

# AMPIGENE<sup>®</sup> qPCR 1- Step Green Kit Hi-ROX

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

AMPIGENE<sup>®</sup> qPCR 1-Step Green 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<sup>®</sup> qPCR 1-Step Green Kits use a proprietary intercalating dye that does not inhibit PCR.

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<sup>®</sup> qPCR 1-Step Green Kits come with either Lo-ROX or Hi-ROX concentration. For Lo-ROX, see Prod. No. [ENZ-NUC108](#). Choose the appropriate ROX concentration for the instrument (see instrument compatibility for reference dye chart below).

## Ordering Information

[Order Online »](#)

ENZ-NUC109-1000	1000Reactions
-----------------	---------------

## Manuals, SDS & CofA

[View Online »](#)

- Convenient one tube reaction
- Thermostable and extremely active RTase
- For dye-based qPCR applications: Green dye does not interfere with PCR
- 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
Analytica Jena	qTower	Yes	Yes
Applied Biosystems	7500, 7500 FAST, Viiia7™	Yes	No
Applied Biosystems	7000, 7300, 7700, 7900, 7900HT, 7900HT FAST, StepOne™, StepOne™ Plus	No	Yes
Bio-Rad®	iCycler®, MyiQ®, iQ™5, Opticon™, Opticon™2, Chromo4™, MiniOpticon™, CFX96™, CFX384™	Yes	No
Cepheid®	Smartcycler®	Yes	Yes
Eppendorf	Mastercycler® ep realplex, Mastercycler® realplex 2S	Yes	Yes
Illumina®	Eco™	Yes	Yes
Qiagen/Corbett	Rotor-Gene™ 3000, 6000, Q	Yes	Yes
Roche Applied Science	Lightcycler®480, Lightcycler®Nano	Yes	Yes
Stratagene (Agilent)	MX 4000P®, MX 3000P®, MX 3005P®	Yes	No
Takara	Cycler Dice®	Yes	Yes
Techne	Quantica®	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 dye-based qPCR in a single tube

Contents 2x AMPIGENE<sup>®</sup> qPCR 1-Step Green Mix Hi-ROX  
(includes dNTPs and polymerases)  
20x RTase (with RNase inhibitor)

Last modified: May 29, 2024



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)