

AMPIGENE[®] qPCR 1-Step Probe Kit Hi-ROX

Optimized and efficient cDNA synthesis and probe-based qPCR in a single tube

AMPIGENE[®] qPCR 1-Step Probe Kits use the latest developments in reverse transcriptase (RTase) technology and buffer chemistry for efficient cDNA synthesis and qPCR in a single tube.

The modified MMLV RTase is blended with RNase inhibitor to prevent degradation of RNA by contaminating RNase. The RTase is thermostable and extremely active.

AMPIGENE[®] qPCR 1-Step Probe Kits are designed for use on a variety of probes including: Taqman[®], Molecular Beacons[®], and Scorpion probes[®].

An optimized buffer system allows efficient amplification and proprietary enhancers prevent the formation of primers/dimers to increase sensitivity and specificity.

ROX is a reference dye which is used to normalize fluorescent signal in qPCR. AMPIGENE[®] qPCR 1-Step Probe Kits come with either Lo-ROX, Hi-ROX, or No-ROX concentration. For Lo-ROX and No-ROX, see Prod. Nos. [ENZ-NUC110](#) and [ENZ-NUC112](#). Choose the appropriate ROX concentration for the instrument (see instrument compatibility for reference dye chart below.)

Ordering Information

[Order Online »](#)

ENZ-NUC111-0200	200Reactions
ENZ-NUC111-1000	1000Reactions

Manuals, SDS & CofA

[View Online »](#)

- Convenient one tube reaction
- Thermostable and extremely active RTase
- For probe-based qPCR applications: Designed for used on a wide range of probe technologies
- Optimized buffer system for efficient amplification
- Enhancers and stabilizers improve reaction sensitivity and specificity

**INSTRUMENT COMPATIBILITY
For Reference Dye**

Manufacturer	Instrument	Lo-ROX	Hi-ROX	No-ROX
Analytica Jena	qTower	Yes	Yes	Yes
Applied Biosystems	7500, 7500 FAST, Viia7™	Yes	No	Yes
Applied Biosystems	7000, 7300, 7700, 7900, 7900HT, 7900HT FAST, StepOne™, StepOne™ Plus	No	Yes	No
Bio-Rad®	iCycler®, MyiQ®, iQ™5, Opticon™, Opticon™2, Chromo4™, MiniOpticon™, CFX96™, CFX384™	Yes	No	Yes
Cepheid®	Smartcycler®	Yes	Yes	Yes
Eppendorf	Mastercycler® ep realplex, Mastercycler® realplex 2S	Yes	Yes	Yes
Illumina®	Eco™	Yes	Yes	Yes
Qiagen/Corbett	Rotor-Gene™ 3000, 6000, Q	Yes	Yes	Yes
Roche Applied Science	Lightcycler® 480, Lightcycler® Nano	Yes	Yes	Yes
Stratagene (Agilent)	MX 4000P®, MX 3000P®, MX 3005P®	Yes	No	Yes
Takara	Cycler Dice®	Yes	Yes	Yes
Techne	Quantica®	Yes	Yes	Yes

Instrument compatibility for ROX dye.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Application Notes For cDNA synthesis and probe-based qPCR in a single tube

Contents 2x AMPIGENE[®] qPCR 1-Step Probe Mix Hi-ROX (includes dNTPs and polymerases)
20x RTase (with RNase inhibitor)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)