

# AMPIGENE<sup>®</sup> Lyo- Probe Mix

Formulated for development of lyophilized  
diagnostic assays targeting viral RNA  
sequences

AMPIGENE<sup>®</sup> Lyo-Probe Mix 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

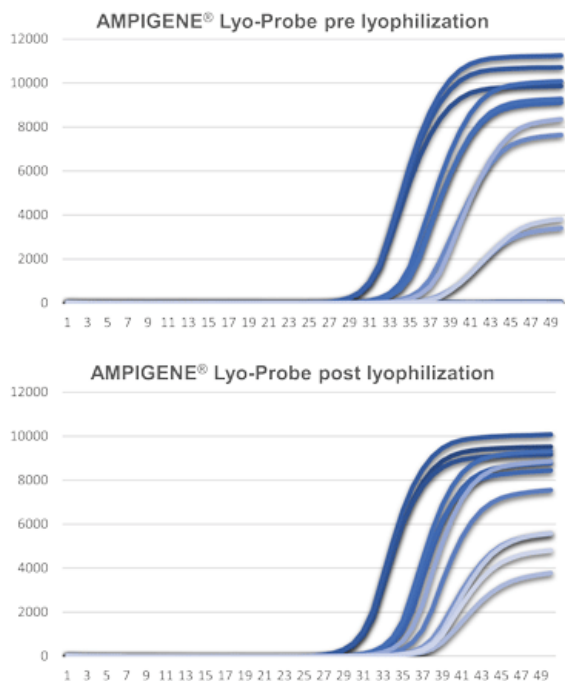
[Order Online »](#)

ENZ-NUC141-0200	200Reactions
ENZ-NUC141-1000	1000Reactions

## Manuals, SDS & CofA

[View Online »](#)

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



### **AMPIGENE® Lyo-Probe Mix Performs Similarly Before Lyophilization and After Reconstitutions**

A singleplex assay on the COVID-19 E-Gene confirmed that there are no major differences in AMPIGENE® Lyo-Probe Mix performance before lyophilization and after reconstitution.

## 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



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)