

# ddPCR Supermix for Residual DNA Quantification

| Catalog # | Description  |
|-----------|--|
| 1864037   | ddPCR Supermix for Residual DNA Quantification, 2 ml (2 x 1 ml vials), 200 x 20 µl reactions     |
| 1864038   | ddPCR Supermix for Residual DNA Quantification, 5 ml (5 x 1 ml vials), 500 x 20 µl reactions     |
| 1864039   | ddPCR Supermix for Residual DNA Quantification, 25 ml (5 x 5 ml vials), 2,500 x 20 µl reactions  |
| 1864040   | ddPCR Supermix for Residual DNA Quantification, 50 ml (10 x 5 ml vials), 5,000 x 20 µl reactions |

For research purposes only.

## Description

ddPCR Supermix for Residual DNA Quantification is a 2x concentrated, ready-to-use reaction cocktail containing all components — except primers, probe(s), and template — required for probe-based Droplet Digital PCR (ddPCR). The mixture delivers maximum target specificity and fluorescence amplitude with minimum droplet variability to ensure precise target quantification for the detection of residual host cell DNA.

The hot-start features of the polymerase enable partitioning of samples into droplets while keeping the enzyme inactive at ambient conditions. The supermix has been optimized to support the amplification and detection of DNA targets using hydrolysis probe-based assays and is compatible with the use of uracil N-glycosylase (UNG) for PCR decontamination. UNG may be purchased from a licensed supplier.

## Storage and Stability

ddPCR Supermix for Residual DNA Quantification is stable at  $-20^{\circ}\text{C}$  through the expiration date printed on the labels. Once thawed, it can be stored at  $4^{\circ}\text{C}$  for up to 2 weeks. Repeated freezing and thawing of the supermix is not recommended.

## Quality Control

ddPCR Supermix for Residual DNA Quantification is free of contaminating DNase and RNase. Stringent specifications are maintained to ensure lot-to-lot consistency.

## Recommendations for Optimal Results

- Follow general guidelines and recommendations for Droplet Digital PCR (refer to the Droplet Digital PCR Applications Guide, bulletin 6407)
- Suggested input quantities of DNA sample are 10 fg–100 ng per 20 µl reaction; add femtogram to picogram amounts of DNA for assays targeting highly repetitive sequences, such as long or short interspersed nuclear elements (LINE or SINE), and picogram to nanogram amounts of DNA for assays targeting single-copy gene sequences
- An assay-specific standard curve is required to convert target copy number to mass concentration for assays targeting an unknown copy number of LINE, SINE, or Alu-like sequences
- A no template control (NTC) should be included to rule out cross-contamination

- Perform DNA extraction for complex samples with high salt ( $\geq 1.0$  M salt) or low pH (buffer  $\leq 3.0$  pH), and perform proteinase K treatment for samples containing a high concentration of protein ( $\geq 0.1$  mg/ml) prior to preparing the ddPCR reaction mix to improve amplification. For less complex samples, pretreatment is not necessary if the sample is diluted at least 25-fold and the protein concentration in the final ddPCR reaction is  $\leq 0.1$  mg/ml

## Required Equipment

The QX200 Droplet Digital PCR System (catalog #1864001), QX200 AutoDG Droplet Digital PCR System (#1864100), or QX ONE Droplet Digital PCR System (#12006536) is required.

Refer to the QX200 Droplet Reader and QuantaSoft Software and QX200 Droplet Generator Instruction Manuals (10031906 and 10031907, respectively), the Automated Droplet Generator Instruction Manual (10043138), or the QX ONE Droplet Digital PCR System Instruction Manual (10000116512) for ordering information about consumables, such as oils, cartridges, gaskets, plates, and seals.

## Reaction Setup

- Thaw all components to room temperature. Mix thoroughly by vortexing the tubes to ensure homogeneity because a concentration gradient may form during  $-20^{\circ}\text{C}$  storage. Centrifuge to collect contents at the bottom of each tube.
- Prepare samples at the desired concentration before setting up the reaction mix.
- Prepare the reaction mix for the number of reactions needed according to the guidelines in Table 1. Assemble all required components except the sample, dispense equal aliquots into each reaction tube, and add the sample to each reaction tube as the final step.

**Table 1. Preparation of the reaction mix.**

| Component   | Volume per Reaction, µl | Final Concentration   |
|---|-------------------------|-----------------------|
| 2x ddPCR Supermix for Residual DNA Quantification | 10                      | 1x                    |
| Target primers/probe                              | Variable                | 900 nM/250 nM         |
| Sample  | Variable                | 10 fg–100 ng/reaction |
| RNase-/DNase-free water                           | Variable                | —                     |
| <b>Total volume*</b>                              | <b>20</b>               | <b>—</b>              |

\* For the Automated Droplet Generator, prepare 22 µl per reaction.

4. Mix thoroughly by vortexing the tubes. Centrifuge briefly to ensure that all components are at the bottom of the reaction tubes.
5. Transfer the reaction mix from the reaction tubes to the appropriate ddPCR Cartridge as follows:
  - For the QX200 Droplet Digital PCR System, load 20 µl of each reaction mix into a sample well of a DG8 Cartridge. Follow subsequent instructions as specified in the QX200 Droplet Generator Instruction Manual (10031907)
  - For the QX200 AutoDG Droplet Digital PCR System, follow instructions in the Automated Droplet Generator Instruction Manual (10043138)
  - For the QX ONE Droplet Digital PCR System, load 20 µl of each reaction mix into a sample well of a GCR96 Cartridge. Follow subsequent instructions as specified in the QX ONE Droplet Digital PCR System Instruction Manual (10000116512)

### Thermal Cycling Conditions

Follow instructions based on the system in use:

- For the QX200 Droplet Digital PCR System, after droplet generation with the QX200 Droplet Generator, carefully transfer droplets into a clean 96-well plate. Seal the plate using the PX1 PCR Plate Sealer at 180°C for 5 sec. Proceed to thermal cycling (see Table 2)
- For the QX200 AutoDG Droplet Digital PCR System, remove the droplet plate containing ddPCR droplets from the Automated Droplet Generator. Seal the plate using the PX1 PCR Plate Sealer at 180°C for 5 sec. Proceed to thermal cycling (see Table 2)
- For the QX ONE Droplet Digital PCR System, thermal cycling is integrated into and sequentially performed by the system itself. Hence, no additional equipment or sample handling is required for this step. Refer to the QX ONE Droplet Digital PCR System Instruction Manual (10000116512) for plate setup instructions. Use appropriate thermal cycling conditions as specified in Table 2

**Table 2. Thermal cycling conditions.\***

| Cycling Step                    |                                | Temperature, °C | Time     | Number of Cycles |
|---------------------------------|--------------------------------|-----------------|----------|------------------|
| Hold (QX ONE ddPCR System only) |                                | 25              | 3 min    | 1                |
| Enzyme activation               |                                | 95              | 10 min   | 1                |
| Denaturation                    |                                | 95              | 30 sec   | 40               |
| Annealing/extension             |                                | 60              | 1 min**  | 40               |
| Enzyme deactivation             |                                | 98              | 10 min   | 1                |
| Hold                            | QX200 ddPCR System (optional)  | 4               | Infinite | 1                |
|                                 | QX ONE ddPCR System (required) | 25              | 1 min    | 1                |

\* For the C1000 Touch Thermal Cycler, use a heated lid set to 105°C and set the sample volume to 40 µl.

\*\* Check/adjust ramp rate settings to ~2°C/sec.

### Data Acquisition and Analysis

1. Follow instructions based on the system in use:
  - For the QX200 Droplet Digital PCR System and the QX200 AutoDG Droplet Digital PCR System, refer to the QX200 Droplet Reader and QuantaSoft Software Instruction Manual (10031906)
  - For the QX ONE Droplet Digital PCR System, refer to the QX ONE Droplet Digital PCR System Instruction Manual (10000116512) and the QX ONE Software Instruction Manual for Standard Edition (10000116655) or Regulatory Edition (10000116656)
2. After data acquisition, select samples in the well selector under Analyze. Examine the automatic thresholding applied to the 1-D amplitude data and, if necessary, set thresholds manually by applying either the single well or multiwell threshold between the positive and negative droplets. Use the NTC well to determine the appropriate threshold for sample wells.
3. The concentration reported is copies/µl of the final 1x ddPCR reaction.

Visit [bio-rad.com/web/ddPCRRDQ](http://bio-rad.com/web/ddPCRRDQ) for more information.

BIO-RAD, DDPCR, DROPLET DIGITAL PCR, and QUANTASOFT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

Purchase of Digital PCR products from Bio-Rad Laboratories includes a limited, non-transferable right under our intellectual property for you to use the products for your internal research purposes only. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the products or to use the products for commercial applications, such as contract services or fee for services. No rights are granted for diagnostic uses. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required. The QX200, QX200 AutoDG, and QX ONE Droplet Digital PCR Systems and/or their use is covered by claims of U.S. patents, and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. See [bio-rad.com/en-us/trademarks](http://bio-rad.com/en-us/trademarks) for details.

All trademarks used herein are the property of their respective owner.