



Takyon[®]
Amplification Technologies

qPCR • RT-qPCR • PCR

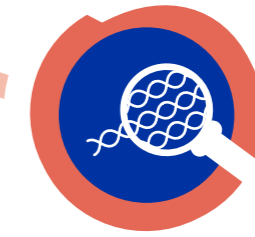
For Life Science Research

- > 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

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 *Positive controls*
- 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.

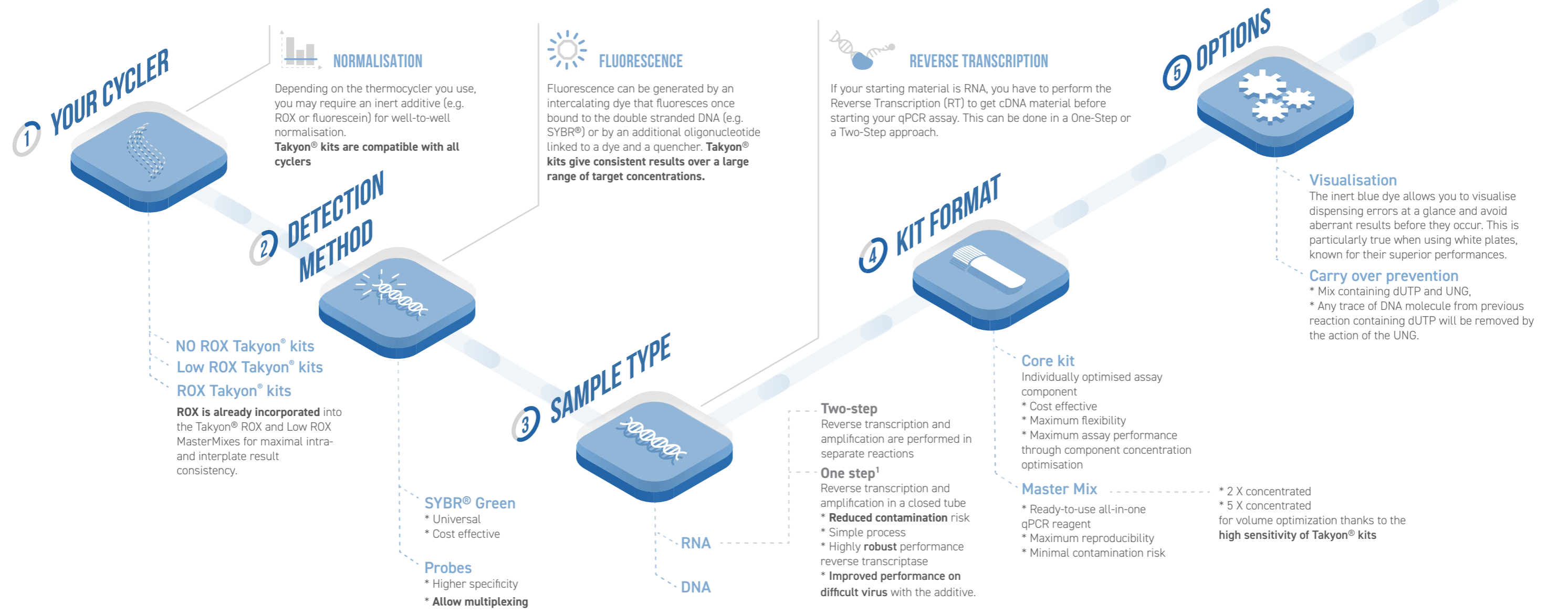
Small reaction volumes
The high sensitivity of the Takyon® kits combined to the highly concentrated 5X format give reliable results even in small volume reaction.

Suitable for multiplex
The Takyon® kits allow the detection of multiple targets in a single reaction tube.

GOOD TO KNOW

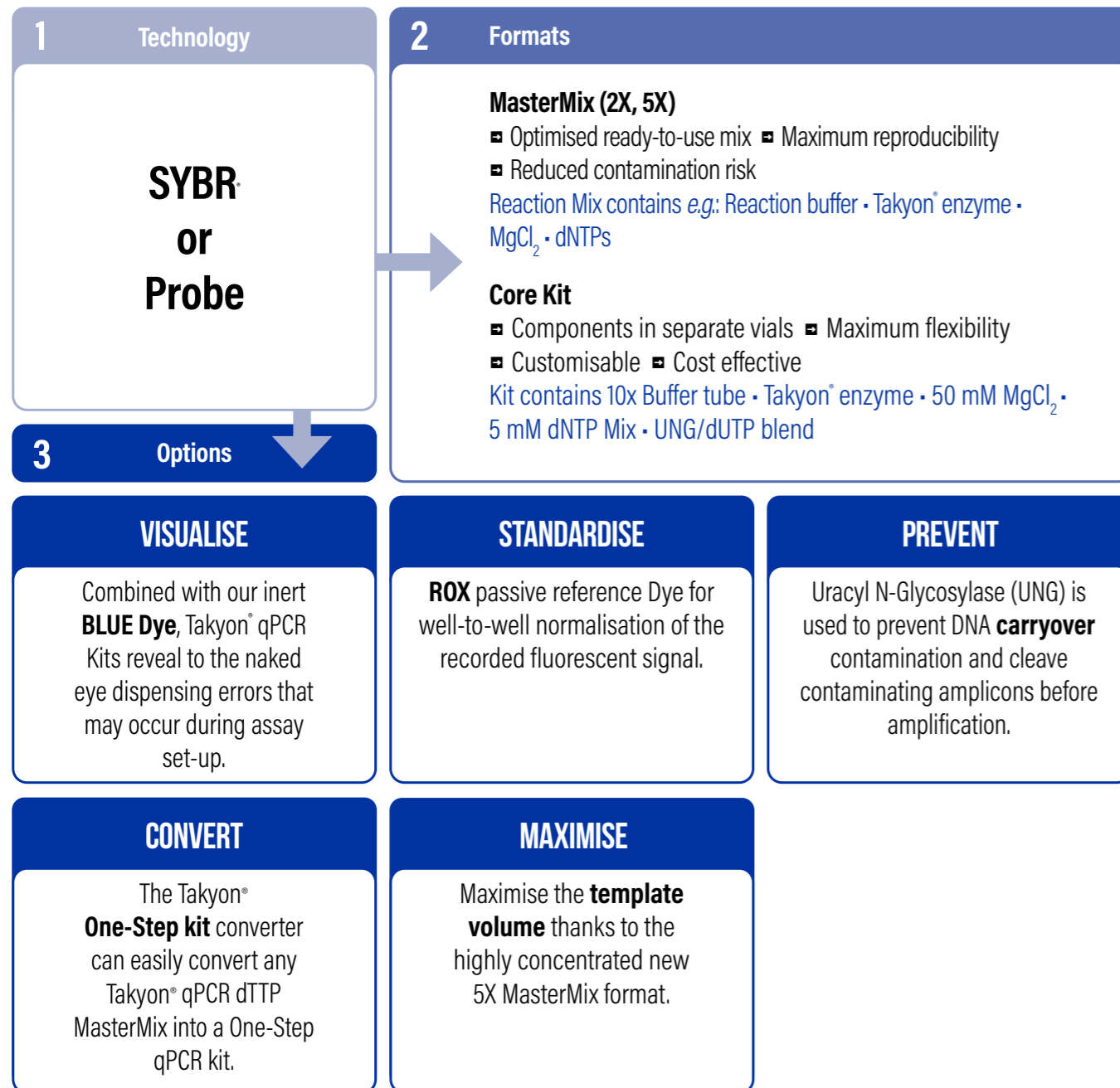
To help you in your kit selection, a convenient **online qPCR selector** is available on <http://www.eurogentec.com/qpcr-selector.html>. You can also evaluate a sample of any MasterMix for free and benefit from a 50% discount on your first kit order.

⇒ 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.

<p>Cycler Group</p> <p>A</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) <p>C</p> <ul style="list-style-type: none"> CFX384 (p.10) CFX96 (p.10) Chromo 4 (p.10) <p>D</p> <ul style="list-style-type: none"> DNA Engine Opticon® 1 & 2 (p.10) 	<p>Cycler Group</p> <p>G</p> <ul style="list-style-type: none"> GeneAmp® SDS 5700 (p.9) <p>I</p> <ul style="list-style-type: none"> iCycler iQ® (p.10) iQ™5 (p.10) <p>L</p> <ul style="list-style-type: none"> LC96 & LC480 (p.10) LightCycler® Nano System (p.10) <p>M</p> <ul style="list-style-type: none"> 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) 	<p>Cycler Group</p> <p>P</p> <ul style="list-style-type: none"> PiKoReal™ (p.10) <p>Q</p> <ul style="list-style-type: none"> Quanta® (p.10) QuantStudio® (p.9) qTower (p.10) <p>R</p> <ul style="list-style-type: none"> Rotor-Gene® 2000/3000/6000/Q (p.10) <p>S</p> <ul style="list-style-type: none"> SmartCycler® 1 & 2 (p.10) Swift Spectrum™ 96 (p.10)
---	--	---

Legend

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

For more precision, use our online selector to help you choose the right kit for your platform.

RT-qPCR kits

The Takyon® One-Step Master Mixes 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).

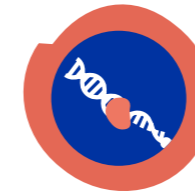
qPCR kits

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 UF-LSMT-B0100
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 UF-FSMT-B0100
	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
NO ROX	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
	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-C0201	UF-NPCT-C0205	UF-NPCT-C0510	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.

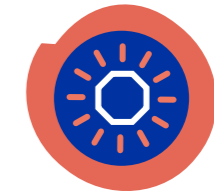
Two-step RT-qPCR kit: Use the Reverse Transcriptase Core Kit prior the qPCR assay to make a two steps reactions with any qPCR kit.



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 UF-RTAD-D0701 UF-RTAD-D0705 UF-RTAD-D0710	150 RXNs 750 RXNs 5x 750 RXNs 10x 750 RXNs
Reverse Transcriptase Core Kit 300	RT-RTCK-03 05-RTCK-03 10-RTCK-03	300 RXNs 5x 300 RXNs 10x 300 RXNs
Reverse Transcriptase Core Kit 500	RT-RTCK-05 05-RTCK-05 10-RTCK-05	500 RXNs 5x 500 RXNs 10x 500 RXNs
dUTP/UNG Additive for 2X dTTP kit conversion	RT-UTPUNG-005	40 µL (150 RXNs)
dUTP/UNG Additive for 2X dTTP kit conversion	RT-UTPUNG-020 05-UTPUNG-020 10-UTPUNG-020	200 µL (750 RXNs) 5x 200 µL (5x 750 RXNs) 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 / iQ5)	RT-FLUO-ADD	1 mL
ROX Passive Reference	RT-PARE-03 05-PARE-03 10-PARE-03 RT-PARE-100	300 µl 5x300 µl 10x300 µl 10 mL
UNG (Uracil-N-Glycosylase)	RT-0610-03 05-0610-03 10-0610-03	300 u 5x300 u 10x300 u
UNG (Uracil-N-Glycosylase)	RT-0610-15 05-0610-15 10-0610-15	1500 u 5x1500 u 10x1500 u
Takyon® Blue Additive	UF-BLUE-ADD	3750 RXN

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
	Sample Processing Control, YY-TAMRA	1000	VIC/YY	RT-SPCY-T10
	Sample Processing Control, YY-BHQ-1	200	VIC/YY	RT-SPCY-B02
	Sample Processing Control, YY-BHQ-1	1000	VIC/YY	RT-SPCY-B12
IPC	Sample Processing Control, Cy ⁵ - QXL 670	200	Cy5	RT-SPCC-Q02
	Sample Processing Control, Cy ⁵ - QXL 670	1000	Cy5	RT-SPCC-Q10
	Universal Exogenous qPCR Positive Control, YY-TAMRA	200	VIC/YY	RT-IPCY-T02
	Universal Exogenous qPCR Positive Control, YY-TAMRA	1000	VIC/YY	RT-IPCY-T10
	Universal Exogenous qPCR Positive Control, YY-BHQ-1	200	VIC/YY	RT-IPCY-B02
	Universal Exogenous qPCR Positive Control, YY-BHQ-1	1000	VIC/YY	RT-IPCY-B10
	Universal Exogenous qPCR Positive Control, Cy ⁵ - QXL 670	200	Cy5	RT-IPCC-Q02
	Universal Exogenous qPCR Positive Control, Cy ⁵ - QXL 670	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

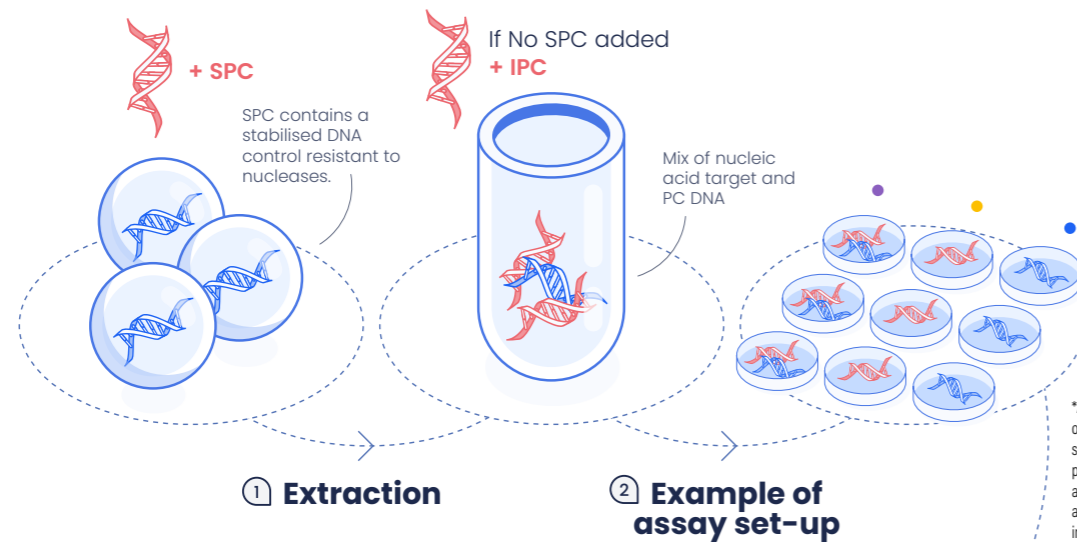
Validated in duplex assays with the target in the FAM channel

Custom labelling is available on request

METHOD

2 controls (SPC or IPC), multiple possibilities

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



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)

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.



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^q polymerase (1 u/25µl), dNTPs (400 µM), MgCl₂ (3 mM) and buffer



ORDERING INFORMATION

Description	Quantity (units)	Product Nbers
GoldStar® Mix	5 x 1 mL	PK-0064-02
Red'y'Gold Mix	5 x 1 mL	PK-0064-02R
HotGoldStar Mix	5 x 1 mL	PK-0073-02
Red'y'Star Mix	5 x 1 mL	PK-0073-02R

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. ■



HIGHLY PURE TAQ POLYMERASES

Eurogentec provides a **wide range of Taq polymerases** to meet your specific requirements, including **classical and hot-start enzymes, high-yield formulations**, and **options with or without loading dye**.

Two classical polymerases are available:

- Diamond Taq®
- Red Diamond Taq®

We also offer **HGS Diamond Taq® DNA Polymerase** for PCR applications that require ultra low levels of bacterial & fungal DNA in Diagnostic and demanding Research fields but also for amplification of templates prone to produce unspecific amplicons.

A kits contains 3 tubes : Taq® polymerase, MgCl₂ and 10X buffer.

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).



ORDERING INFORMATION

Description	Quantity	#Ref
HGS Diamond Taq®	1000 U	TAQ-I011-1000+
	5000 U	TAQ-I011-5000+
Diamond Taq®	1000 U	TAQ-I021-1000
	5000 U	TAQ-I021-5000
Red Diamond Taq®	25000 U	TAQ-I021-25000-
	1000 U	TAQ-I041-1000
Red Diamond Taq®	5000 U	TAQ-I041-5000
	25000 U	TAQ-I041-25000-



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.

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



IP-FREE OFFER

BHQ® Quenchers

Label your custom oligos with original, license-free BHQ® quenchers—available at Eurogentec for research, diagnostics, and commercial use. Compatible with most fluorescent dyes and all qPCR channels.

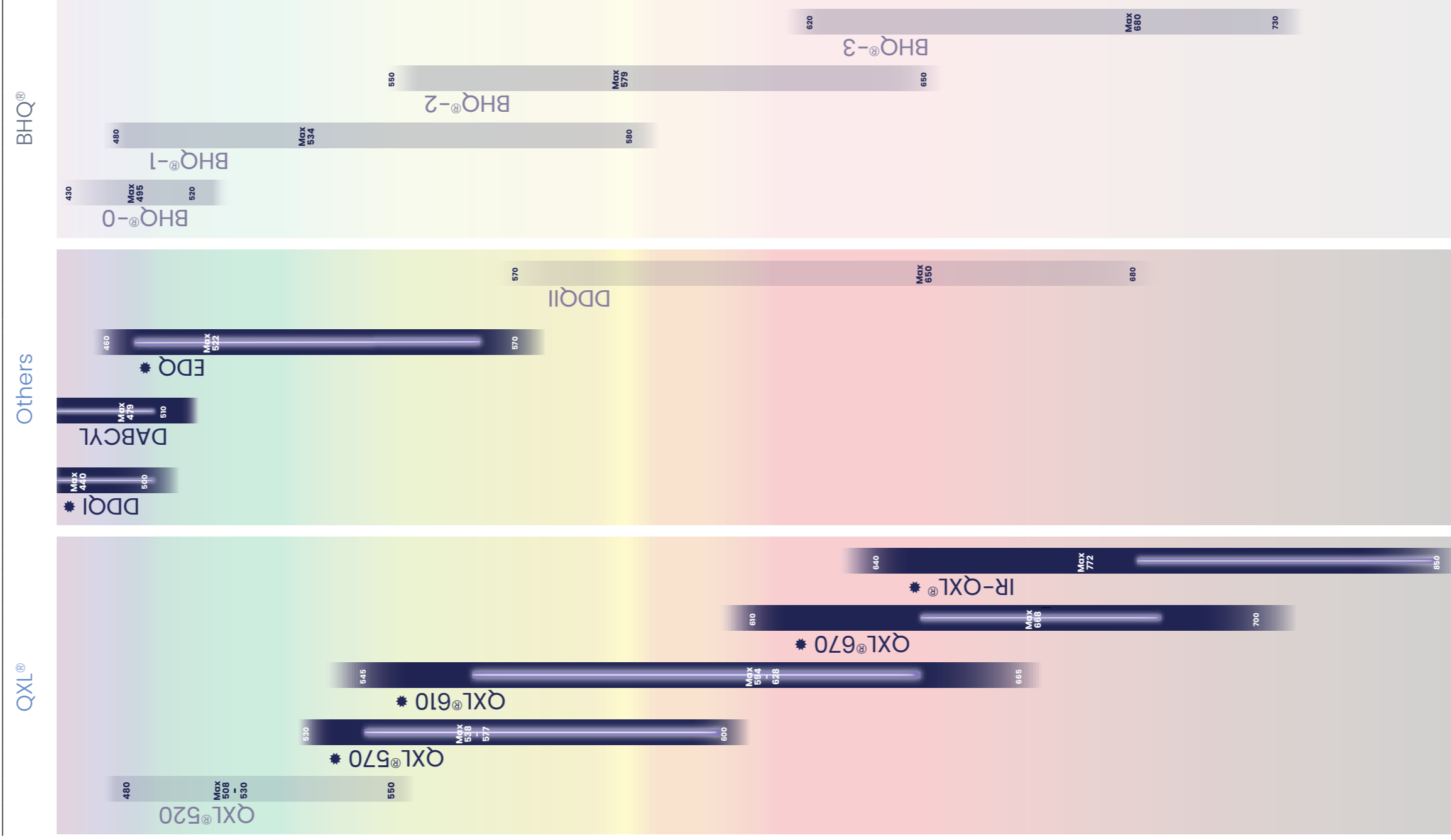
ATTO Dyes

Explore our broad selection of IP-friendly ATTO dyes with strong absorption, high quantum yield, and full-spectrum coverage—ideal for bioanalytical applications.

* 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	ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
		ATTO 520	520	538
700	UV	ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
		ATTO 568	568	586
800	UV	ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628
		ATTO 610	610	628



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 indicative quenching range of recommended quenchers
Grey bands represent the indicative 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

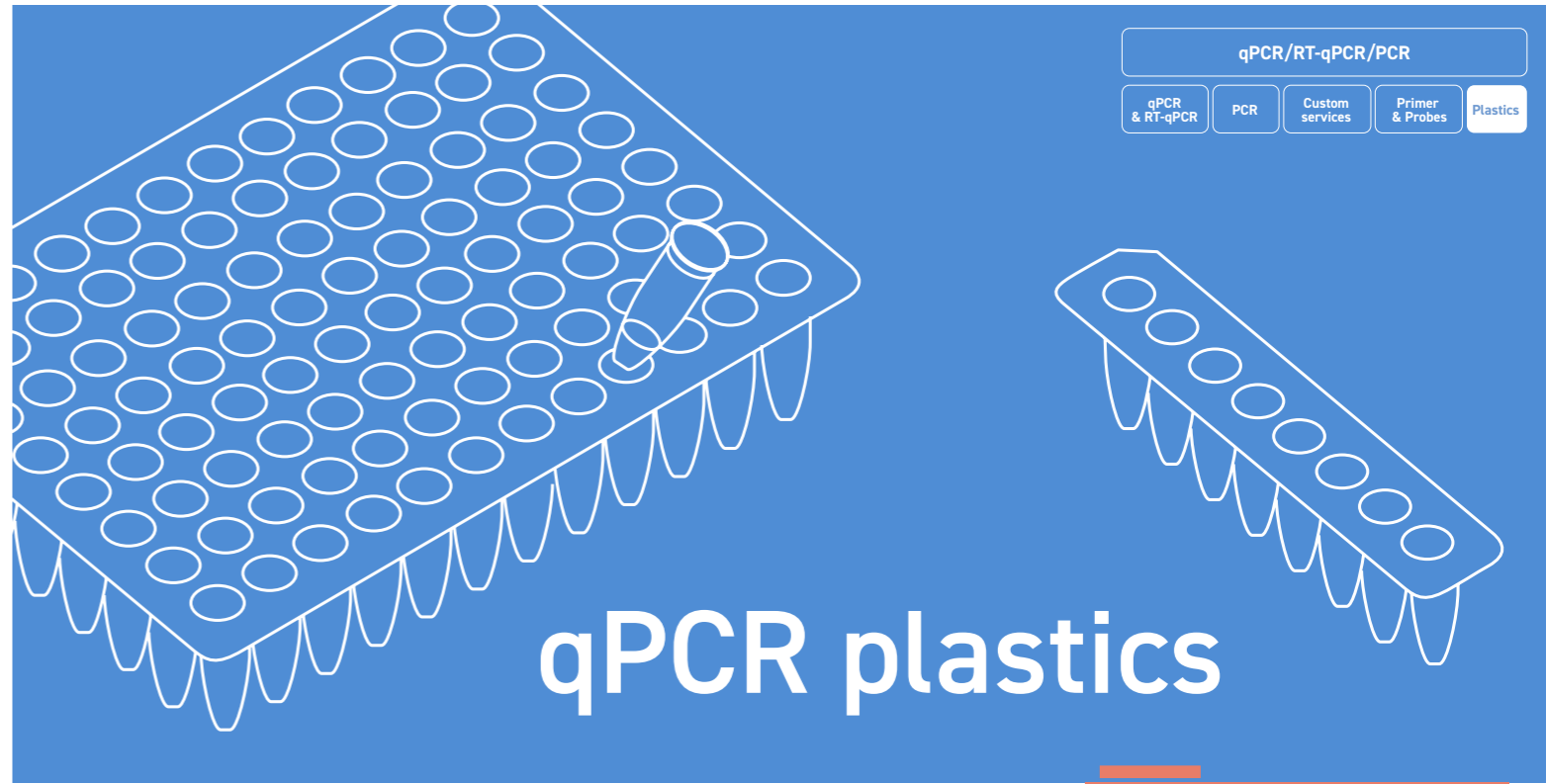
* Perform a dye calibration for optimal results
For complementary information, please refer to instrument manufacturer technical guide or contact us at scientific.support@eurogentec.com
YV = Yokima Yellow™
DFO = Dragonfly Orange™
TR = Texas Red™
In blue = Recommended by Eurogentec
In black = Available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

Thermocycler	UV- / Blue Channel	Channel 1	Channel 2	Channel 3	Channel 4	Channel 5	Channel 6	Channel 7
Real-Time PCR Thermocyclers								
AridMX & AridX (up to 6 filters)	ATTO 425	FAM	YY*/HEX*/JOE*/ATTO 532*	ATTO 550*/TAMRA/Cy 3	ATTO 590*/ROX/TR	ATTO 647N*/Cy5		
qTOWER*		FAM	YY*/HEX*/JOE*/VIC	DFO*/TAMRA/Alexa546/NED	ROX/TR/Cy 3.5	ATTO 647N*/Alexa633/Cy5	Cy 5.5	
qTOWERiris	ATTO 425/ATTO 390*	FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/TAMRA/DFO*	ATTO 590/ROX/TR/Cy 3.5	ATTO 633*/Cy5	Cy 5.5/ATTO 665	
ABI Prism® SDS 7500		FAM	YY*/HEX*/JOE*/TET*/VIC	DFO*/TAMRA/Cy 3/NED	ROX/TR	ATTO 647N*/HL647/Cy5		
ABI Prism® SDS 7900HT		FAM	YY*/JOE*/TET*/VIC	DFO*/TAMRA/NED	ROX			
QuantStudio™ 1		FAM	YY*/HEX*/JOE*/TET*/VIC	ROX/TR/JUN				
QuantStudio™ 3		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy3*	ROX/ATTO 590*/TR*			
QuantStudio™ 5 (384-W.) & 6 Flex/Pro		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy3*	ROX/ATTO 590*/TR*	ATTO 647N*/HL647*/Cy5		
QuantStudio™ 5 (96-W.) & 7 Flex/Pro & 12k Flex		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy3*	ROX/ATTO 590*/TR*	ATTO 647N*/HL647*/Cy5	ATTO 680*/Cy5.5*	
StepOnePlus		FAM	YY*/HEX*/JOE*/TET*/VIC	DFO*/TAMRA/Cy 3*/NED	ROX			
VIA7		FAM	YY*/HEX*/JOE*/TET*/VIC	ATTO 550*/DFO*/TAMRA/Cy3*	ROX/ATTO 590*/TR*	ATTO 647N*/HL647*/Cy5	ATTO 680*/Cy5.5*	
CFX Connect & CFX Duet		FAM	YY*/HEX*/TET*/JOE*					
CFX Opus & CFX96 Touch		FAM	YY*/HEX*/TET*/JOE*	TR/ROX*/ATTO 590*	ATTO 647N*/Cy5	ATTO 700*/Cy5.5*		
CFX384 Touch		FAM	YY*/HEX*/TET*/JOE*	TR/ROX*/ATTO 590*	ATTO 647N*/Cy5			
iCycler iQ™		FAM	YY*/HEX*/Cy 3*/JOE*/TET	ATTO 550*/DFO*/TAMRA/Cy3	ROX/TR/ATTO 590*	ATTO 647N*/Cy5		
iQ™5		FAM	YY*/HEX*/Cy3*/JOE*/TET	ATTO 550*/DFO*/TAMRA/Cy3	ROX/TR/ATTO 590*	ATTO 647N*/Cy5		
QIAquant 96 & 384 Splex		FAM	YY*/HEX*/JOE*/VIC	ROX/TR/ATTO 590*	ATTO 647N*/Cy5	Cy 5.5/ATTO 680*		
QIAquant 96 2plex		FAM	YY*/HEX*/JOE*/VIC					
Rotor-Gene® 6000 & Q	ATTO 390*	FAM	YY*/HEX*/JOE*/TET*/VIC	ROX/TR/Cy 3.5/ATTO 590*	ATTO 647N*/Cy5/Alexa633	ATTO 680*/Alexa680		
LightCycler® 480 I	ATTO 425*/Cyan 500	FAM	YY*/HEX*/JOE*/VIC	ROX*/TR*/LC610	ATTO 620/LC640	ATTO 633*/Cy5		
LightCycler® 480 II	ATTO 425*/Cyan 500	FAM	YY*/HEX*/JOE*/VIC	ROX*/TR*/LC610	ATTO 620/LC640	ATTO 633*/Cy5/Cy5.5		
LightCycler® 96		FAM	YY*/HEX*/VIC	TR/ROX*/ATTO 590*/LC610	ATTO 647N*/Cy5			
LightCycler® PRO	ATTO 425*/Cyan 500	FAM	YY*/HEX*/JOE/R6G*/VIC	ROX*/TR*/ATTO 590*/LC610	ATTO 633*/LC640/JA 270	ATTO 647N*/Cy5	Cy 5.5/ATTO 700*	
Mic		FAM	YY*/HEX*/JOE*/VIC	ATTO 590/ROX/TR	ATTO 647N/HL647/Cy5			
Digital PCR Platforms								
Naica® Prism3		FAM	HEX/Cy 3*/VIC	ATTO 647N*/Cy5				
Naica® Prism6		FAM	YY	ATTO 550/TAMRA	ROX	ATTO 647N*/Cy5	ATTO 700/Cy5.5	
Nio & Nio+		FAM	YY*/HEX*/ATTO 532*/JOE*	ATTO 550/TAMRA/Cy3	ROX/TR*/ATTO 590*	ATTO 647N*/Cy5	Cy 5.5/ATTO 700*	DY-521-XL
QX One ddPCR		FAM	YY*/HEX*/VIC	ATTO 647N*/Cy5	Cy 5.5			
QX100 & QX200 ddPCR		FAM	YY*/HEX*/VIC					
QX600 ddPCR		FAM	YY*/HEX*/VIC	ATTO 647N*/Cy5	Cy 5.5	ROX	ATTO 590	
QuantStudio™ Absolute Q		FAM	HEX/YY*/VIC	ATTO 550*/TAMRA*/ABY	ROX/ATTO 590*/TR*	ATTO 633*/Cy5/JUN		
QIAcquity Four/Eight (Splex)		FAM	YY*/HEX*/JOE*/VIC	TAMRA/ATTO 550	ROX/TR/ATTO 590*	ATTO 647N*/Cy5		
QIAcquity One (2plex)		FAM	YY*/HEX*/JOE*/VIC					
Digital LightCycler®	ATTO 425*/Cyan 500	FAM	HEX*/VIC	TR/CFR610/LC610	ATTO 647N*/Cy5/JA270/LC640	Cy 5.5		
DQ24 Plus	ATTO 425	FAM	YY*/HEX*/JOE*/VIC/SUN	ROX/TR	ATTO 647N*/Cy5/Q670	Cy 5.5/Quasar 705		

i For complementary information, please refer to instrument manufacturer technical guide or contact us at scientific.support@eurogentec.com

* For optimal results in multiplex assays, a dye calibration may be required to avoid fluorescence cross-talk

in bold = Recommended by Eurogentec
 in regular black = Available at Eurogentec
 in grey = Not 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.

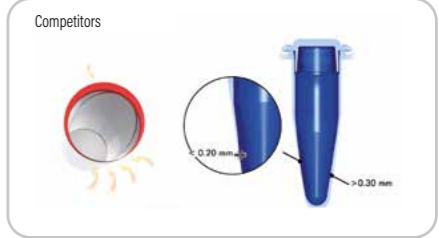
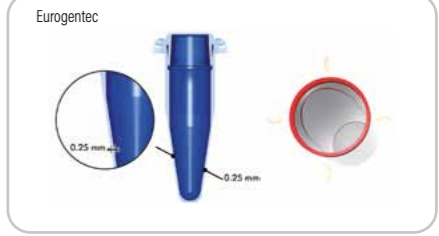
Please refer to the annexes to know which plastics is the most suited for your thermocycler. (pages 28 → 30) ■

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, DNase, RNase, 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 a well.

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-MQ	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

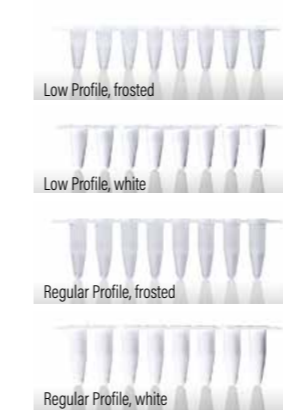
The extremely clear glass-grade area of the cap ensures very good transmission of the fluorescence and equals the performance properties of optical seals. When combined with our qPCR plates they seal better than other brands, thus less evaporation.

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

Compatibility of plastic plates and tubes with Real-Time PCR thermocyclers

qPCR 96-well plate type	Regular block		FAST block		Regular block		FAST block										
	96-well blocks (1, 3, 5, 6&7 Flex/PRO, 12k Flex)	QuantStudio w/ standard (0.2 ml) 12k Flex)	QuantStudio w/ FAST (0.1 ml) 96-well blocks (3, 5, 6&7 Flex/PRO, 12k Flex)	QuantStudio w/ 384-well blocks (5, 6&7 Flex/PRO, 12k Flex)	Agilent AriaIIX & AriaIIX	Analytik Jena qTOWER ³ & qTOWER ³ HiS	Bio-rad iCycler iQ-105 & MyiQ ³	Bio-rad CFX96 & CFX Opus/Connect 96 & CFX Duet	Bio-rad CFX384 & CFX Opus/Connect 384	Qiagen QIAquant (2- & 5-plex)	Roche LightCycler [®] 480 (I & II)	Roche LightCycler [®] 96	ABI Prism [®] SDS 7500	ABI Prism [®] SDS 7900HT	ABI StepOne [™] Plus	ABI ViiA7	
Cover																	
Sub-skirted, high profile, frosted or white type (ABI systems) RT-PL96-AB or RT-PL96-ABW	*	*											*	*		*	*
Sub-skirted, low profile, frosted or white type (ABI systems) RT-PL96-AF or RT-PL96-AFW			*			*									*		*
Non-skirted, high profile, natural or white type RT-PL96-MQ or RT-PL96-MCW	✓	✓				*							✓	✓			
Non-skirted, low profile, frosted or white type RT-PL96-OP or RT-PL96-OPW			✓			*											
Semi-skirted, low profile, white RT-PL96-LC480							*	*	*	*	*	*					
qPCR 384-well plate type																	
Skirted, low profile, white plate for ABI RT-PL384-ABW				*				*	*	*	*	*	*	*	*	*	*
Skirted, low profile, white plate for LC RT-PL384-LCW																	
qPCR 8-tube strip type																	
Extra Robust, Regular Profile, Frosted, 0.2 ml RT-STR8-RH20	✓	✓				*							✓	✓		✓	✓
Extra Robust, Regular Profile, White, 0.2 ml RT-STR8-RW120	✓	✓				*							✓	✓		✓	✓
Extra Robust, Low Profile, Frosted, 0.1 ml RT-STR8-LF120																	
Extra Robust, Low Profile, White, 0.1 ml RT-STR8-LW120																	

* recommended for optimal compatibility / ▲ adaptor required (ref: RT-PL96-LCAD) / ✓ compatible plastics / ✓ Refer toycler manufacturer instructions for required adaptor
 qPCR optical semi-domed 8-cap strips are also available (ref: RT-OCAP-300; 300 cap strips) and can be used on several machines.

Need help finding the right plastic vessel?
 Contact us at scientific.support@eurogentec.com

HOW TO ORDER

ONLINE

ON 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.

TRADEMARKS/LABELS

ABI Prism[®] 5700, 7000, 7500, 7700 and 7900 are registered trademarks of Applied Biosystems, Inc.
Acrydite[®] is a registered trademark of Mosaic Technologies Corp.
Alexa Fluor[®] is a registered trademark of Molecular Probes, Inc.
AnaSpec[®] is a registered trademark of Anaspec, Inc.
AnaTag[™] is a trademark of Anaspec, Inc.
Benzonase[®] is a registered trademark of Merck KGaA, Darmstadt, Germany.
BHQ[®] is a registered trademark of Biosearch Technologies, Inc. (Novato, California, USA)
BlackBerry[®] is a registered trademark of Berry & Associates, Inc.
Black Hole Quencher[®] is a registered trademark of Biosearch Technologies, Inc. (Novato, California, USA)
BODIPY[®] is a registered trademark of Life Technologies Corp.
CAL Fluor[®] is a registered trademark of Kaneka Eurogentec S.A.
Cascade Blue[®] is a registered trademark of Life Technologies Corp.
CDP Star[®] is a registered trademark of Applera Corp.
Cy[®] and CyDye[®] are registered trademarks of Cytiva.
CyLyte Fluor[™] is a trademark of Anaspec, Inc.
DHL[™] is a trademark of Anaspec, Inc.
Diamond Tag[®] is a registered trademark of Kaneka Eurogentec S.A.
DNA Engine Option[®] is a registered trademark of Bio-Rad Laboratories, Inc.
Dragonfly Orange[™] is a trademark of Kaneka Eurogentec S.A.
DyLight[®] is a registered trademark of Thermo Fisher Scientific Inc. and its subsidiaries.
Eclipse[™] Dark Quencher is a trademark of ElitechGroup, Inc.
Eclipse[™] Quencher is a trademark of ElitechGroup, Inc.
EGT[®] is a registered trademark of Kaneka

Eurogentec S.A.

Eppendorf[®] is a registered trademark of Eppendorf AG.
Excel[™] is a trademark of Microsoft Corp.
Eurogentec[®] is a trademark of Kaneka Eurogentec S.A.
Exiqon[®] is a registered trademark of Exiqon A/S.
FLAG[®] is a registered trademark of Sigma-Aldrich Co.
GeneAmp[®] is a registered trademark of Roche Molecular Systems, Inc.
GenePix[®] is a registered trademark of Astro Acquisition Sub II.
GenePix 4000[®] is a registered trademark of Astro Acquisition Sub II.
Gold'N Oligo[®] is a registered trademark of Kaneka Eurogentec S.A.
GO[®] is a registered trademark of Anaspec, Inc.
GoldStar[®] is a registered trademark of Kaneka Eurogentec S.A.
Hilyte[™] is a trademark of Anaspec, Inc.
Hilyte fluor[™] is a trademark of Anaspec, Inc.
iCycler iQ[®] is a registered trademark of Bio-Rad Laboratories, Inc.
iCycler[®] is a registered trademark of Bio-Rad Laboratories, Inc.
I-MUPID[®] is a registered trademark of Advance Co., Ltd.
iQ[™]5 is a trademark of Bio-Rad Laboratories, Inc.
iQ-CHECK[®] is a registered trademark of Bio-Rad Laboratories, Inc.
LightCycler[®] is a registered trademark of Roche Diagnostics GmbH.
LNA[®] is a registered trademark of Exiqon A/S.
Lumavidin[®] is a registered trademark of Luminex Corp.
Luminex[®] is a registered trademark of Luminex Corp.
Luminex[®] 200[™] is a trademark of Luminex Corp.
Luminex[®] XYP[™] is a trademark of

Luminex Corp.

Luminex[®] SD[™] is a trademark of Luminex Corp.
Luminex[®] SeroMAP[™] is a trademark of Luminex Corp.
MagPlex[™] is a trademark of Luminex Corp.
MapMarker[®] is a registered trademark of BioVentures, Inc.
Marina Blue[®] is a registered trademark of Life Technologies Corp.
MasterCycler[®] is a registered trademark of Eppendorf, Inc.
MGB[™] is a trademark of ElitechGroup, Inc.
MGB Pleiades[™] is a trademark of ElitechGroup, Inc.
MGB Eclipse[®] is a registered trademark of ElitechGroup, Inc.
MicroPlex[™] is a trademark of Luminex Corp.
Mupid[®] is a registered trademark of Advance Co., Ltd.
Mx3000[®] is a registered trademark of Stratagene
Mx3000P[®] is a registered trademark of Stratagene
Mx3005P[®] is a registered trademark of Stratagene
Mx4000[®] is a registered trademark of Stratagene
Onco-Mir[®] is a registered trademark of Ambion Inc.
Opticon[®] is a registered trademark of MJ Research, Inc.
Oregon Green[®] is a registered trademark of Life Technologies Corp.
Oyster[®] is a registered trademark of Denovo Biolabels
Pacific Blue[™] is a trademark of Life Technologies Corp.
Pulsar[®] is a trademark of Biosearch Technologies, Inc.
Quantica[®] is a registered trademark of Techné Ltd and Techné, Inc.
Quasar[®] is a trademark of Biosearch Technologies, Inc.

Qiagen[®] and **QiaQuick[®]** are trademarks of Qiagen, Inc.
QSY[®] is a registered trademark of Life Technologies Corp.
QXL[®] is a registered trademark of Luminex Corp.
Rhodamine Red[™] is a trademark of Life Technologies Corp.
Rotor-Gene[®] is a registered trademark of Corbett Research Pty Ltd
ROX[™] is a trademark of Applied Biosystems
RP-Cartridge Gold[™] is a trademark of Eurogentec S.A.
SELEX[™] is a trademark of Gilead Sciences, Inc.
Sensolyte[®] is a registered trademark of Anaspec, Inc.
Smartcyber[®] is a registered trademark of ElitechGroup, Inc.
StepOne[®] is a registered trademark of Applied Biosystems
SuperScript[®] is a registered trademark of Invitrogen Corp.
SYBR[®] Green is a registered trademark of Life Technologies Corp.
Takyon[®] is a registered trademark of Kaneka Eurogentec S.A.
Taqman[®] is a registered trademark of Roche Diagnostics GmbH.
Texas Red[®] is a registered trademark of Life Technologies Corp.
TRIZOL[®] is a registered trademark of Molecular Research Center Inc.
VersaArray[®] is a registered trademark of Bio-Rad Laboratories, Inc.
Vic[®] is a registered trademark of Applied Biosystems
Word[™] is a trademark of Microsoft Corp.
xPONENT[®] is a registered trademark of Luminex Corp.
xMAP[®] is a registered trademark of Luminex Corp.
Yakima Yellow[®] is a registered trademark of ElitechGroup, Inc.
Z-Fish[®] is a registered trademark of Anaspec, Inc.

BE AWARE OF

EUROGENTEC LICENSE STATEMENT



SCAN AND READ THEM

www.eurogentec.com

PCR/qPCR

Contact:

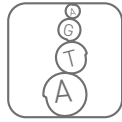
For more information,
please contact your Sales
representative or contact us at:
isales@eurogentec.com

Kaneka Eurogentec S.A. 2025

Follow us on



Life Science Research



Oligos



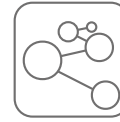
PCR/qPCR



Genes/DNA



Antibodies



Peptides



Proteins