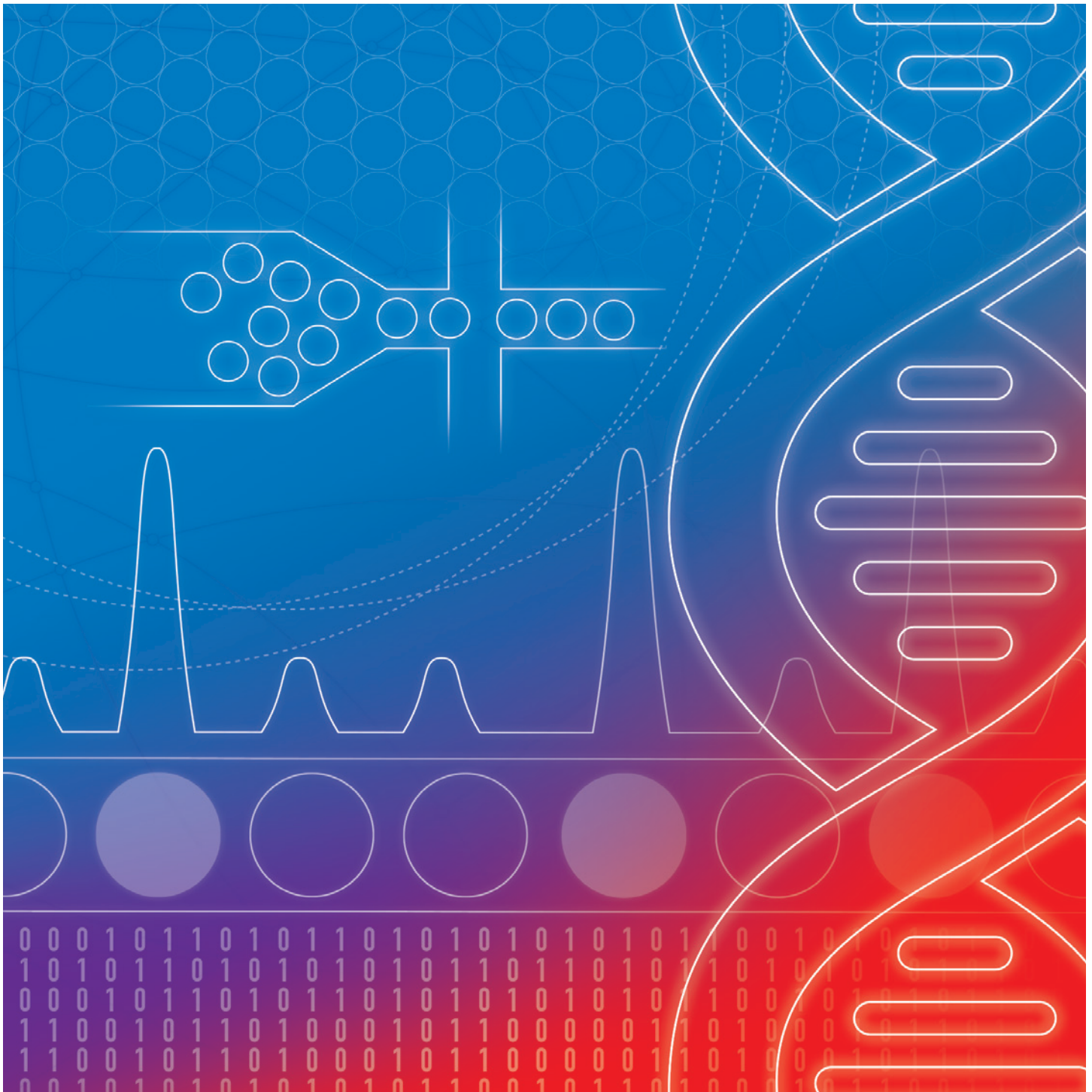


QX600™ Droplet Digital™ PCR System



MULTIPLEXING SIMPLIFIED — GET MORE DATA FROM YOUR PRECIOUS SAMPLES

The QX600 Droplet Digital PCR (ddPCR™) System provides advanced, six-color multiplexing to maximize the number of biomarkers interrogated per well, allowing the preservation of precious samples and reducing laboratory burden. For maximum flexibility, the QX200™ Droplet Digital PCR System assays, consumables, and reagents can be used on the QX600 ddPCR System.



FEATURES AND BENEFITS



SENSITIVE MULTIPLEXING

- 6-color detection capability (FAM/EvaGreen®, HEX/VIC, Cy5, Cy5.5, ROX, and ATTO 590)
- Quantification of 12 targets in a single well
- Absolute quantification with 0.1% or better sensitivity



ACCURATE AND REPRODUCIBLE RESULTS

- Confidence in your results — the system has been qualified by traceable material from the National Measurement Institute of Australia (NMIA)
- Reproducible results — market-leading ddPCR technology has been cited in reproducibility studies



PORTFOLIO COMPATIBILITY

- Same assays, reagents, and consumables as QX200 Droplet Digital PCR System
- Extensive menu of existing ddPCR Assays for EvaGreen® and hydrolysis probe chemistry

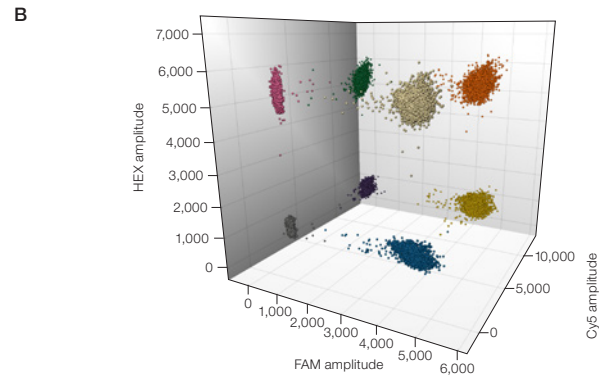
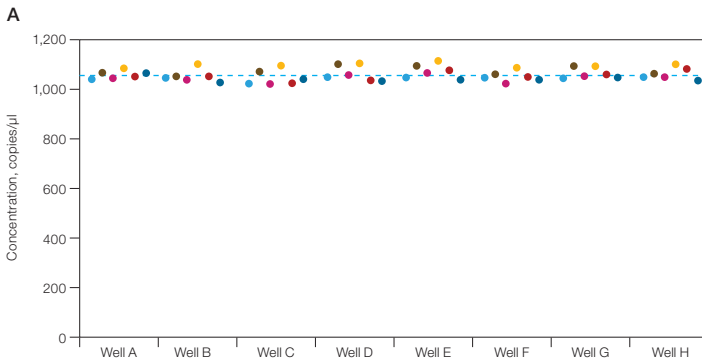


POWERFUL SOFTWARE

- Results that are easy to analyze
- Premium edition available featuring audit trails and tools for U.S. FDA 21 CFR Part 11 compliance
- Premium edition supports assay protocol files for automated data analysis and reporting

SIX-COLOR DETECTION CAPABILITY WITH HIGH ACCURACY, PRECISION, AND REPRODUCIBILITY

The QX600 ddPCR System offers six-color multiplexing for the clear discrimination of multiple targets and more accurate and reproducible quantification with six-color detection capability (FAM/EvaGreen®, HEX/VIC, Cy5, Cy5.5, ROX, and ATTO 590). It has been qualified by traceable material from the National Measurement Institute of Australia.



Metric	QX600 ddPCR System, %	Other Digital PCR Platform, %
Accuracy, difference from NMIA target concentration	0.7	3.2 ^{**}
Precision, % CV, mean concentration across a plate	1.7	3.5–3.8 ^{***}

^{*} NMIA standard, data shown, 2023.

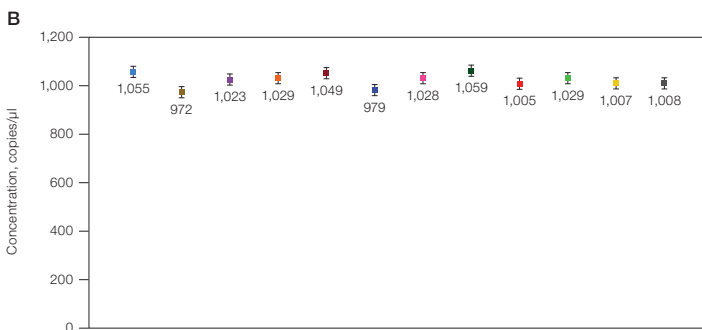
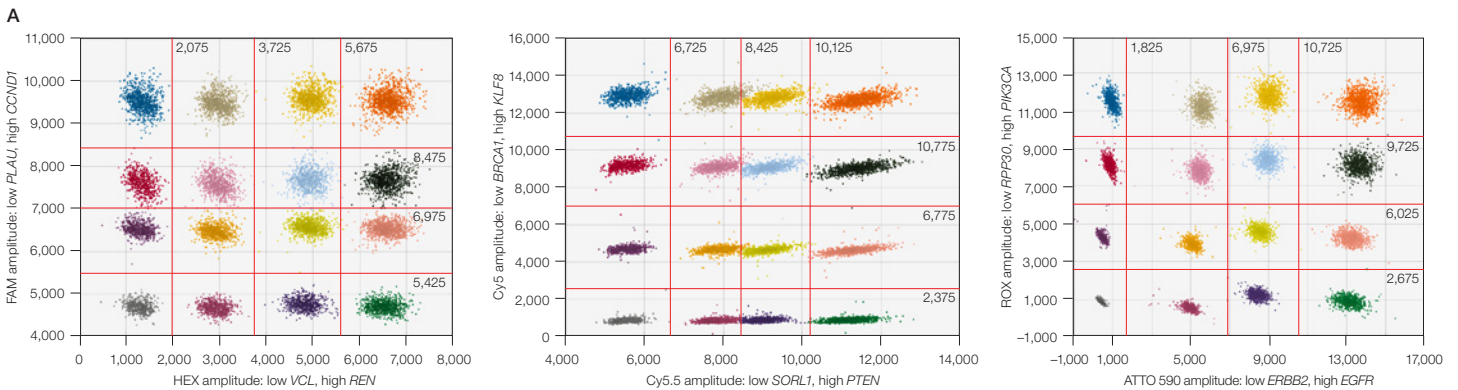
^{**} NIST standard, as published by competitor, 2023.

^{***} *ERBB2* assay, as published by competitor, 2022.

Accurate, precise, and reproducible data of a known metrology standard for six targets. **A**, concentration plot shows accurate, precise, and reproducible values across eight wells of the 6-plex assay. NMIA standard target concentration is 1,060 copies/μl (---). FAM (■), HEX (■), Cy5 (■), Cy5.5 (■), ROX (■), ATTO 590 (■). **B**, 3-D fluorescence amplitude plots of a 6-plex NMIA standard. **C**, the concentration calls for the NMIA standard were accurate and precise on the QX600 Droplet Reader, with no greater than 1% difference compared to the NMIA standard concentration and with CV <2.0%. Data were generated on a QX600 Droplet Reader using ddPCR Multiplex Supermix and a Bio-Rad performance qualification test. CV, coefficient of variation; NMIA, National Measurement Institute of Australia.

QUANTIFICATION OF 12 TARGETS IN A SINGLE WELL

The QX600 ddPCR System provides unprecedented multiplexing capability for 12 copy number targets in a single well using amplitude multiplexing.



Detection of 12 copy number targets in a single well. **A**, 2-D fluorescence amplitude plots show a single well with 12 targets across six fluorescence channels using amplitude multiplexing; **B**, concentration plot shows measured concentration of 12 genes (*PLAU*, *CCND1*, *VCL*, *REN*, *BRCA1*, *KLF8*, *SORL1*, *PTEN*, *RPP30*, *PIK3CA*, *ERBB2*, and *EGFR*) for which the expected concentration was 1,000 copies/μl. Data were generated on a QX600 Droplet Reader using ddPCR Multiplex Supermix and ddPCR Copy Number Assays (available at [bio-rad.com/digital-assays](https://www.bio-rad.com/digital-assays)).

WHY BIO-RAD?

Bio-Rad has more than 70 years' experience supporting life science research and biopharma. We are a leader in digital PCR and other 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.

KEY APPLICATIONS

The QX600 ddPCR System's six-color detection capabilities set a new standard in user friendliness when developing highly multiplexed assays in the following key applications.



Mutation Detection

The QX600 ddPCR System enables six-color mutation detection with ultrasensitive, cost-effective circulating tumor DNA analysis and a rapid turnaround time — unlike next-generation sequencing and other digital PCR platforms.



Viral Titer and Vector Identity

The QX600 ddPCR System enables rapid viral vector titer characterization of multiple critical targets like inverted terminal repeat, target gene of interest, control elements, and terminator, all in the same well, saving time and cost.



Wastewater Surveillance

The QX600 ddPCR System can quantify multiple pathogens in wastewater with high precision and sensitivity.

DIGITAL ASSAYS

Bio-Rad has a large menu of off-the-shelf assays, kits, and design services. The digital assays website allows ddPCR users to find and buy assays across a variety of applications. You can search curated lists of predesigned and validated assays, or design and customize your own assay by using one of the assay design engines. Built by ddPCR experts, the design engines customize a ready-to-use assay optimized for Droplet Digital PCR.

Go to [bio-rad.com/digital-assays](https://www.bio-rad.com/digital-assays) to find and buy ddPCR Assays.

SPECIFICATIONS

Parameter	QX600 Droplet Reader	Automated Droplet Generator	QX200 Droplet Generator
Dimensions (W x D x H)	66 x 52 x 29 cm (26 x 20 x 11 in.)	66 x 56 x 66 cm (26 x 22 x 26 in.)	28 x 36 x 13 cm (11 x 14 x 5 in.)
Weight	26 kg (56.6 lb)	45.5 kg (100 lb)	4.5 kg (10 lb)
Capacity	1–96 samples/run	1–96 samples/run	1–8 samples/run
Droplets per 20 µl sample	20,000	20,000	20,000
Detection channels	FAM/EvaGreen®, HEX/VIC, Cy5, Cy5.5, ROX, and ATTO 590	—	—
Sample Illumination	Light-emitting diodes	—	—
Sample detection	Avalanche photodiodes	—	—

ORDERING INFORMATION

Catalog # Description

QX600 Droplet Digital PCR Systems

17007769	QX600 Droplet Digital PCR System
17008371	QX600 AutoDG™ Droplet Digital PCR System
12013328	QX600 Droplet Reader
1864002	QX200 Droplet Generator
1864101	Automated Droplet Generator

Visit [bio-rad.com/QX600](https://www.bio-rad.com/QX600) for more information.

BIO-RAD, AUTODG, DDPCR, DROPLET DIGITAL, DROPLET DIGITAL PCR, QX200, and QX600 are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.

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 and/or Single-Cell NGS Sample Preparation products (the "Products") from Bio-Rad Laboratories is subject to 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>.

BIO-RAD

**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 1 905 364 3435 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

