EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00 • Fax: +32 4 372 75 00 Toll-free: +800 666 00 123 • info@eurogentec.com • www.eurogentec.com



NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560 Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • service@anaspec.com www.anaspec.com

Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

Technical Data Sheet

ROX Passive Reference RT-PARE-03 • RT-PARE-100 • 05-PARE-03 • 10-PARE-03

[300 μ L] [10 mL] [5 x 300 μ L] [10 x 300 μ L]

Eurogentec products are sold for research or laboratory use only and are not to be administrated to humans or used for medical diagnostics.

Description

The ROX passive reference consists of an inert dye whose fluorescence is stable along the qPCR reaction, allowing the normalization of well-to-well variations that may result from pipetting errors or instrument limitations.

Eurogentec's ROX passive reference has been specifically formulated for use on ABI, Stratagene and Eppendorf Real-Time thermocyclers. Any of our No ROX and Low ROX kits can be used in combination with a separate tube of the ROX passive reference.

Package content

RT-PARE-03: 300 µl of ROX passive reference. RT-PARE-100: 10 mL of ROX passive reference. 05-PARE-03: 5 x 300 µl of ROX passive reference. 10-PARE-03: 10 x 300 µl of ROX passive reference.

Storage and shipping condition

- For long term storage the ROX passive reference should be stored at -15°C to -25°C in a constant temperature freezer. When stored under these conditions the reagents are stable for 2 years. Avoid repetitive freeze-thaw cycles.
- ROX passive reference is stable at 4°C for one month.
- Store in the dark. Minimizing exposure to direct light avoid loss of fluorescent signal intensity.
- Shipped on dry ice.

Protocol

The following table indicates the appropriate quantity of ROX passive reference to add in Takyon™ MasterMix depending on the thermocycler used, the Mix volume and concentration.

Starting Takyon™ MasterMix	No ROX		Low ROX	
Platform used	2 x	5 x	2 x	5 x
Mx3000P® Mx3005P® Mx4000®	8 μL / mL or 60 μL / 7.5 mL	20 μL / mL	N.A.	N.A.
ABI Prism® 7500 ABI Prism® FAST 7500 ViiA7™ QuantStudio™	2.8 μL / mL or 21 μL / 7.5 mL	7μL/mL	N.A	N.A
ABI Prism® 5700 ABI Prism® 7000 ABI Prism® 7300 ABI Prism® 7700 ABI Prism® 7900 ABI Prism® FAST 7900 ABI Step One ABI Step One Plus MasterCycler® ep realplex I & II (Firmware 2.1 or below)	40 μL / mL or 300 μL / 7.5 mL	100 μL / mL	37.2 μl / ml or 279 μl / 7.5 ml	93 μl / ml or 697.5 μl / 7.5 ml

For further information please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com

Tel: +32 4 372 76 65 • Toll-free: +800 666 00 123

For USA:

E-mail: service@anaspec.com

Tel.: +1-510-791-9560 • Toll-free: +1-800-452-5530

ABI Prism® 5700, 7000, 7500, 7700 and 7900 are registered trademarks of Applied Biosystems, Inc. iCycler iQ® is a registered trademark of Bio-Rad Laboratories, Inc. iCycler® is a registered trademark of Bio-Rad Laboratories, Inc. iQ¹™S is a trademark of Bio-Rad Laboratories, Inc.

Mx3000P°, Mx3005P° and Mx4000° are registered trademarks of Stratagene

MasterCycler® is a registered trademark of Eppendorf, Inc

UDG Use of UDG employs patent rights licensed from Invitrogen Corporation.

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.