

AMPIGENE[®] cDNA

Synthesis Kit

Efficient and sensitive cDNA synthesis for downstream qPCR

AMPIGENE[®] cDNA Synthesis Kit uses the latest developments in reverse transcriptase (RTase) technology and buffer chemistry for efficient and accurate cDNA synthesis. 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 contains oligo(dT), random hexamers, enhancers, dNTPs and MgCl₂, allowing efficient and sensitive cDNA synthesis.

Citations: 18

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-KIT106-0050	50Reactions
ENZ-KIT106-0200	200Reactions

Manuals, SDS & CofA

[View Online »](#)

- Unbiased, efficient, sensitive cDNA synthesis
- Thermostable and extremely active RTase
- Enhanced buffer system for efficient synthesis
- Optimized random hexamer and anchored oligo(dT) concentrations for generation of cDNA for qPCR

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

For cDNA synthesis for downstream qPCR reactions

Contents

5x cDNA synthesis mix (anchored oligo (dT), random hexamers, 15mM MgCl₂, 5mM dNTPs, enhancers, 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