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

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

with RNase inhibitor

Last modified: May 29, 2024



info-

eu@enzolifesciences.com