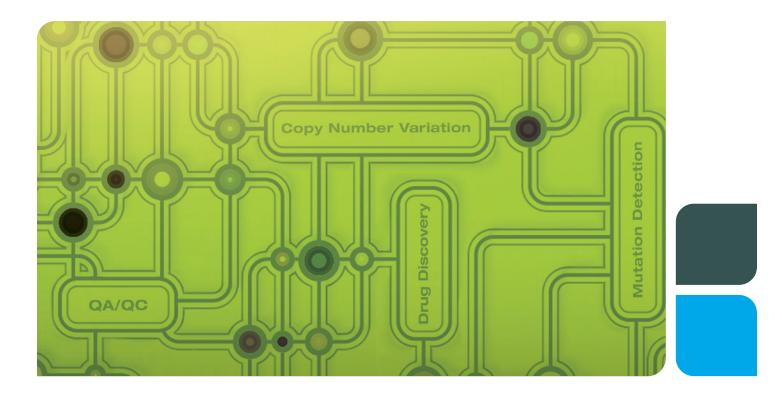
Droplet Digital PCR: QX200 AutoDG Droplet Digital PCR System



Automated Droplet Generator

In a Complicated World, the Future Just Got Simpler

Bio-Rad's QX200 Droplet Digital PCR (ddPCR) System gives scientists the power to unveil new discoveries through precise and absolute nucleic acid quantification. The Automated Droplet Generator (AutoDG) simplifies the ddPCR workflow, making digital PCR both scalable and practical.

Simplifies the ddPCR Workflow and Minimizes Hands-On Time

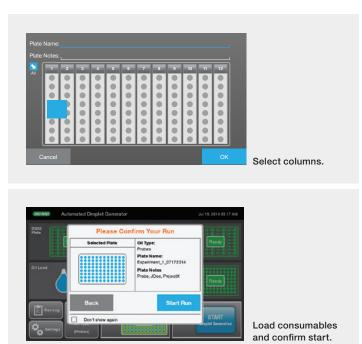
- Generates droplets for 96 ddPCR reactions in less than 45 minutes
- Guides setup in less than 5 minutes with large, color touch-screen interface and light-emitting diode (LED) lighting
- Requires no supervision or hands-on activity during droplet generation

Eliminates User-to-User Variability and Offers Flexibility

- More consistent droplet counts across users and plates
- User-definable plate layout with increments of 8 wells
- A single AutoDG can keep multiple QX200 Droplet Readers running
- HEPA-filtered enclosure reduces contamination and allows the system to be placed in a standard laboratory









Countdown timer indicates time remaining until the plate is ready.



time elapsed since droplet generation.

Simply Set Up and Go

The Automated Droplet Generator's large, color touch-screen interface guides users to droplet generation in just three easy steps.

- Select desired columns across a 96-well plate
- Load only the consumables you need, as prompted by the system
- Confirm and start the run

A countdown timer will display the time remaining so users can maximize their time in the laboratory. When the plate is ready, users are notified and a time-elapsed counter begins. Every run is stored in an exportable log file for future reference.

Bio-Rad's QX200 AutoDG Droplet Digital PCR System combines workflow simplification with established leadership in absolute quantification.

Ordering Information

Catalog #	Description
1864100	QX200 AutoDG Droplet Digital PCR System, includes Automated Droplet Generator, QX200 Droplet Reader, laptop computer
1864101	Automated Droplet Generator
1864108	DG32 Automated Droplet Generator Cartridges, pkg of 30, enough for 10 x 96-well ddPCR Plates
1864109	DG32 Automated Droplet Generator Cartridges, pkg of 60, enough for 20 x 96-well ddPCR Plates
1864110	Automated Droplet Generation Oil for Probes, 140 ml, enough for 20 x 96-well ddPCR Plates
1864112	Automated Droplet Generation Oil for EvaGreen [®] , 140 ml, enough for 20 x 96-well ddPCR Plates
1864120	Pipet Tips for AutoDG System , pkg of 20, enough for 10 x 96-well ddPCR Plates
1864121	Pipet Tips for AutoDG System , pkg of 40, enough for 20 x 96-well ddPCR Plates
1864125	Pipet Tip Waste Bins for AutoDG System , pkg of 10, enough for 10 x 96-well ddPCR Plates

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

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

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen Dye for use in real-time PCR, for research purposes only.

Purchase of Digital PCR products from Bio-Rad Laboratories includes a limited, nontransferable 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 and QX200 AutoDG 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/trademark for details.

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



Bio-Rad Laboratories, Inc.

Life Science Group

Website 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
 Website
 bio-rad.com
 USA
 1800
 424
 6723
 Austral 00 800
 00 24
 6723
 Beigrum 100 800
 00 24
 6723
 Finand 100 800
 02 46
 723
 Finand 100 800
 02 46

23