

ddPCR Expert Design Assays

| Catalog # | Description |
|--|---|
| 12008183, 12008184, 12008185, 12008186, 12008187, 12008188, 12008189, 12008190, 12008201, 12008202, 12008203, 12008204, 12008205, 12008206, 12008207, 12008208, 12008209, 12008210, 12008211, 12008212, 12008213, 12008214, 12008215, 12008216, 12008217, 12008218, 12008219 | ddPCR Expert Design Assay, 200 x 20 µL reactions |
| 12003898, 12003899, 12003900, 12003901, 12003902, 12003903, 12003904, 12003905, 12003906, 12003907, 12003908, 12003909, 12003910, 12003923, 12004111, 12004112, 12004113, 12004114, 12004115, 12004116, 12004117, 12004118, 12004119, 12004120, 12004121, 12004122, 12004123 | ddPCR Expert Design Assay, 1000 x 20 µL reactions |

For research purposes only.

Description

ddPCR Expert Design Assays are designed specifically for Droplet Digital™ PCR (ddPCR) applications by the experts who developed the technology. The assays can be used for a wide range of experiments, including translocations, mutation detection, multiplex mutation screening, and more. Each one has a unique assay ID that corresponds to a specific formulation of primers and probes. These assays should be validated prior to use by the end user.

Ordering Information

The ddPCR assays can be ordered online at bio-rad.com/digital-assays. Required ddPCR Supermixes can be also be ordered at bio-rad.com/web/ddPCRmultiplexsmx.

Storage and Stability

The ddPCR assays are stable for 12 months when stored at 4°C in the dark. The assay can be kept at -20°C for longterm storage.

Kit Contents

The ddPCR Expert Design Assay is a concentrated, ready-to-use primer-probe mix. Each tube has either 200 or 1,000 µL of the selected assay.

See Table 1 for a list of fluorophores compatible with your system.

Visit bio-rad.com/ddPCRconsumables to find ordering information on consumables such as oils, cartridges, gaskets, plates, and seals.

Table 1. Fluorophore compatibility.

| QX200 Droplet Reader | QX ONE ddPCR System | QX Continuum ddPCR System | QX600 Droplet Reader | QX700 ddPCR System |
|----------------------|---------------------|---------------------------|----------------------|--------------------|
| FAM | FAM | FAM | FAM | FAM |
| HEX | HEX | HEX | HEX | HEX |
| | Cy5 | ROX | ROX | ATTO 550 |
| | Cy5.5 | Cy5 | ATTO 590 | ROX |
| | | | Cy5 | ATTO 590 |
| | | | Cy5.5 | Cy5 |
| | | | | Cy5.5 |

Required Reagents and Equipment

For assays using the QX200™ Droplet Generator (catalog #1864002) or Automated Droplet Generator (AutoDG™, #1864101):

- iScript Advanced cDNA Synthesis Kit (#1725037, 1725038)
- For 1–2 targets, ddPCR Supermix for Probes (No dUTP) (#1863023, 1863024, 1863025)
- For >2 targets, ddPCR Multiplex Supermix (#12005909, 12005910, 12005911)
- QX100, QX200, or QX600™ Droplet Reader (#1863003, 1864003, or 12013328, respectively)
- PX1 PCR Plate Sealer (#1814000)

For assays using the QX ONE™ Droplet Digital PCR System (#12006536):

- iScript Advanced cDNA Synthesis Kit (#1725037, 1725038)
- ddPCR Multiplex Supermix (#12005909, 12005910, 12005911)
- PX1 PCR Plate Sealer (#1814000)

For assays using the QX Continuum™ Droplet Digital PCR System (#12019613):

- QX Continuum ddPCR Supermix for Probes (#12019028, 12019141, 12019133)
- PX1 PCR Plate Sealer (#1814000)

For assays using QX700™ Droplet Digital PCR Systems (#17011036, #17010638, #17010628):

- naica™ 5X Multiplex ddPCR Mix (#12025253, #12025254)
- naica 5X ddPCR Mix (#12025257)
- naica 10X Multiplex ddPCR Mix (#12025255, #12025256, #12025258)
- PX1 PCR Plate Sealer (#1814000)

Refer to the relevant instrument manual for operation (Table 2).



Table 2. Instrument manuals.

| Bulletin Number | Document |
|-----------------|---|
| 10031906 | QX200 Droplet Reader and QuantaSoft™ Software Instruction Manual |
| 10031907 | QX200 Droplet Generator Instruction Manual |
| 10000116512 | QX ONE Droplet Digital PCR System and QX ONE Software Instrument Guide |
| 10000153877 | QX600 Droplet Reader and QX Manager Software Standard Edition User Guide |
| 10043138 | Automated Droplet Generator Instruction Manual (for ordering information on consumables such as oils, cartridges, gaskets, plates, and seals) |
| 10000170603 | QX Continuum Droplet Digital PCR System Instruction Manual |
| 10000171494 | QX700 System Analysis Software User Guide |
| 10000171493 | QX700 Droplet Digital PCR System Instruction Manual |

Bio-Rad, AutoDG, ddPCR, Droplet Digital, Droplet Digital PCR, QuantaSoft, QX Continuum, QX ONE, QX200, QX600, and QX700 are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2026 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 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>.

Instructions for Use

Visit the Assay Details page at [bio-rad.com/digital-assays](https://www.bio-rad.com/digital-assays) for assay-specific instructions, including recommended supermix, restriction enzyme digestion, reaction setup, thermal cycling conditions, and data acquisition and analysis. Additional guidelines can be found on the product inserts for any ddPCR supermix.



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

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 1 800 361 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