

# Preparing a GCR96 Cartridge for a QX ONE™ Droplet Digital™ PCR System Run

## Quick Start Guide

### Transferring Droplet Digital PCR (ddPCR™) Reactions to a GCR96 Cartridge

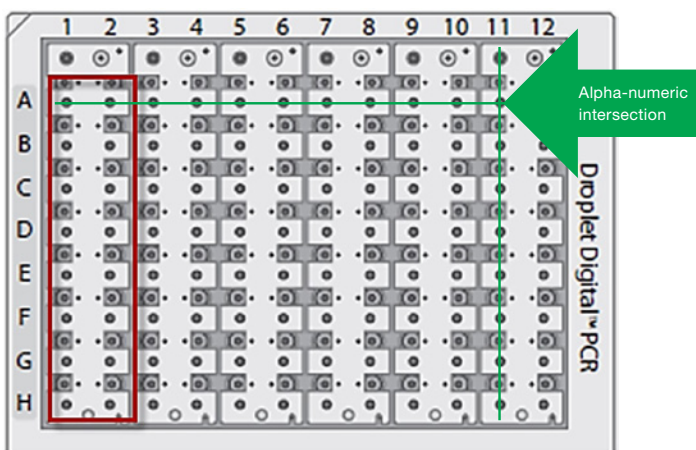
Using a Rainin LTS P20 Multichannel Pipettor, transfer 20 µl of ddPCR reactions, consisting of RNA, DNA, or cDNA template, sample, supermix, assay(s), and nuclease-free water, into the wells of a GCR96 Cartridge that are designated by the labeled alpha-numeric intersections (the wells that align with both a letter and number, as shown). Insert the pipet tips into the wells at approximately a 15° angle, contacting the bottom of the wells. Dispense the ddPCR reactions slowly (approximately 5 sec). Do not push the pipet plunger past the first stop.

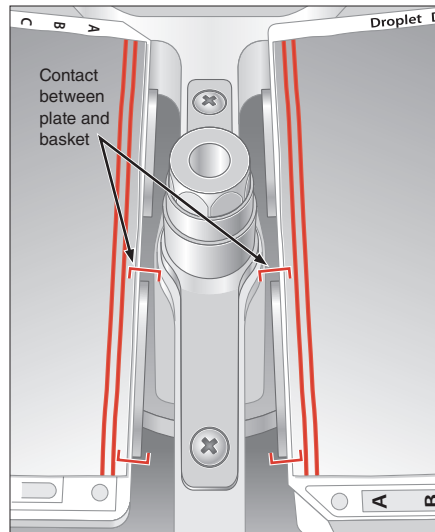
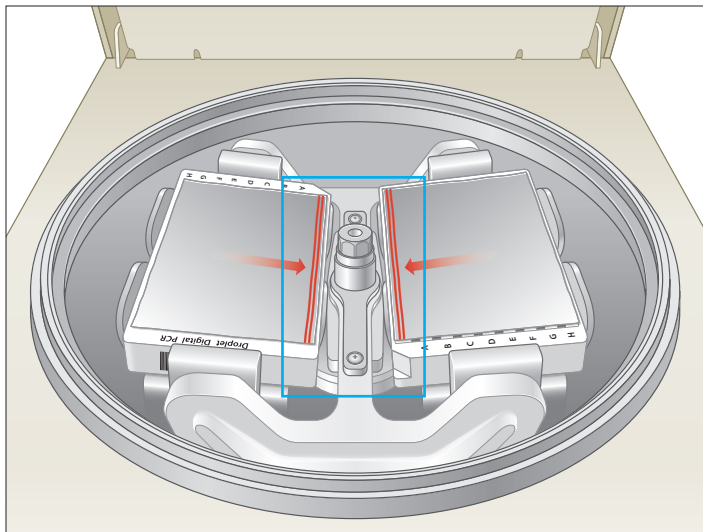
**Note:** Before transferring ddPCR reactions to a GCR96 Cartridge, **thoroughly mix reagents** after thawing them at room temperature (vortex supermix for 30 sec after thawing). Upon generating master mix(es) and after adding the sample/template to the master mix, vortex once more (approximately 15 sec). Centrifuge the ddPCR reactions after vortexing to collect the volume at the bottom of each well (30 sec, 1,150 rcf). Centrifugation eliminates air bubbles and ensures that the ddPCR reactions are at the bottom of each well. **Mixing or vortexing is critical to obtain homogeneity of reagents within droplets.**

**Note:** GCR96 Cartridges contain six strips of 16 wells, as shown. Each strip of 16 is called a DG16 Cartridge. All DG16 Cartridges must be filled with ddPCR reactions or a 1:1 ratio of ddPCR Buffer Control for Probes (Bio-Rad™ Laboratories, Inc., catalog #1863052) and water. Failure to do so will affect the performance of the QX ONE ddPCR System and subsequent experimental data.

#### Sealing a GCR96 Cartridge

- 1 Check the PX1 PCR Plate Sealer to ensure that the correct settings are programmed for sealing the GCR96 Cartridge (Temp.: 180°C; Time: 0.5 sec).
- 2 Securely load the GCR96 Cartridge on the silver heat block (Z-shaped support block; bottom protruding edge of heat block should face the operator). Orient the GCR96 Cartridge so that well A1 is at the upper left corner of the plate.
- 3 Place the pierceable foil heat seal with the double red stripes face up (facing toward the operator) on the GCR96 Cartridge. Ensure the foil seal covers the entire surface area of the cartridge, with the exception of the cartridge labels identifying rows A–H (left-hand side of cartridge) and the Droplet Digital PCR label (right-hand side of cartridge), and that the seal remains flush with the top and bottom of the cartridge (ensure no foil overhangs on any side).





- 4 Select **Eject** from the PX1 PCR Plate Sealer touch screen to eject the drawer. Load the heat block, cartridge, and seal into the PX1 Sealer. Close the drawer by selecting **Close** in the touch-screen menu on the PX1 Sealer (check to make sure seal is still aligned correctly before closing the drawer).
- 5 Select **Seal** (Seal will be green when the PX1 Sealer has reached the sealing temperature of 180°C). The drawer will automatically eject when the seal is completed.
- 6 When the PX1 Sealer ejects the drawer, carefully rotate the semisealed cartridge 180° (do not rotate the heat block from its original orientation) and repeat steps 5–6. Then remove the sealed cartridge and heat block from the PX1 Sealer. Store the heat block on top of the PX1 Sealer until further use and execute steps 7–9 with the sealed cartridge.
- 7 Place the GCR96 Cartridge in a swinging bucket rotor centrifuge, placing row A closest to the rotor spindle (refer to images shown).
- 8 Load a counterbalance cartridge on the opposite rotor bucket.
- 9 Push the individual DG16 Cartridges in the GCR96 frame toward the numbered edge as far as it will go to prevent breaking the tabs of the cartridge. Centrifuge cartridges for 30 sec at 1,150 rcf. Proceed with loading your cartridge into the QX ONE ddPCR System. Refer to Initiating and Managing a QX ONE Droplet Digital PCR System Run Quick Start Guide (bulletin 3431).

Refer to the QX ONE Droplet Digital PCR System and QX ONE Software User Guide (10000116512) for additional details.

Visit [bio-rad.com/QXONE](https://bio-rad.com/QXONE) for more information.

BIO-RAD, DDP, DROPLET DIGITAL, and QX ONE are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2022 Bio-Rad Laboratories, Inc.

The QX ONE Droplet Digital PCR System, and the consumables and reagents designed to work with this system, 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](https://bio-rad.com/en-us/trademarks) for details. Purchase of the product includes a limited, non-transferable right under such intellectual property for use of the product for internal research purposes in the field of digital PCR only. No rights are granted for diagnostic uses. No rights are granted for use of the product for commercial applications of any kind, including but not limited to manufacturing, quality control, or commercial services, such as contract services or fee for services. 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.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Website [bio-rad.com](https://bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399  
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23  
France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050  
Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670  
The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23  
Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23  
Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

