



QX600 Droplet Reader

The QX600™ Droplet Reader enables six-color detection for multiplexing, allowing the clear discrimination of multiple targets with assays that can port over and grow from the QX200™ Droplet Digital™ PCR (ddPCR™) System. The QX600 Droplet Reader is designed for researchers who need to quantify multiple targets with high accuracy, reproducibility, and sensitivity.

The QX600 Droplet Reader offers users:

- Sensitive multiplexing
- Accurate and reproducible results
- Portfolio compatibility
- Powerful software



HIGHER MULTIPLEXING, UPGRADED PERFORMANCE



The QX600 Droplet Reader is a new six-color Droplet Digital PCR reader that delivers the same best-in-class performance you've come to expect from Bio-Rad Droplet Digital PCR Systems, with the addition of more channels for multiplexing. The QX600 Droplet Reader has a wide variety of applications, including rare mutation detection, gene expression analysis, genome edit detection, copy number variation studies, multiplex mutation screening, viral titer measurement, contamination testing, and research into infectious diseases such as coronavirus disease 2019 (COVID-19). For maximum flexibility, QX200 System assays, consumables, and reagents can be used on the QX600 Droplet Digital PCR System.

FEATURES AND BENEFITS

SENSITIVE
MULTIPLEXING

- 6-color detection capability (FAM, HEX, Cy5, Cy5.5, ROX, and ATTO 590)
- Quantification of 12 targets in a single well with 6-color detection
- Absolute quantification with 0.1% or better sensitivity
- Sensitive and precise gene expression multiplexing

ACCURATE AND
REPRODUCIBLE
RESULTS

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

PORTFOLIO
COMPATIBILITY

- Same reagents and consumables as QX200 System
- Extensive menu of existing ddPCR assays

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

Ordering Information

Catalog #	Description
12013328	QX600 Droplet Reader

BIO-RAD, DDPCR, DROPLET DIGITAL, QX200, and QX600 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 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. 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
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
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 2 3473 4460 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

