

ddPCR™ Assay Design Service Assays

Catalog #	Description	
12008226	ddPCR Assay Design Service FAM Assay	1,000 x 20 µl reactions
12008227	ddPCR Assay Design Service HEX Assay	1,000 x 20 µl reactions
12008228	ddPCR Assay Design Service Cy5 Assay	1,000 x 20 µl reactions
12008229	ddPCR Assay Design Service Cy5.5 Assay	1,000 x 20 µl reactions
12008230	ddPCR Assay Design Service 450 µM Primer	50 µl
12008231	ddPCR Assay Design Service 125 µM FAM Probe	50 µl
12008232	ddPCR Assay Design Service 125 µM HEX Probe	50 µl
12008233	ddPCR Assay Design Service 125 µM Cy5 Probe	50 µl
12008234	ddPCR Assay Design Service 125 µM Cy5.5 Probe	50 µl
12008235	ddPCR Assay Design Service 375 µM Dark Probe with Iowa Black FQ	50 µl
12008236	ddPCR Assay Design Service 375 µM Dark Probe with Iowa Black RQ-Sp	50 µl

For research purposes only.

Description

ddPCR Assay Design Service (ADS) Assays are designed specifically for Droplet Digital™ PCR (ddPCR) based on customer requests. ddPCR ADS Assays can encompass a wide variety of assays (for example, translocations, mutation detection, multiplex mutation screening, etc). An assay design may require a combination of multiple catalog numbers and unique assay IDs. Information required to formulate an assay for a particular assay design request can be obtained from the assay information pages available 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 also be ordered at bio-rad.com. Assays 12008226, 12008227, 12008228, and 12008229 each contain 9 µM primers and 5 µM probe.

Adding 4 µl of the required number of components of 12008230, 12008231, 12008232, 12008233, 12008234, 12008235, and 12008236, and water to a final volume of 100 µl results in a concentration of 18 µM primer, 5 µM probe, and 15 µM dark probe. A further 20x dilution in the ddPCR reaction results in 900 nM primer, 250 nM probe, and 750 nM probe.

Storage and Stability

The ddPCR Assays are stable through the expiration date printed on the label. Assays are stable for 12 months when stored at 4°C protected from light. The assays can be kept at -20°C for long-term storage.

Kit Contents

The ddPCR ADS Assay is a concentrated, ready-to-use primer, probe, or primer-probe mix.

Required Reagents and Equipment

- QX100™ or QX200™ Droplet Generator (catalog #1864002) or Automated Droplet Generator (catalog #1864101)
- QX100 or QX200 Droplet Reader (catalog #1864003)
- 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 Manual (#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.

For more information on assay design and optimization for ddPCR, please refer to the Droplet Digital PCR Applications Guide (bulletin 6407) and the Rare Mutation Detection Best Practices Guide (bulletin 6628).



FAM and HEX are trademarks of Life Technologies Corporation. Cy is a trademark of GE Healthcare. Iowa Black is a trademark of Integrated DNA Technologies, Inc. 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.