

Automated Droplet Generator

In a Complicated World, the Future Just Got Simpler

Bio-Rad's QX100™ and QX200 Droplet Digital PCR (ddPCR™) Systems gave scientists the power to unveil new discoveries through precise and absolute nucleic acid quantification. The new Automated Droplet Generator (AutoDG) simplifies the ddPCR workflow, making digital PCR both scalable and practical.

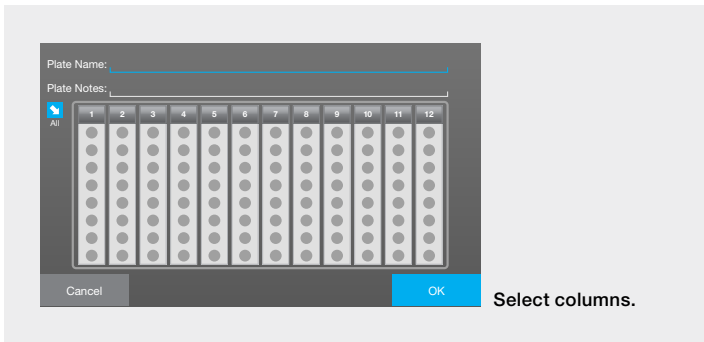
Simplifies the ddPCR Workflow and Minimizes Hands-On Time

- 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

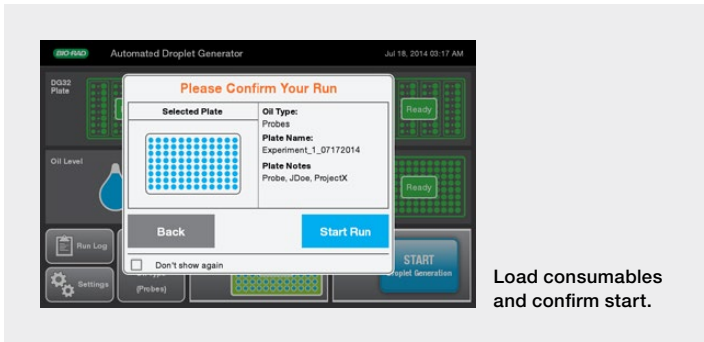
Eliminates User-to-User Variability and Offers Flexibility

- More consistent droplet counts across users and plates
- User-definable plate layout with increments of 8 wells
- A single AutoDG can keep multiple QX100 or QX200 Droplet Readers running
- HEPA-filtered enclosure reduces contamination and allows the system to be placed in a standard laboratory

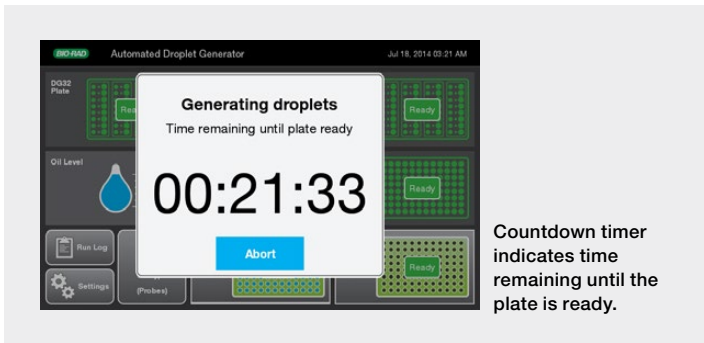




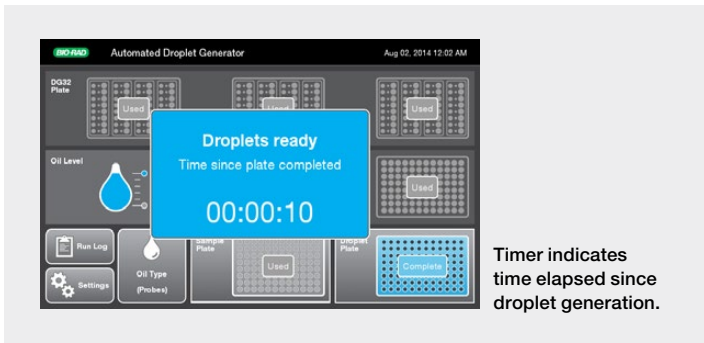
Select columns.



Load consumables and confirm start.



Countdown timer indicates time remaining until the plate is ready.



Timer indicates 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
186-4100	QX200 AutoDG Droplet Digital PCR System , includes Automated Droplet Generator, QX200 Droplet Reader, laptop computer
186-4101	Automated Droplet Generator
186-4108	DG32™ Automated Droplet Generator Cartridges , pkg of 30, enough for 10 x 96-well ddPCR plates
186-4109	DG32 Automated Droplet Generator Cartridges , pkg of 60, enough for 20 x 96-well ddPCR plates
186-4110	Automated Droplet Generation Oil for Probes , 140 ml, enough for 20 x 96-well ddPCR plates
186-4112	Automated Droplet Generation Oil for EvaGreen , 140 ml, enough for 20 x 96-well ddPCR plates
186-4120	Pipet Tips for AutoDG System , pkg of 20, enough for 10 x 96-well ddPCR plates
186-4121	Pipet Tips for AutoDG System , pkg of 40, enough for 20 x 96-well ddPCR plates
186-4125	Pipet Tip Waste Bins for AutoDG System , pkg of 10, enough for 10 x 96-well ddPCR plates

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

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.

The QX100 or QX200 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 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

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000