

ddPCR™ Expert Design Assays

Catalog #	Description	
12003898,12003899,12003900	ddPCR™ Expert Design Assay C1_0, C2_0, C3_0	1,000 x 20 µl reactions
12003907,12003908,12003909	ddPCR Expert Design Assay C1_5, C2_5, C3_5	1,000 x 20 µl reactions
12004115,12004116,12004117	ddPCR Expert Design Assay-C1_10, C2_10, C3_10	1,000 x 20 µl reactions
12003901,12003902,12003903	ddPCR Expert Design Assay-G1_0, G2_0, G3_0	1,000 x 20 µl reactions
12003910,12003923,12004111	ddPCR Expert Design Assay-G1_5, G2_5, G3_5	1,000 x 20 µl reactions
12004118,12004119,12004120	ddPCR Expert Design Assay-G1_10, G2_10, G3_10	1,000 x 20 µl reactions
12003904,12003905,12003906	ddPCR Expert Design Assay-Q1_0, Q2_0, Q3_0	1,000 x 20 µl reactions
12004112,12004113,12004114	ddPCR Expert Design Assay-Q1_5, Q2_5, Q3_5	1,000 x 20 µl reactions
12004121,12004122,12004123	ddPCR Expert Design Assay-Q1_10, Q2_10, Q3_10	1,000 x 20 µl reactions

For research purposes only.

Description

ddPCR Expert Design Assays are designed specifically for Droplet Digital™ PCR (ddPCR) by the experts that developed the technology. ddPCR Expert Design Assays encompass a wide variety of assays (for example, translocations, mutation detection, multiplex mutation screening, etc). Each assay has a Unique Assay ID that corresponds to a specific formulation of primers and probes. Details about each assay can be obtained online at bio-rad.com/digital-assays. 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 on bio-rad.com.

Storage and Stability

The ddPCR assays are stable for 12 months when stored at 4°C protected from light. The assay can be kept at -20°C for long-term storage.

Kit Contents

The ddPCR Expert Design Assay is a concentrated, ready-to-use primer-probe mix. Each tube of assay comes with 1,000 µl of the assay mix.

Required Reagents and Equipment

- QX100™ or QX200™ Droplet Generator (catalog #1863002 or 1864002, respectively) or Automated Droplet Generator (catalog #1864101)
- QX100 or QX200 Droplet Reader (catalog #1863003 or 1864003, respectively)
- C1000 Touch™ Thermal Cycler with 96-Deep Well Reaction Module (catalog #1851197)
- PX1™ PCR Plate Sealer (catalog #1814000)

Please refer to the QX100 or QX200 Instruction Manuals (#10026321 and 10026322 or 10031906 and 10031907, respectively) or the Automated Droplet Generator Instruction Manual (#10043138) for ordering information on consumables such as oils, cartridges, gaskets, plates, and seals.

Instructions for Use

Assay-specific instructions, including recommended supermix, restriction enzyme digestion, reaction setup, thermal cycling conditions, and data acquisition and analysis are provided on the Assay Details Page on bio-rad.com/digital-assays. Additional guidelines can be found on the product inserts for any ddPCR Supermix.



FAM and HEX are trademarks of Life Technologies Corporation.

Bio-Rad's thermal cyclers and real-time thermal cyclers are 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.

The QX100 and QX200 Droplet Digital PCR Systems 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 purposes in the field of digital PCR only. No rights are granted for diagnostic uses. 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.

Bio-Rad Laboratories, Inc.
2000 Alfred Nobel Drive, Hercules, CA 94547
1-800-424-6723