

## Droplet Digital™ PCR: One-Step RT-ddPCR Kit for Probes



## One-Step RT-ddPCR Kit for Probes

The Bio-Rad one-step RT-ddPCR kit for probes creates a new paradigm for the precise quantitation of RNA by combining reverse transcription with Droplet Digital™ PCR (ddPCR™). The one-step RT-ddPCR kit for probes provides an absolute measure of target RNA molecules with unrivaled precision and sensitivity.

In combination with the QX100™ Droplet Digital™ PCR system, the one-step RT-ddPCR kit for probes lets you:

- Enrich for rare target RNA sequences
- Detect small differences in gene expression levels
- Determine the number of copies of an RNA molecule without a standard curve

For more information, visit us on the Web at [www.bio-rad.com/one-stepddPCR](http://www.bio-rad.com/one-stepddPCR).

**BIO-RAD**

## A Breakthrough in Quantitative PCR

With the one-step RT-ddPCR kit for probes, the sample is partitioned into 20,000 droplets, with target and background RNA randomly distributed among the droplets. An RNase inhibitor included in the formulation minimizes template degradation during reaction setup and droplet generation. After reverse transcription, the resulting cDNA is amplified for target detection using TaqMan hydrolysis probes. After PCR amplification, each droplet provides a fluorescent positive or negative signal indicating the target RNA was present or not present after partitioning. Each droplet provides an independent digital measurement. Positive and negative droplets are counted and software calculates the concentration of target RNA as copies per microliter.

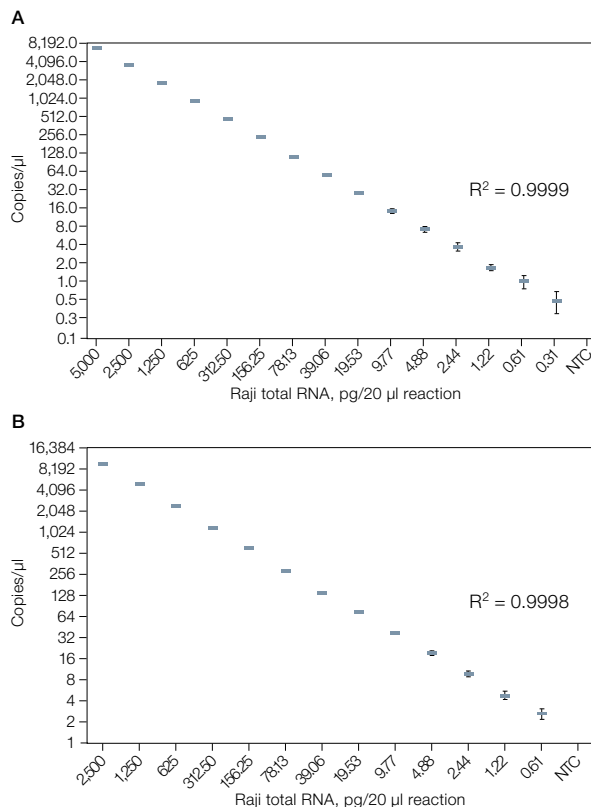
The one-step RT-ddPCR kit for probes is formulated for efficient and sensitive reverse transcription over a wide linear dynamic range of input RNA for Droplet Digital PCR. Our unique hot-start reverse transcriptase enables convenient reaction setup. The reverse transcription reaction is performed at 55–60°C, enhancing the specificity and efficiency of primer-mediated cDNA conversion. The thermostable enzymes allow RNA template to be reverse transcribed and subsequently amplified in the same reaction tube. Template-specific primer annealing at elevated temperatures significantly improves stringency and melting of secondary structures.

## Ordering Information

| Catalog # | Description  |
|-----------|--|
| 186-3021  | <b>One-Step RT-ddPCR Kit for Probes</b> , 2 x 1 ml, 200 x 20 µl reactions, 2x RT-ddPCR mix, contains manganese acetate |
| 186-3022  | <b>One-Step RT-ddPCR Kit for Probes</b> , 5 x 1 ml, 500 x 20 µl reactions, 2x RT-ddPCR mix, contains manganese acetate |

TaqMan is a trademark of Roche Molecular Systems, Inc.

Purchase of this product includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.



**ddPCR absolute quantification of *EEF2* (A) and *GAPDH* (B) gene transcripts using the one-step RT-ddPCR kit for probes.** Raji total RNA was serially diluted twofold from 5 ng to 310 fg per 20 µl reaction. The data demonstrate the precision and sensitivity of the QX100 Droplet Digital PCR system in detecting thousands of copies down to a single copy per microliter. N = 6. NTC, no template control.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000