

QX200™ Droplet Generation Oil for EvaGreen®

Catalog #	Description
1864005	QX200 Droplet Generation Oil for EvaGreen® , 14 ml (2 x 7 ml vials), enough for 192 x 20 µl reactions
1864006	QX200 Droplet Generation Oil for EvaGreen® , 70 ml (10 x 7 ml vials), enough for 960 x 20 µl reactions

For research purposes only.

Description

QX200 Droplet Generation Oil for EvaGreen® is used to create water (sample) in oil droplets required for Droplet Digital™ PCR (ddPCR™) analysis and is formulated for use exclusively with QX200 ddPCR EvaGreen® Supermix (catalog #1864033, 1864034, 1864035, and 1864036).

Following reaction preparation using QX200 ddPCR EvaGreen® Supermix, 20 µl each of up to eight prepared samples (or blanks) and QX200 Droplet Generation Oil for EvaGreen® are transferred to a DG8 Droplet Generator Cartridge. The loaded cartridge is covered with a DG8 Gasket and placed in the QX200 Droplet Generator. There, the samples and oil are combined within the microchannels of the cartridge to create an emulsion of ~20,000 monodispersed, nanoliter-sized droplets for each of the samples.

Additional requirements for droplet generation are listed in Table 1. For complete system requirements, refer to the QX600™ or QX200 Droplet Reader and QX Manager Software Standard Edition User Guide (10000153877 or 10000107223, respectively).

Storage and Stability

QX200 Droplet Generation Oil for EvaGreen® is stable for 12 months when stored at a constant temperature of 25°C (room temperature).

Sample Preparation

Prepare the PCR reaction as described in the QX200 ddPCR EvaGreen® Supermix Product Insert (10028376).

Droplet Generation

The QX200 Droplet Generator prepares droplets for up to eight samples at a time. Droplet generation takes ~2 min for each set of eight samples (~45 min for a full 96-well plate). The detailed protocol for droplet generation can be found in the QX200 Droplet Generator Instruction Manual (10031907).

Note: QX200 Droplet Generation Oil for EvaGreen® should be used only with QX200 ddPCR EvaGreen® Supermix. The use of other supermixes can lead to poor results and potential damage to the QX600 or QX200 Droplet Readers.

Table 1. QX200 droplet generation components.

Component	Description	Catalog #
QX200 Droplet Generator	Instrument used for droplet generation	1864002
Droplet generator cartridges and gaskets	Microfluidic DG8 Cartridge used to mix sample and oil to generate droplets; DG8 Gaskets seal the cartridge to prevent evaporation and apply the pressure required for droplet formation	1864007
DG8 Cartridge Holder	Positions and holds the DG8 Cartridge in the instrument for droplet generation	1863051
QX200 ddPCR EvaGreen® Supermix	PCR supermix	1864033, 1864034, 1864035, 1864036
QX200 Buffer Control for EvaGreen®	Control	1864052
Pipets	20 µl pipet for sample loading 50 µl pipet for droplet transfer 8-channel, 200 µl pipet for oil	Rainin L-20 Rainin L50, L8-50 Rainin L8-200
Pipet tips	Filtered	Rainin GP-L10F, GP-L200F
ddPCR 96-well plates	Pkg of 25, clear well/clear shell semi-skirted plates, for use with QX600/QX200 Droplet Digital PCR Systems	12001925
Reagent reservoir	Reservoirs for dispensing reagent with multichannel pipets	2244872
Foil seal	Pierceable foil heat seal	1814040
PX1 PCR Plate Sealer	Plate sealer	1814000
8-cap strips		Any

1. Insert the DG8 Cartridge into the holder, aligning the notch in the cartridge with the upper left corner of the holder.
2. Transfer 20 µl of each prepared sample to the sample wells (middle row) of the DG8 Cartridge.
3. Dispense the QX200 Droplet Generation Oil for EvaGreen® into a reagent trough (see Table 2 for volumes required).
4. Using a multichannel pipet, fill each oil well (bottom row) with 70 µl of QX200 Droplet Generation Oil for EvaGreen® from the reagent reservoir.
5. Hook the gasket over the cartridge holder using the holes on both sides. The gasket must be securely hooked on both ends of the holder to ensure pressure is sufficient to generate droplets.
6. Open the QX200 Droplet Generator by pressing the button on the green top and place the cartridge holder into the instrument.
7. Press the button on the top again to close the door and initiate droplet generation.
8. When droplet generation is complete, all three indicator lights are solid green and the door can be opened by pressing the button. The holder, with the DG8 Cartridge still in place, can be removed.
9. Pipet 40 µl of the contents of the top wells (the droplets) into a single column of a 96-well PCR plate. Repeat for each cartridge, filling up each column across the 96-well plate.
10. To avoid evaporation, seal the PCR plate with foil immediately after transferring all the columns of droplets.
11. Begin thermal cycling (PCR) within 30 min of sealing the plate, or store the plate at 4°C for up to 4 hr prior to thermal cycling. Refer to the QX200 ddPCR EvaGreen® Supermix Product Insert for cycling conditions.

BIO-RAD, DDPCR, DROPLET DIGITAL, DROPLET DIGITAL PCR, QX600, and QX200 are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen Dye for use in real-time PCR, for research purposes only.

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>.

Table 2. QX200 Droplet Generation Oil requirements.

Number of Wells	Volume of Oil, µl
8	700
24	1,820
48	3,500
96	6,860

When PCR amplification is complete, remove the 96-well plate from the thermal cycler and read the droplets using the QX600 or the QX200 Droplet Readers (follow the directions in the corresponding system user guide).

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