

AMPIGENE[®] Pre Lyo- Probe 1-Step Evaluation Kit

Version of AMPIGENE[®] Lyo-Probe 1-Step Kit (ENZ-NUC143)

AMPIGENE[®] Lyo-Probe 1-Step Evaluation 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-NUC142-0200	200Reactions
-----------------	--------------

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, 20x AMPIGENE[®] RT with RNase inhibitor



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)