

AMPIGENE® 1-Step RT-PCR Kit

Optimized and efficient cDNA synthesis and PCR in a single tube

AMPIGENE® 1-Step RT-PCR Kit uses the latest developments in reverse transcriptase (RTase) technology and buffer chemistry for efficient cDNA synthesis and PCR 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.

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-KIT105-0050	50Reactions
ENZ-KIT105-0200	200Reactions
ENZ-KIT105-1000	1000Reactions

Manuals, SDS & CofA

[View Online »](#)

- Convenient one tube reaction
- Thermostable and extremely active RTase
- Optimized buffer system for efficient amplification
- Enhancers and stabilizers improve reaction sensitivity and specificity

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 PCR in a single tube

Contents 2x 1-Step Mix (AMPIGENE[®] HS Taq Polymerase, 6mM MgCl₂, 2mM dNTPs, enhancer, and stabilizers)
20x RTase (with RNase inhibitor)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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