Perform a dye calibration for optimal results
For complementary information, please refer to instrument manufacturer technical guide or contact us at scientific.support@eurogentec.com
YV = Yokima Yellow™
DFO = Dragonfly Orange™
TR = Texas Red™
In blue = Recommended by Eurogentec
In black = Available at Eurogentec
In grey = Not available at Eurogentec

Thermocycler	Blue channel	Channel 1	Channel 2	Channel 3	Channel 4	Channel 5	Channel 6	Channel 7
GeneAmp SDS 9700	-	FAM	AP5/Y/Y/JOE/IC/TET*	DFO*/TAMRA/RED	-	-	-	-
ABI Prisma SDS 7000	-	FAM	AP5/Y/Y/JOE/IC/TET*	DFO*/TAMRA/RED	ROX	-	-	-
ABI Prisma SDS 7700	-	FAM	AP5/Y/Y/JOE/IC/TET*	DFO*/TAMRA/RED	ROX	-	-	-
ABI Prisma SDS 7900 HT	-	FAM	AP5/Y/Y/JOE/IC/TET*	DFO*/TAMRA/RED	ROX	-	-	-
ABI Prisma SDS 7300	-	FAM	AP5/Y/Y/JOE/IC/TET*	DFO*/TAMRA/RED	ROX	-	-	-
ABI Prisma SDS 7500	-	FAM	AP5/Y/Y/JOE/IC/TET*	DFO*/TAMRA/Cy3/RED	ROX/TR	Cy5/HL647	-	-
QuantStudio 3.5, 6.7	-	FAM	AP5/Y/Y/HEX*/JOE/IC/TET*	DFO*/TAMRA/Cy3/RED	ROX/TR	Cy5/HL647	-	-
QuantStudio 7	-	FAM	AP5/Y/Y/HEX*/JOE/IC/TET*	DFO*/TAMRA/Cy3/RED	ROX/TR	Cy5/HL647	Cy5.5/ATTO 700	-
Vii7	-	FAM	AP5/Y/Y/HEX*/JOE/IC/TET	DFO*/TAMRA/RED	ROX/TR*	ATTO 633/LIZ	ATTO 680/Alexa Fluor 680	-
StepOne	-	FAM	AP5/Y/Y/HEX*/JOE/IC/TET*	ROX	-	-	-	-
StepOnePlus	-	FAM	AP5/Y/Y/HEX*/JOE/IC/TET*	DFO*/TAMRA/Cy3/RED	ROX	-	-	-
Cycler iQ	-	FAM	AP5/Y/Y/HEX/Cy3/TET	DFO*/TAMRA/Cy3/RED	ROX/TR	Cy5	-	-
My iQ	-	FAM	-	-	-	-	-	-
iQ5	-	FAM	AP5/Y/Y/HEX/JOE/TET	DFO*/TAMRA/Cy3/RED	ROX/TR	Cy5	-	-
CFX 96	-	FAM	AP5/Y/Y/HEX/JOE/TET	ROX/TR	Cy5	Cy5.5	-	-
MiniAptari	-	FAM	AP5/Y/Y/HEX/TET*	-	-	-	-	-
DNA Engine Opticon 1	-	FAM	-	-	-	-	-	-
DNA Engine Opticon 2	-	FAM	AP5/Y/Y/HEX/TAMRA/IC/TET	-	-	-	-	-
Chromo 4	-	FAM	AP5/Y/Y/HEX/TAMRA/Cy3/JOE/IC/TET	ROX/TR	Cy5	-	-	-
Mx3005P (choice of 4 filters)	-	FAM	TET	YY*/HEX/JOE/IC	Cy3	DFO/TAMRA/RED	TR/ROX	Cy5
Mx3005P (choice of 5 filters)	-	FAM	TET	YY*/HEX/JOE/IC	Cy3	DFO/TAMRA/RED	TR/ROX	Cy5
Mx4000 (choice of 4 filters)	-	FAM	TET	YY*/HEX/JOE/IC	Cy3	DFO/TAMRA/RED	TR/ROX	Cy5
AriaMx (choice of up to 6 filters)	-	FAM	HEX	Cy3	ROX	Cy5	-	-
Mastecycler ep resiplex2	-	FAM	AP5/Y/Y/HEX/JOE/IC/TET	-	-	-	-	-
Mastecycler ep resiplex4	-	FAM	AP5/Y/Y/HEX/T/JOE/IC	DFO*/TAMRA	ROX	-	-	-
LightCycler 15	-	FAM	AP5/Y/Y/HEX/JOE/IC/TET	Cy5	-	-	-	-
LightCycler 2	-	FAM	AP5/HEX/YY*/JOE*/IC	TR/LC - Red 610	ATTO 620/LC Red 640	Cy5/LC Red 670	ATTO 680 / LC Red 705	-
LightCycler 480 I	-	FAM	AP5/Y/Y/HEX/JOE/IC	ROX/TR/LC Red 610	ATTO 620/LC Red 640	Cy5	-	-
LightCycler 480 II	-	FAM	AP5/Y/Y/HEX/JOE/IC	ROX/TR/LC Red 610	ATTO 620/LC Red 640	Cy5/Cy 5.5	-	-
LightCycler 96	-	FAM	AP5/Y/Y/HEX/IC	TR/LC - Red 610	Cy5	-	-	-
SmartCycler1	-	FAM	AP5/Y/Y/Cy3/JOE/IC/TET	TR	Cy5	-	-	-
SmartCycler2	-	FAM	AP5/Cy3/YY*/JOE*/TET	TR/ROX	Cy5	-	-	-
Rotor-Gene 2000 / 3000	-	FAM	AP5/Y/Y/JOE/IC/TET	TAMRA/ROX/Cy3.5/TR	Cy5	-	-	-
Rotor-Gene 6000	-	FAM	AP5/Y/Y/HEX/JOE/IC/TET	TAMRA/ROX/Cy3.5/TR	Cy5	ATT 0880	-	-
PhicoReal	-	FAM	AP5/HEX/YY	ROX/TR	Cy5	-	-	-
Clower	-	FAM	AP5/Y/Y/HEX/JOE/IC	DFO/TAMRA/RED	ROX/TR/Cy3.5	Cy5	Cy5.5	-

* perform a dye calibration for optimal results YY = Yakima Yellow® DFO = Dragonfly Orange™ TR = Texas Red® DFO = Dragonfly Orange™ TR = Texas Red® In grey = Recommended by Eurogentec In blue = Available at Eurogentec



Our virgin optical quality polypropylene blend is optimised for qPCR and causes no background fluorescence. The wall thickness has been designed for an optimal heat transfer and to minimise evaporation of the reaction mix.

The antistatic properties of our plastics reduce reagents adsorption on the tube walls, increase the reaction yield and allow the use of small reaction volumes. The plates are perfectly flat, thin-walled and rigid, and show extreme uniformity to avoid edge artefacts.

Natural, white or frosted plates are available. White plates significantly increase signal-to-noise ratio compared to the natural plates but it may be difficult to see through the content. The best compromise remains the frosted plates which allow a direct view of the content and a high signal-to-noise ratio.

The super-tight closure and sealing properties reduce evaporation, leading to more reproducible results. The optical seals do not interfere with the signal, contrary to many other brands.

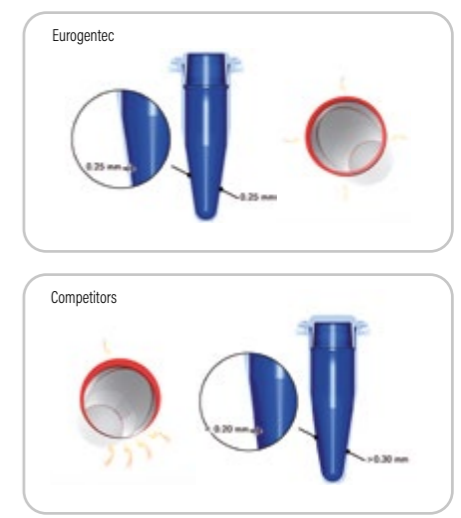
Please refer to the annexes to know which plastics is the most suited for your thermocycler. (pages 28 → 30) ■

FEATURES

- No background fluorescence to increase the signal to noise ratio
- Optimal heat transfer to increase the reaction yield
- Extreme uniformity to avoid edge artefacts
- High purity, DNA(se), RNA(se), pyrogen-free to avoid sample contamination and degradation

BENEFITS

- More accurate and reproducible results
- Increased signal intensity



Eurogentec's wells show an extreme uniformity of wall thickness avoiding edge artefacts. Moreover, the wall is thin at the bottom for an optimal heat transfer and thicker at the top to minimise evaporation. Red circles represent a top view of the wells.

96-well plates

Our well plates are available for all thermocyclers and can be combined with either optical flat 8-caps strips as well as with qPCR optical seals. All non-skirted plates can be easily cut into 16, 24, 32 or 48-well pieces.

qPCR 96-well plates for ABI GeneAmp® 5700 SDS, ABI Prism® SDS 7000, 7300, 7500 Regular block, 7700, 7900 Regular block and ViiA™7 Regular block, Quanstudio regular (0,2 ml block)

- High profile
- Thin-walled
- Sub-skirted
- Optical quality
- Frosted or white



qPCR 96-well plates for iCycler iQ®, iQ5, MyiQ, Mx4000®, Mx3000P®, Mx3005P® and MasterCycler® ep realplex

- Also compatible with most regular 96-well ABI Real-Time thermocyclers
- High profile
- Thin-walled
- Non-skirted
- Natural or white
- Easy to cut into 16, 24, 32, 48-well pieces



Description	Reference	Quantity
qPCR 96-well plate sub-skirted, high profile, frosted (ABI systems)	RT-PL96-AB	5 x 5 plates
qPCR 96-well plate sub-skirted, high profile, frosted (ABI systems), 200 pack	RT-PL96-AB-P200	200 plates + seals
qPCR 96-well plate sub-skirted, high profile, frosted (ABI systems), 500 pack	RT-PL96-AB-P500	500 plates + seals
qPCR 96-well plate sub-skirted, high profile, white (ABI systems)	RT-PL96-ABW	5 x 5 plates
qPCR 96-well plate sub-skirted, high profile, white (ABI systems), 200 pack	RT-PL96-ABW-P200	200 plates + seals
qPCR 96-well plate sub-skirted, high profile, white (ABI systems), 500 pack	RT-PL96-ABW-P500	500 plates + seals

Description	Reference	Quantity
qPCR 96-well plate non-skirted, high profile, natural	RT-PL96-MQ	5 x 5 plates
qPCR 96-well plate non-skirted, high profile, natural, 200 pack	RT-PL96-MQ-P200	200 plates + seals
qPCR 96-well plate non-skirted, high profile, natural, 500 pack	RT-PL96-MQ-P500	500 plates + seals
qPCR 96-well plate non-skirted, high profile, white	RT-PL96-MQW	5 x 5 plates
qPCR 96-well plate non-skirted, high profile, white, 200 pack	RT-PL96-MQW-P200	200 plates + seals
qPCR 96-well plate non-skirted, high profile, white, 500 pack	RT-PL96-MQW-P500	500 plates + seals

qPCR 96-well plates for ABI Prism® SDS 7500 Fast block, 7900 Fast block, StepOne Plus and ViiA™7 Fast block, Quanstudio regular (0,1 ml block)

- Low profile
- Thin-walled
- Sub-skirted
- Optical quality
- Frosted or white



qPCR 96-well plates for DNA Engine Opticon® 1 and 2, MiniOpticon, Chromo4, StepOne and Quantica®, Biorad CFX 96

- Low profile
- Thin-walled
- Non-skirted
- Frosted or white
- Easy to cut into 16, 24, 32, 48-well pieces



Description	Reference	Quantity
qPCR 96-well plate sub-skirted, low profile, frosted (ABI FAST systems)	RT-PL96-AF	5 x 5 plates
qPCR 96-well plate sub-skirted, low profile, frosted (ABI FAST systems), 200 pack	RT-PL96-AF-P200	200 plates + seals
qPCR 96-well plate sub-skirted, low profile, frosted (ABI FAST systems), 500 pack	RT-PL96-AF-P500	500 plates + seals
qPCR 96-well plate sub-skirted, low profile, white (ABI FAST systems)	RT-PL96-AFW	5 x 5 plates
qPCR 96-well plate sub-skirted, low profile, white (ABI FAST systems), 200 pack	RT-PL96-AFW-P200	200 plates + seals
qPCR 96-well plate sub-skirted, low profile, white (ABI FAST systems), 500 pack	RT-PL96-AFW-P500	500 plates + seals

Description	Reference	Quantity
qPCR 96-well plate non-skirted, low profile, frosted	RT-PL96-OP	5 x 5 plates
qPCR 96-well plate non-skirted, low profile, frosted, 200 pack	RT-PL96-OP-P200	200 plates + seals
qPCR 96-well plate non-skirted, low profile, frosted, 500 pack	RT-PL96-OP-P500	500 plates + seals
qPCR 96-well plate non-skirted, low profile, white	RT-PL96-OPW	5 x 5 plates
qPCR 96-well plate non-skirted, low profile, white, 200 pack	RT-PL96-OPW-P200	200 plates + seals
qPCR 96-well plate non-skirted, low profile, white, 500 pack	RT-PL96-OPW-P500	500 plates + seals

qPCR 96-well plates for LC480

- Developed for LC480, no adaptor needed.
- Low profile
- Thin-walled
- Semi-skirted
- White



It is necessary to use our qPCR optical caps (flat 8-cap strips) or seals, to seal our plates

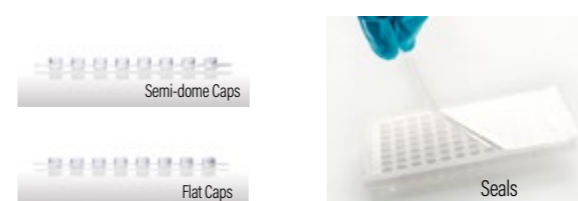
Description	Reference	Quantity
qPCR 96-well plate for LC480, low profile, thin-walled, semi-skirted, white	RT-PL96-LC480	5 x 5 plates
qPCR 96-well plate for LC480, low profile, thin-walled, semi-skirted, white, 200 pack	RT-PL96-LC480-P200	200 plates + seals
qPCR 96-well plate for LC480, low profile, thin-walled, semi-skirted, white, 500 pack	RT-PL96-LC480-P500	500 plates + seals

Discontinued: the previous model RT-PL96-LC which must be combined with the adaptor has been discontinued.

qPCR optical caps & seals

The extremely clear glass-grade area of the cap ensures very good transmission of the fluorescence and equals the performance properties of optical seals. When combined with our qPCR plates they seal better than other brands, thus less evaporation.

Our adhesive optical seals have excellent optical quality properties and will not interfere with the signal, leading to higher signal-to-noise ratio compared to other brands. The glue that we use has been specifically developed for the type of plastics that we use, thus it sticks to the plate very well and leads to better sealing and less evaporation. The seal itself is thicker than usual films, which avoids potential tears; moreover it will not curl and stick to itself.



Description	Reference	Quantity
qPCR optical semi-dome 8-cap strip, natural	RT-OCAP-300	300 strips
qPCR optical "wide indented" flat 8-cap strip, natural	RT-FLAT-120, RT-FLAT-300	120 strips or 300 strips
qPCR opti-seals (optical, disposable, adhesive)	RT-OPSL-25, RT-OPSL-100	25 or 100

qPCR 8-tube strips



These tube-strips are designed for PCR and qPCR applications. For the highest S/N ratios in qPCR the use of white or frosted products is recommended. qPCR 8-tube strips are compatible with almost all PCR and qPCR cyclers which accept regular or low profile products. Please refer to the compatibility table to select the 8-tube strip adapted to your thermocycler (p.28).

Description	Reference	Quantity
0.1 mL Low Profile		
qPCR 8-tube strip, frosted	RT-STR8-LF120	1 bag of 120 strips
qPCR 8-tube strip, frosted + flat caps	RT-STR8-LF-P600	600x strips + 600 caps
qPCR 8-tube strip, white	RT-STR8-LW120	1 bag of 120 strips
qPCR 8-tube strip, white + flat caps	RT-STR8-LW-P600	600x strips + 600 caps
0.2 mL Regular Profile		
qPCR 8-tube strip, frosted	RT-STR8-RF120	1 bag of 120 strips
qPCR 8-tube strip, frosted + flat caps	RT-STR8-RF-P600	600x strips + 600 caps
qPCR 8-tube strip, white	RT-STR8-RW120	1 bag of 120 strips
qPCR 8-tube strip, white+ flat caps	RT-STR8-RW-P600	600x strips + 600 caps
LC480 96-well plate adaptor*	RT-PL96-LCAD	1 piece

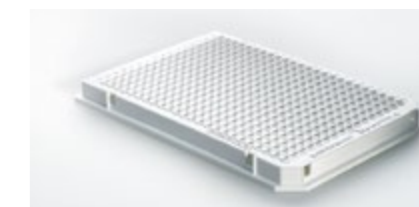
*This adaptor is mandatory for use in LC480

384-well plates

For Roche, ABI, Bio-Rad & other systems

Our skirted 384-well plate types perfectly fit to Roche, ABI, Bio-Rad and Eppendorf qPCR platforms, as well as many regular PCR cyclers. The plates are fully optimised for robotic applications used for high-throughput, low volume processing. The plates are perfectly suited for diagnostic & clinical procedures. Each plate is lasermarked with a unique code.

- Perfect fit to Roche, ABI, Bio Rad and Eppendorf systems
- Fully skirted, low profile, white plate
- Low evaporation, very robust, due to clever design
- Extreme Uniform wells lead to reproducible results
- Robotic friendly, easily stackable and flat
- High S/N Ratio



Description	Reference	Quantity
For Roche LC480	RT-PL384-LCW	8 x 5 plates
For ABI, Bio-Rad & Eppendorf systems	RT-PL384-ABW	8 x 5 plates

Plastics Samples

Reference	Product name	Quantity
qPCR 96-well plate type		
RT-PL96-ABSA	qPCR 96-well plate sub-skirted, high profile, frosted (ABI systems) sample	2
RT-PL96-ABWSA	qPCR 96-well plate sub-skirted, high profile, white (ABI systems) sample	2
RT-PL96-AFSA	qPCR 96-well plate sub-skirted, low profile, frosted (ABI FAST systems) sample	2
RT-PL96-AFWSA	qPCR 96-well plate sub-skirted, low profile, white (ABI FAST systems) sample	2
RT-PL96-MQSA	qPCR 96-well plate non-skirted, high profile, natural sample	2
RT-PL96-MQWSA	qPCR 96-well plate non-skirted, high profile, white sample	2
RT-PL96-OPSA	qPCR 96-well plate non-skirted, low profile, frosted sample	2
RT-PL96-OPWSA	qPCR 96-well plate non-skirted, low profile, white sample	2
RT-PL96-LC480SA	qPCR 96-well plate semi-skirted, low profile, white + seals for LC480 sample	2
qPCR 384-well plate type		
RT-PL384-LCWSA	qPCR 384-well white plate, skirted, low profile for LC480 sample	2
RT-PL384-ABWSA	qPCR 384-well white plate, skirted, low profile sample	2
qPCR 8-tube Strips		
RT-STR8-RFSA	qPCR 8-tube strip, regular profile - 0.2 mL -frosted sample	6
RT-STR8-RWSA	qPCR 8-tube strip, regular profile - 0.2 mL -white sample	6
RT-STR8-LFSA	qPCR 8-tube strip, low profile - 0.1 mL -frosted sample	6
RT-STR8-LWSA	qPCR 8-tube strip, low profile - 0.1 mL -white sample	6
Covers		
RT-FLAT-SA	qPCR optical flat 8-cap strip, natural sample	6
RT-OCAP-SA	qPCR optical semi-domed 8-cap strip, natural sample	24
RT-OPSL-SA	qPCR optical seals (disposable, adhesive) sample	2

96-WELL PLATE COMPATIBILITY WITH REAL-TIME THERMOCYCLERS

qPCR 96-well plate type	Cover	Cycler Models																								
		ABI GeneAmp 5700	ABI Prism SDS 7000-7300-7500-7700	ABI Prism SDS 7500 FAST	ABI Prism SDS 7900HT	ABI Prism SDS 7900HT FAST	ABI Step One	ABI Step One Plus	ABI ViiA7 qPCR system	ABI ViiA7 FAST	ABI QuantStudio 6-7-12K Flex	ABI QuantStudio 6-7-12K Flex FAST	Bio-Rad MyiQ-iQ-iQ5	Analytik Jena/Biometra TOptical-flower	Agilent/Stratagene Mx3000P-Mx3005P-Mx4000	Eppendorf Mastercycler ep realplex Silver	Bio-Rad DNA Engine (Dyad-Tetrad 2) Chromo4	Bio-Rad DNA Engine Opticon 2	MiniOpticon	Bio-Rad CFX 96	Bio-Rad CFX 384	Roche LightCycler 480	Roche COBAS z 480	Roche LightCycler 96	Thermo Fisher/Finzymes PikoReal 24-well	
Sub-skirted, high profile, frosted or white type (ABI systems) RT-PL96-AB or RT-PL96-ABW	Flat caps RT-FLAT-120	V	V	V			V	V	C	V	C	C														
Sub-skirted, low profile, frosted or white type (ABI systems) RT-PL96-AF or RT-PL96-AFW	Adhesive seals RT-OPSL-100	V	V	V			V	V	C	V	C	C														
Non-skirted, high profile, natural or white type RT-PL96-MQ or RT-PL96-MQW	Flat caps RT-FLAT-120	C	C				C	C	V	V	V	V						X								
Non-skirted, low profile, frosted or white type RT-PL96-OP or RT-PL96-OPW	Adhesive seals RT-OPSL-100	C	C				C	C	V	V	V	V						X								
Semi-skirted, low profile, white RT-PL96-LC480	Flat caps RT-FLAT-120			C	C	X	C				C					V	V						V	V	V	
Semi-skirted, low profile, white RT-PL96-LC480	Adhesive seals RT-OPSL-100			C	C	X	C				C					V	V						V	V	V	
Skirted, low profile, white plate for ABI RT-PL384-ABW	Adhesive seals RT-OPSL-100			V	V		V	V														V				
Skirted, low profile, white plate for LC RT-PL384-LCW	Adhesive seals RT-OPSL-100																				V	V				
qPCR 8-tube Strips	Cover	Refers to cycler manufacturer instructions for required adaptor																								
Extra Robust, Regular Profile, Frosted, 0.2 ml RT-STR8-RF120	Flat caps RT-FLAT-120	C	C	C			C	C	C	C	C	C						C								
Extra Robust, Regular Profile, White, 0.2 ml RT-STR8-RW120	Flat caps RT-FLAT-120	C	C	C			C	C	C	C	C	C						C								
Extra Robust, Low Profile, Frosted, 0.1 ml RT-STR8-LF120	Flat caps RT-FLAT-120			C	C	C	C	C								C	C						C**	C	V	
Extra Robust, Low Profile, White, 0.1 ml RT-STR8-LW120	Flat caps RT-FLAT-120			C	C	C	C	C								C	C						C**	C**	C	V

V for best recommended plastics.
 C for compatible plastics. qPCR optical semi-dome 8 cap strips are also available (RT-OCAP-300: 300 strips). They can be used on several machines.
 X plate must be cut to size (48 well) to fit Step One or MiniOpticon. ** adaptor needed (Ref: RT-PL96-LCAD).
 Please refer to cycler manufacturer instruction for required adaptor or contact us at scientific.support@eurogentec.com

PCR PLATE COMPATIBILITY TABLE (PART 1/2)

Manufacturer	Model	# Wells	Sub-skirted, high profile, frosted	Sub-skirted, low profile, frosted	Non-skirted, high profile, natural	Non-skirted, low profile, frosted	Skirted, low profile, white
			RT-PL96-AB	RT-PL96-AF	RT-PL96-MQ	RT-PL96-OP	RT-PL384-LCW
ABI	GeneAmp PCR System 2400	24	x	x	X	x	x
ABI	GeneAmp PCR System 9600	96	V	x	C	x	x
ABI	2700 Thermal Cycler	96	V	x	C	x	x
ABI	2720 Thermal Cycler	96	V	x	C	x	x
ABI	GeneAmp PCR System 9700	96/384	C	x	C	x	C
ABI	9800 Fast Thermal Cycler	96 fast	x	V	x	C	x
ABI	Verity	96/384	C	x	C	x	C
Bioer	XP Cycler	30/48/96/384	x	x	C	C	x
Bioer	Little Genius	25	x	x	X	X	x
Bioer	Life Express	96	x	x	C	C	x
Biometra	Uno	24/48/96	x	x	V	C	x
Biometra	Uno II	24/48/96/384	x	x	V	C	x
Biometra	Tpersonal	20/48	x	x	X	X	x
Biometra	T1 (plus)	48/96/384	x	x	V	C	C
Biometra	TGradient	48/96/384	x	x	V	C	x
Biometra	T3/T3000	20/48	x	x	X	X	x
Biometra	TProfessional	96/384	x	x	V	C	C
Biometra	TRobot	96/384	x	x	V	C	C
Bio-Rad	Gene Cycler	24	x	x	X	x	x
Bio-Rad	iCycler	48/96/384	x	x	C	x	C
Bio-Rad	MyCycler	96	x	x	C	x	x
Bio-Rad	C1000	96/384	x	x	C	x	C
CLP	ATC 201	96	x	x	C	C	x
CLP	ATC 401	96	x	x	C	C	x
Corbett Research	Palm-Cycler	60/96/384	x	x	C	C	C
Eppendorf	MasterCycler	96/384	x	x	C	C	C
Eppendorf	MasterCycler gradient	96/384	x	x	C	C	C
Eppendorf	MasterCycler ep gradient	96/384	x	x	C	C	C
Eppendorf	MasterCycler ep gradient S	96/384	x	x	C	C	C
GRI	G-storm GSI	48 0.2ml/0.5ml	x	x	C	C	x
GRI	G-storm GSI	96/384	x	x	X	X	x
GRI	G-storm GSX	48 0.2ml/0.5ml	x	x	C	C	x
GRI	G-storm GSX	96/384	x	x	X	X	x
GRI	G-storm GS4	48 0.2ml/0.5ml	x	x	C	C	x
GRI	G-storm GS4	96/384	x	x	X	X	x
MJ Research/Bio-Rad	Mini	48	x	x	X	X	x
MJ Research/Bio-Rad	PTC-100	60/96	x	x	C	C	x
MJ Research/Bio-Rad	PTC-200	30/48/96/384	x	x	C	C	C

C Fits this Cycler
 V Most optimal plastic for this thermocycler
 X The plate must be cut
 x Not compatible

PCR PLATE COMPATIBILITY TABLE (PART 2/2)

Manufacturer	Model	# Wells	Sub-skirted, high profile, frosted	Sub-skirted, low profile, frosted	Non-skirted, high profile, natural	Non-skirted, low profile, frosted	Skirted, low profile, white
			RT-PL96-AB	RT-PL96-AF	RT-PL96-MQ	RT-PL96-OP	RT-PL384-LCW
MJ Research/Bio-Rad	PTC-221	30/48/96/384	x	x	C	C	C
MJ Research/Bio-Rad	PTC-220	30/48/96/384	x	x	C	C	C
MJ Research/Bio-Rad	PTC-225	30/48/96/384	x	x	C	C	C
MWG	Primus 96	96	x	x	C	C	x
MWG	Primus HT Dualblock	96/384	x	x	C	C	C
MWG	Primus Multiblock System	96/384	x	x	C	C	C
Stratagene	Robo-Cycler	96	x	x	C	C	x
Takara	TP 600	96	x	x	C	C	x
Takara	TP 650	96	x	x	C	C	x
Takara	TP 3000	96	x	x	C	C	x
Techne	Genius	40/96/384	x	x	C	C	C
Techne	TouchGene	40/96	x	x	C	C	x
Techne	FlexiGene/FlexiQuad	96/384	x	x	C	C	C
Techne	TechGene	25/96	x	x	C	C	x
Techne	Gene-E	96	x	x	C	C	x
Techne	TC-412	96	x	x	C	C	x
Techne	TC-512	96/384	x	x	C	C	C
ThermoHybaid	PxE	96	x	x	C	C	x
ThermoHybaid	OmniGene	96	x	x	C	C	x
ThermoHybaid	PCR Express	48/96/384	x	x	C	C	C
ThermoHybaid	Omn-E	96	x	x	C	C	x
ThermoHybaid	TouchDown	48/96	x	x	C	C	x
ThermoHybaid	Px2	96/384	x	x	C	C	C
ThermoHybaid	PCR sprint	24	x	x	⊂	⊂	x
ThermoHybaid	MBS	96/384	x	x	C	C	C
ThermoHybaid	MBSR	96/384	x	x	C	C	C

SEQUENCER COMPATIBILITY

Manufacturer	Model	Sub-skirted, high profile, frosted	Sub-skirted, low profile, frosted	Non-skirted, high profile, natural	Non-skirted, low profile, frosted
		RT-PL96-AB	RT-PL96-AF	RT-PL96-MQ	RT-PL96-OP
ABI	310	x	x	C	x
ABI	3100	x	x	C	x
ABI	3130	x	x	C	x
ABI	3700	C	x	C	x
ABI	3730	C	x	C	x
Beckman	CEQ 8000	C	x	x	x
Beckman	CEQ 8800	C	x	x	x
Transgenomics	Wave MD System 4000	x	x	C	x

Models not mentioned within this list are not compatible with our plates.

Any type of cap/seal can be combined with these machines. Semi-domed caps are recommended for ABI machines.

C Fits this Cycler

V Most optimal plastic for this thermocycler

⊂ The plate must be cut

x Not compatible

ABI Prism® is a registered trademarks of The Perkin-Elmer Corp. • Chromo4® is a registered trademark of Bio-Rad Laboratories, Inc. • GeneAmp® is a registered trademark of Roche Molecular Systems, Inc. • iCycler iQ® is a registered trademark of Bio-Rad Laboratories, Inc. • LightCycler® is a registered trademark of Roche Diagnostics GmbH. • MasterCycler® is a registered trademark of Eppendorf, Inc. • Mx3000®, Mx3000P®, Mx4000® are registered trademarks of Stratagene • Opticon® is a registered trademark of Bio-Rad Laboratories, Inc. • QuantStudio® is a trademark of Life Technologies. • Quantica® is a registered trademark of Techne Ltd and Techne, Inc. • StepOne® and StepOne Plus® are trademarks of Applied Biosystems. • ViiA™7 is a trademark of Life Technologies.

HOW TO ORDER

ONLINE

WWW.EUROGENTEC.COM

FAST AND SIMPLE

1. Register to or log in www.eurogentec.com
2. Select your product(s)
3. Fill in your cart
4. Finalise your order

FULL OF ADVANTAGES

1. Manage your profile
2. Fast order and favourite products
3. Request quotes

Discounted prices are automatically applied and tagged with a four leaves clover on the related product web pages.

CUSTOM PRODUCTS

Contact us at info@eurogentec.com to discuss your project.

RELATED PRODUCTS

Oligonucleotides

- Custom and catalogue oligonucleotides and probes
- RNAi oligonucleotides

Smart gene Analysis

- Nucleic acids purification and extraction kits
- Electrophoresis Systems and Reagents

TRADEMARKS/LABELS

ABI Prism® 5700, 7000, 7500, 7700 and 7900 are registered trademarks of Applied Biosystems, Inc.

Alexa Fluor® is a registered trademark of Life Technologies Corp.

Black Hole Quencher® is a registered trademark of Biosearch Technologies, Inc.

Cy® and CyDye® are registered trademarks of Cytiva.

Diamond Taq® is a registered trademark of Kaneka Eurogentec S.A.

DNA Engine Opticon® is a registered trademark of Bio-Rad Laboratories, Inc.

Dragonfly Orange™ is a trademark of Kaneka Eurogentec S.A.

Eclipse™ Quencher and Dark Quencher are a trademark of ElitechGroup, Inc.

GeneAmp® is a registered trademark of Roche Molecular Systems, Inc.

GoldStar® is a registered trademark of Kaneka Eurogentec S.A.

HiLyte™ is a trademark of Anaspec, Inc.

iCycler iQ® is a registered trademark of Bio-Rad Laboratories, Inc.

iQ™5 is a trademark of Bio-Rad Laboratories, Inc.

LightCycler® is a registered trademark of Roche Diagnostics GmbH.

LNA® is a registered trademark of Exiqon A/S.

MasterCycler® is a registered trademark of Eppendorf, Inc.

Mx3000®, Mx3000P®, Mx 3005P®, Mx4000® are registered trademarks of Stratagene.

Opticon® is a registered trademark of MJ Research, Inc.

Quantica® is a registered trademark of Techne Ltd and Techne, Inc.

QXL® is a registered trademark of Anaspec, Inc.

Rotor-Gene® is a registered trademark of Corbett Research Pty Ltd.

SmartCycler® is a registered trademark

of Cepheid, Inc.

StepOne™ is a registered trademark of Applied Biosystems.

SYBR® Green is a registered trademark of Life Technologies Corp.

Taqman® is a registered trademark of Roche Diagnostics GmbH.

Takyon® is a trademark of Kaneka Eurogentec S.A.

Texas Red® is a registered trademark of Life Technologies Corp.

Vic® is a registered trademark of Applied Biosystems.

Yakima Yellow® is a registered trademark of ElitechGroup, Inc.

BE AWARE OF

EUROGENTEC LICENSE STATEMENT

Sept_2022

1. Elitech Dyes
 - Use of this product (including but not limited to Yakima Yellow®) is covered by patents owned by ElitechGroup, Inc., including but not limited to the following US patents and corresponding patent claims outside the US: US 6,972,339, 7,541,454, 7,671,218, 7,767,834, 8,163,910. This product, including both Research Grade Product and IVD Grade Product, is sold under a licensing arrangement with ElitechGroup, Inc. for the purchaser's own internal research. Commercial use of this product requires a separate license from Kaneka Eurogentec S.A. Further information on purchasing licenses may be obtained by contacting licensing@eurogentec.com.
2. SYBR® Green I nucleic acid gel stain
 - This product is provided under a technology agreement from Life Technologies Corporation to Kaneka Eurogentec S.A. The sale of this product is conditioned on the buyer using the product only in research conducted by the buyer and not to sell or otherwise commercialise this product or its components (1) in manufacturing; (2) to provide a service, information, or data in exchange for payment; (3) for therapeutic, diagnostic or prophylactic purposes; or (4) for resale, not even if resold for use in research. For information on purchasing a license to this product for purposes other than research, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, CA 92008 USA or outlicensing@lifetech.com.
3. QXL® Quenchers
 - Use of this product is covered by one or more of the following US patents owned by AnaSpec, Inc.: USP 7,910,753, USP 8,093,411, and divisionals, continuations, continuations-in-part, reissues, substitutes, and extensions thereof. The purchase of this product conveys to the buyer a limited, non-exclusive, nontransferable right (without the right to sell, repackage, or further sublicense) under these patent rights to use only the amount of product purchased for the purchaser's own internal research. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of this product does not include nor carry any right or license to use, develop, or otherwise exploit this product commercially, and no rights are conveyed to the buyer to use the product or components of the product for any other purposes, including without limitation, provision of services to a third party, generation of commercial databases, or clinical diagnostics or therapeutics. This product is for research use only. For information on purchasing a license to this product for purposes other than research, contact licensing@eurogentec.com.
4. UDG
 - Use of UDG employs patent rights licensed from Invitrogen Corporation.
5. MGB Pleiades Probes
 - Use of this product is covered by patents owned by ElitechGroup, Inc., including but not limited to the following US patents and corresponding patent claims outside the US: US-2009-0048427-A1 and US 7,381,818. This product, including both Research Grade Product and IVD Grade Product, is sold under a licensing arrangement with ElitechGroup, Inc. for the purchaser's own internal research. Commercial use of this product requires a separate license from Kaneka Eurogentec S.A. Further information on purchasing licenses may be obtained by contacting licensing@eurogentec.com.

www.eurogentec.com

PCR/qPCR

Contact:

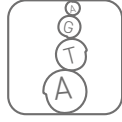
For more information,
please contact your Sales
representative or contact us at:
isales@eurogentec.com

Kaneka Eurogentec S.A. 2019

Follow us on



Life Science Research



Oligos



PCR/qPCR



Genes/DNA



Antibodies



Peptides



Proteins