

Technical Product Specification

Quant Nova® Probe RT-PCR Kit 208352 (100), 208354 (500), or 208356 (2500) reactions



Product description

For 100 x 20 µL reactions: 1 mL Quant Nova Probe RT-PCR Master Mix, 20 µL Quant Nova Probe RT Mix, 20 µL Internal Control RNA, 500 µL Yellow Template Dilution Buffer, 250 µL ROX Reference Dye, 1.9 µL RNase-Free Water

For 500 x 20 µL reactions: 3 x 1.7 mL Quant Nova Probe RT-PCR Master Mix, 100 µL Quant Nova Probe RT Mix, 100 µL Internal Control RNA, 500 µL Yellow Template Dilution Buffer, 1 mL ROX Reference Dye, 2 x 1.9 µL RNase-Free Water

For 2500 x 20 µL reactions: 15 x 1.7 mL Quant Nova Probe RT-PCR Master Mix, 5 x 100 µL Quant Nova Probe RT Mix, 5 x 100 µL Internal Control RNA, 5 x 500 µL Yellow Template Dilution buffer, 5 x 1 mL ROX Reference Dye, 10 x 1.9 µL RNase-Free Water

Applications

For highly sensitive and specific one-step qRT-PCR using sequence-specific probes:

- From fresh or frozen adherent, semi-adherent or suspension cells using the supplementary protocol directly from cells
- Also from total RNA or mRNA purified from various starting materials used in gene expression analysis
- Accurate RNA quantification over several logs of template in single or duplex reactions
- two-phase hot-start mechanism keeping both reverse transcriptase and Taq polymerase inactive at room temperature

Required sample amount

The Quant Nova One-Step Probe RT-PCR Kits have been optimized for the following:

- As few as single cells up to 2000 cells using the supplementary protocol directly from cells
- 100 fg to 400 ng of purified template RNA input

To be used with the Quant Nova RT-PCR Kit

- All commonly used real-time cyclers or assays are applicable for use with the Quant Nova kits
- We recommend use of Quant Nova LNA Probe PCR Assays for gene expression analysis, conveniently available from GeneGlobe (www.qiagen.com/QuantNova-LNA)

Process duration

Cells need to be washed, and the desired cell concentration needs to be adjusted before use in One-Step RT-qPCR.

- Transcript verification from washed cells possible within 1.5 hours
- No RNA preparation required
- Cell lysis within 5 minutes at 74°C in PCR vessel
- Integrated reverse transcription step takes 10 minutes at 45°C
- PCR amplification within 45 minutes depending on ramping speed and data acquisition of real time cycler

Ordering webpage: www.qiagen.com/Cell-Line-Solutions

For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at www.qiagen.com or can be requested from QIAGEN Technical Services or your local distributor.

Trademarks: QIAGEN®, Samples to Insights®, Quant Nova® (QIAGEN Group). Registered names, trademarks, etc., used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

05/23 PROM-22183-001 © 2023 QIAGEN, all rights reserved.