

AMPIGENE[®] Lyo- Probe 1-Step kit

AMPIGENE[®] Lyo-Probe 1-Step Kit
formulated for reliable detection of viral RNA
sequences and development of lyophilized
diagnostic assays

AMPIGENE[®] Lyo-Probe 1-Step Kit was designed for development of highly sensitive RT-qPCR detection of RNA sequences. Lyophilization offers advantages such as ambient storage, extended shelf-life, and versatility in sample volume.

Manuals, SDS & CofA

[View Online »](#)

- Same high performance as our standard kits
- Separate, high concentration AMPIGENE[®] Reverse Transcriptase (with RNase inhibitor) included
- Compatible with high-throughput and high-multiplexed assays

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Contents 4x AMPIGENE[®] Lyo-Probe Mix, 3200x AMPIGENE[®] RT 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