

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.

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

