



QX ONE Droplet Digital PCR System



BUILT FOR AUTOMATED, WALK-AWAY DIGITAL PCR OPERATION



DELIVERING ddPCR FOR BIOPHARMA AND CLINICAL RESEARCH

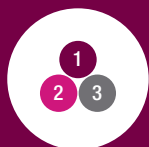
Bio-Rad's unique Droplet Digital PCR (ddPCR) technology has long provided absolute quantification of target DNA and RNA via the generation and thermal cycling of thousands of nanoliter droplets. Now, joining the family alongside the QX200 ddPCR System, the new QX ONE ddPCR System seamlessly integrates the components of a ddPCR workflow — all in one instrument — making it an ideal lab partner for optimizing processes, scaling production, and growing businesses.

FEATURES AND BENEFITS



WALK-AWAY OPERATION

- The QX ONE System automates a fully integrated ddPCR workflow, allowing for walk-away operation
- A 5-plate capacity allows 480 samples to be analyzed in a single day



EASY TO USE

- Simple, intuitive run setup using a large onboard touch screen
- Optimized plate setup means fewer handling steps, fewer operator errors, and faster onboarding



ADVANCED MULTIPLEXING

- Four color channels combined with amplitude multiplexing and/or radial multiplexing allow for multiple unique targets to be detected in a single reaction
- Specially formulated ddPCR Multiplex Supermix is optimized to empower advanced multiplexing in probe-based assays



REGULATORY COMPLIANT

- QX ONE Software, Regulatory Edition is a robust package that enables U.S. FDA 21 CFR Part 11 compliance and offers audit trails with tracked protocol changes for regulatory requirements
- RFID barcoded consumables allow lot management and traceability

WHY BIO-RAD?

Bio-Rad has more than 60 years' experience supporting life science research and biopharma. We are a leader in digital PCR and other analytical technologies. Our innovative products are available as stand-alone solutions or combined into workflows that provide the right answer the first time, every time with minimized downtime.

SPECIFICATIONS

| | |
|-----------------------------|--|
| Dimensions (W x D x H) | 122 cm (48 in) x 66 cm (26 in) x 38 cm (15 in) [76 cm (30 in) to top of monitor] |
| Weight | 100 kg (220 lb) [includes monitor, no reagents] |
| Detection channels | 4 (FAM, HEX, Cy5, and Cy5.5) plus scatter channel |
| Capacity | 5 plates (480 samples) per run |
| Electrical requirements | 100–240 VAC, 50/60 Hz, 10–4.6 A, plugs into standard AC receptacle (similar to the QX200 ddPCR System) Fuse: 10 A, 250 V slow blow |
| Environment | Standard laboratory, 18–30°C, relative humidity 20–80% (noncondensing), altitude between 0 and 2,500 meters (0 to 8,202 feet) |
| Benchtop space requirements | W x D x H: 193 cm (76 in) x 76 cm (30 in) x 126 cm (50 in); 35 cm (14 in) on each side and 10 cm (4 in) behind the machine for proper ventilation Single benchtop rated for more than 114 kg (250 lb) Level within approx. 1° about each axis |
| Touch screen | 61 cm (24 in), Windows 10 Professional 64-Bit, 1 TB SSD hard drive, Intel Core i5-6300U processor, 16 GB RAM, 1,920 x 1,080 display resolution |

| Catalog # | Description |
|---------------------|---|
| Reagents | |
| 1863023 | ddPCR Supermix for Probes (No dUTP) , 2 ml (2 x 1 ml), 200 x 20 µl reactions |
| 1863024 | ddPCR Supermix for Probes (No dUTP) , 5 ml (5 x 1 ml), 500 x 20 µl reactions |
| 1863025 | ddPCR Supermix for Probes (No dUTP) , 25 ml (5 x 5 ml), 2,500 x 20 µl reactions |
| 1863026 | ddPCR Supermix for Probes , 2 ml (2 x 1 ml), 200 x 20 µl reactions |
| 1863010 | ddPCR Supermix for Probes , 5 ml (5 x 1 ml), 500 x 20 µl reactions |
| 1863027 | ddPCR Supermix for Probes , 25 ml (5 x 5 ml), 2,500 x 20 µl reactions |
| 1863028 | ddPCR Supermix for Probes , 50 ml (10 x 5 ml), 5,000 x 20 µl reactions |
| 12005909 | ddPCR Multiplex Supermix , 1 ml (2 x 0.5 ml), 200 x 20 µl reactions |
| 12005910 | ddPCR Multiplex Supermix , 2.5 ml (5 x 0.5 ml), 500 x 20 µl reactions |
| 12005911 | ddPCR Multiplex Supermix , 12.5 ml (5 x 2.5 ml), 2,500 x 20 µl reactions |
| 1864037 | ddPCR Supermix for Residual DNA Quantification , 2 ml (2 x 1 ml), 200 x 20 µl reactions |
| 1864038 | ddPCR Supermix for Residual DNA Quantification , 5 ml (5 x 1 ml), 500 x 20 µl reactions |
| 1864039 | ddPCR Supermix for Residual DNA Quantification , 25 ml (5 x 5 ml), 2,500 x 20 µl reactions |
| 1864040 | ddPCR Supermix for Residual DNA Quantification , 50 ml (10 x 5 ml), 5,000 x 20 µl reactions |
| 1864021 | One-Step RT-ddPCR Advanced Kit for Probes , 200 x 20 µl reactions |
| 1864022 | One-Step RT-ddPCR Advanced Kit for Probes , 500 x 20 µl reactions |
| 1863052 | ddPCR Buffer Control for Probes , 9 ml (2 x 4.5 ml) |
| Plate Sealer | |
| 1814000 | PX1 PCR Plate Sealer , includes heat sealing instrument, plate support block that holds 96-well and 384-well plates, sealing frame, power cord |

ORDERING INFORMATION

| Catalog # | Description |
|--|--|
| QX ONE Droplet Digital PCR (ddPCR) System | |
| 12006536 | QX ONE Droplet Digital PCR System , integrated system used for sequential droplet generation, thermal cycling, and droplet reading; includes instrument, touch screen, and QX ONE Accessory Kit |
| — | QX ONE Software (Standard Edition) , connects to the QX ONE Droplet Digital PCR System for instrument control, assay setup, and data collection and analysis |
| 12012078 | QX ONE Software (Regulatory Edition) , connects to the QX ONE Droplet Digital PCR System for instrument control, assay setup, and data collection and analysis; offers audit trails |
| 17005930 | QX ONE Droplet Digital PCR System with PX1 PCR Plate Sealer , includes QX ONE Droplet Digital PCR instrument, touch screen, QX ONE Accessory Kit, and PX1 PCR Plate Sealer |

Consumables

| | |
|-------------|--|
| 12006859 | GCR96 Cartridges , pkg of 10 |
| 12006858 | GCR96 Cartridge , pkg of 1 |
| 12006843 | GCR96 Foil Heat Seal , pkg of 100 |
| 12006060 | QX ONE ddPCR System Waste Bottle |
| Oils | |
| 12006058 | QX ONE Droplet Generation Oil for Probes , enough for 18 x 96-well plates |
| 12006057 | QX ONE Droplet Reader Oil , enough for 10 x 96-well plates |

Visit bio-rad.com/QXONE for more information.

BIO-RAD, DDPCR, DROPLET DIGITAL PCR, and QX ONE are trademarks of Bio-Rad Laboratories 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 QX ONE and QX200 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 for details.

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

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

