



## QX ONE Droplet Digital PCR (ddPCR) System

The QX ONE ddPCR System is designed to deliver accurate, precise, and multiplex digital PCR results. This system seamlessly integrates a standard ddPCR workflow of droplet generation, thermal cycling, droplet reading, and analysis into an automated precision platform.

The QX ONE offers users:

- Walk-away operation (5-plate capacity)
- Advanced multiplexing capabilities
- A new multiplex supermix and compatibility with existing supermixes
- Easy-to-use QX ONE Software

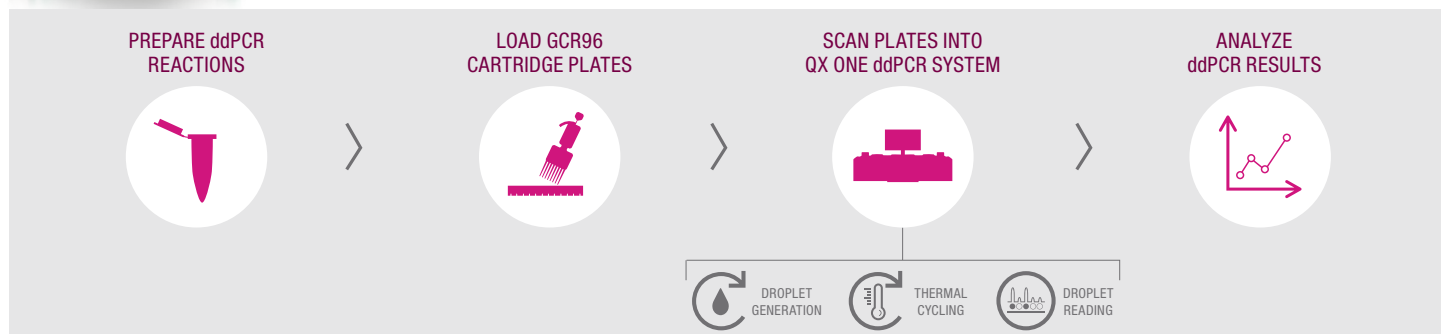
Visit [bio-rad.com/QXONE](http://bio-rad.com/QXONE) for more information.




## THE WORKFLOW




The QX ONE ddPCR System combines partitioning, amplification, and droplet reading into a single, completely automated process. Once a protocol is selected, plates are scanned and loaded into the QX ONE ddPCR System.




## 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

## Ordering Information

Catalog #	Description
12006536	<b>QX ONE Droplet Digital PCR System</b>

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

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 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. See [bio-rad.com/en-us/trademarks](http://bio-rad.com/en-us/trademarks) for details.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](http://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

