

# ddPCR™ 96-Well Plates

| Catalog # | Description   |
|-----------|---|
| 12001925  | <b>ddPCR™ 96-Well Plates</b> , pkg of 25, clear well/clear shell, semi-skirted plates. For use with QX100™, QX200™, and QX200™ AutoDG™ Droplet Digital™ PCR Systems |

For research purposes only.

## Description

Designed and validated for Bio-Rad's Droplet Digital PCR workflow, ddPCR 96-Well Plates are for use with the QX100, QX200, and AutoDG Droplet Digital PCR Systems. These 96-well semi-skirted PCR plates are optimized for maximum ddPCR performance. Sample wells utilize an ultrathin-wall polypropylene construction for maximal thermal transfer. The robust polycarbonate frame design is highly resistant against warping and torque.

## Compatible Equipment

- Automated Droplet Generator (catalog #1864101)
- QX100 or QX200 Droplet Reader (catalog #1863003 or 1864003, respectively)
- C1000 Touch™ Thermal Cycler with 96–Deep Well Reaction Module (catalog #1851197)\*
- C1000 Touch™ Thermal Cycler with 96-Well Fast Reaction Module (catalog #1851196)
- PX1™ PCR Plate Sealer (catalog #1814000) with Pierceable Foil Heat Seals (catalog #1814040)

\* Recommended for optimal ddPCR performance.

## Quality Control

Each lot of ddPCR 96-Well Plates is tested with Droplet Digital PCR to ensure quality. Strict manufacturing processes ensure that the plates are free of detectable RNase and DNase.

## Key Specifications

|                                   |                                |
|-----------------------------------|--------------------------------|
| Skirt format                      | Semi-skirted                   |
| Well/tube color                   | Clear                          |
| Maximum well volume               | 250 µl                         |
| Maximum centrifuge speed          | 2,250 x g in a swing-out rotor |
| Thermal cycling temperature range | 4–105°C                        |

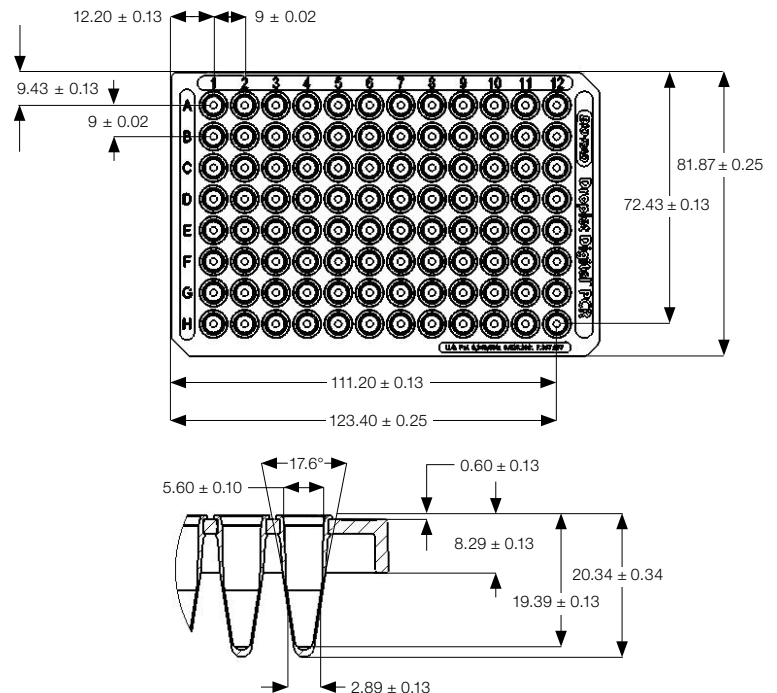
## Sealing Protocol

Set the plate sealer temperature to 180°C and the time to 5 sec. Follow the instructions in the QX100, QX200, or Automated Droplet Generator Instruction Manual (catalog #10026322, 10031907, or 10043138, respectively). For additional operation information, see the PX1 PCR Plate Sealer Instruction Manual (catalog #10023997).

## Plate Dimensions

|  |           |
|--|-----------|
| Length at base plane                       | 123.40 mm |
| Width at base plane                        | 81.87 mm  |
| Height overall                             | 20.68 mm  |
| Well depth                                 | 19.39 mm  |
| Well diameter at opening                   | 5.60 mm   |
| Well diameter at bottom of conical section | 2.89 mm   |
| Well volume                                | 250 µl    |
| Well spacing                               | 9.00 mm   |
| Well angle                                 | 17.6°     |
| Left edge to well A1                       | 12.20 mm  |
| Top edge to well A1                        | 9.43 mm   |
| Left edge to well H12                      | 111.20 mm |
| Top edge to well H12                       | 72.43 mm  |

## Plate Diagram



ddPCR 96-Well Plates are covered by one or more U.S. patents or their foreign counterparts owned by Eppendorf AG.

Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

The QX100, QX200, or QX200 AutoDG Droplet Digital PCR System and/or its 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. 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.