



QXDx™ Universal Kit for AutoDG™ ddPCR™ System

Droplet Digital™ PCR (ddPCR) provides key advantages, including precision, absolute quantification, high sensitivity, and low cost, making it the perfect choice when quick and iterative decisions are essential. The launch of the QXDx Universal Kit for AutoDG ddPCR System with GMP manufactured general purpose reagents now allows for a more streamlined workflow and even greater reproducibility.

The kit includes all of the non-analyte-specific reagents and consumables that are required for the QXDx AutoDG ddPCR System. It allows use of the QXDx AutoDG ddPCR System in Open Mode and is designed and packaged to:

- Simplify Droplet Digital PCR by eliminating the need to determine which consumables are required
- Provide a validated kit with verified performance standards manufactured under GMP
- Deliver a complete digital PCR solution that requires design of only primers and probe(s)
- Achieve high assay precision and accuracy and consistent results through absolute quantification

Visit [bio-rad.com/web/CE-IVD/AutoDGUniversalKit](https://www.bio-rad.com/web/CE-IVD/AutoDGUniversalKit) for more information.



The Complete Workflow

To simplify digital PCR we provide a complete workflow, from reagents and consumables to instruments and analysis software.



The QXDx Universal Kit for AutoDG ddPCR System

The QXDx Universal Kit for AutoDG ddPCR System consists of the following:

- QXDx AutoDG Supermix Pack
- QXDx AutoDG Consumable Pack
- QXDx Droplet Reader Oil Pack

The only reagents not included are the primers and probe(s) that are unique to each assay.

Ordering Information

Catalog #	Description
17001378*	QXDx Universal Kit for AutoDG ddPCR System , 480 reaction kit for QXDx AutoDG ddPCR System
12003031*	QXDx AutoDG Supermix Pack , 480 reaction Universal Supermix Pack
12001922*	QXDx AutoDG Consumable Pack , 480 reaction consumable pack consisting of 96-well plates (3 x 5 pack); Foil Heat Seals (1 x 50); Droplet Generation Oil (1 x 50 ml); DG32™ Cartridges (1 x 15 pack); AutoDG Pipette Tips (1 x 10 pack)
12002526*	QXDx Droplet Reader Oil Pack , 1 L of ready-to-use Droplet Reader Oil for QXDx AutoDG ddPCR System
Material not included in the kit:	
17002229*	QXDx AutoDG ddPCR System , QXDx Automated Droplet Generator; QXDx Droplet Reader
1814000	PX1 PCR Plate Sealer
—	Thermal cycler , with the following specifications: <ul style="list-style-type: none">– Accuracy: ± 0.2°C– Uniformity: ± 0.4°C well-to-well within 10 sec– Adjustable ramp capability with required ramp rate: up to 2°C/sec– Temperature range: 0–100°C

* These products are CE Marked IVDs. They have not been submitted to U.S. FDA and are not available in the U.S. market.

Bio-Rad is in the process of updating IVD product names to include QXDx. Visit bio-rad.com/web/QXDxNameUpdate for more detailed information about product name updates. Current product is labeled as QX200.

Droplet Digital PCR products 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. Purchase of the product includes a limited, non-transferable right under such intellectual property for use of the product for internal research and diagnostic purposes only. 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.

The C1000 Touch Thermal Cycler is covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00
France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 972 03 963 6050
Italy 39 02 216091 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 (0)318 540 666 New Zealand 64 9 415 2280
Norway 47 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 (0) 861 246 723
Spain 34 91 590 5200 Sweden 46 08 555 12700 Switzerland 41 026 674 55 05 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311
United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

