AMPIGENE[®] Lyo- Probe Mix

Formulated for development of lyophilized diagnostic assays targeting viral RNA sequences

AMPIGENE® 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

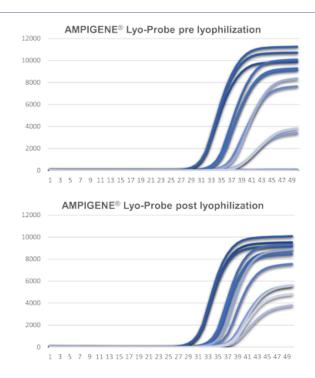
Order Online »

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

Manuals, SDS & CofA

View Online »

- Same high performance as our standard kits
- Separate, high concentration AMPIGENE[®]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

4x AMPIGENE® Lyo-Probe Mix **Contents**

Last modified: May 29, 2024



info-

eu@enzolifesciences.com