

QX700 HT Droplet Digital PCR System

Bio-Rad's Simplest High-Throughput Droplet Digital PCR (ddPCR) Powerhouse

When Droplet Digital™ PCR (ddPCR™) precision is required for your high-throughput research applications, the QX700™ HT System* is ready to deliver. QX700 HT System integrates the complete ddPCR workflow into a single instrument, capable of processing 384 samples every 8 hours — allowing your lab to scale analysis with confidence and reliability. Dynamically load samples at any time during the workday to maximize throughput and operational flexibility.



*For Research Use Only. Not for use in diagnostic procedures.

Features and Benefits

Simple Workflow

Droplet generation, thermal cycling, and analysis — all in one instrument. Run everything from a large touchscreen with intuitive, plug-and-play software

High Throughput

Seamless scale from bench-top experiments to production volumes. Process 96 samples in under 2.5 hours and up to 384 samples every 8 hours

Dynamic Loading

Queue up to 384 samples at once or load on demand with random-access capability. Independent thermal control for each 16-well cartridge enables true sample scheduling flexibility

Flexible Options for Dynamic or Full Instrument Loading

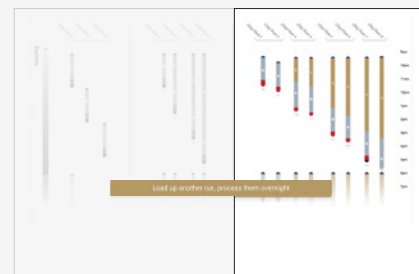
Random Access — Great for Core Facility, Production Scale Testing



Scalable



Flexible Options for Dynamic or Full Instrument Loading



Product Specifications

Channels	7
Input volume	5 µL
Partitions	~17,000
Automation	Full
Throughput (RDG48 plates)	8
Throughput (# wells)	384
Time to full result	8 hours

Visit [bio-rad.com/QX700HT](https://www.bio-rad.com/QX700HT) for more information.

Bio-Rad, QX700, ddPCR, Droplet Digital, and Droplet Digital PCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

© 2025 Bio-Rad Laboratories, Inc.

Purchase of Digital PCR and/or Single-Cell NGS Sample Preparation products (the "Products") from Bio-Rad Laboratories is subject to the Bio-Rad Laboratories, Inc. Standard Terms and Conditions of Sale, which can be accessed at <https://www.bio-rad.com/en-us/terms-conditions>. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the Products. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for the development or commercialization of diagnostic assays for use with the Products without a license from Bio-Rad. It is the user's obligation to obtain a commercial license from Bio-Rad for (i) all commercial uses (not just diagnostic uses) and (ii) sale of assays for use on Bio-Rad's dPCR and ddSEQ instruments. The Products and/or their use are covered by U.S. and foreign patents and/or pending patent applications owned by or under license to Bio-Rad Laboratories, Inc. See <https://www.bio-rad.com/en-us/trademarks>.

*For Research Use Only. Not for use in diagnostic procedures.

Ordering Information

Catalog #	Description
17010628	QX700 HT Droplet Digital PCR System
12025252	RDG16 Cartridges, Pack of 12
12025262	RDG16 Cartridge Holder
12025253	naica® 5X Multiplex ddPCR Mix, 1 × 0.75 mL
12025254	naica® 5X Multiplex ddPCR Mix, 2 × 0.75 mL
12025257	naica® 5X ddPCR Mix, 2 × 0.75 mL
12025255	naica® 10X Multiplex ddPCR Mix, 1 × 0.375 mL
12025256	naica® 10X Multiplex ddPCR Mix, 2 × 0.375 mL
12025258	naica® 10X ddPCR Mix, 2 × 0.375 mL



**Bio-Rad
Laboratories, Inc.**

Life Science Group

Website [bio-rad.com](https://www.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 **Canada** 1 800 361 1808 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23