

# AMPIGENE<sup>®</sup> qPCR 1-Step Green Kit Lo-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 Hi-ROX, see Prod. No. [ENZ-NUC109](#). Choose the appropriate ROX concentration for the instrument (see instrument compatibility for reference dye chart below.)

## Ordering Information

[Order Online »](#)

|                 |               |
|-----------------|---------------|
| ENZ-NUC108-0200 | 200Reactions  |
| ENZ-NUC108-1000 | 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, Viia7™                                                             | 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 Lo-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)