

# 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.

## Ordering Information

[Order Online »](#)

ENZ-NUC143-2000	2000Reactions
-----------------	---------------

## 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<sup>®</sup> Lyo-Probe Mix, 3200x AMPIGENE<sup>®</sup> RT 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)