

QX200™ ddPCR™ EvaGreen® Supermix



Bio-Rad's QX200 ddPCR EvaGreen supermix is a ready-to-use 2x supermix that contains a dsDNA binding dye that enables double-stranded DNA detection following PCR amplification. It is the only dye chemistry optimized for digital PCR.

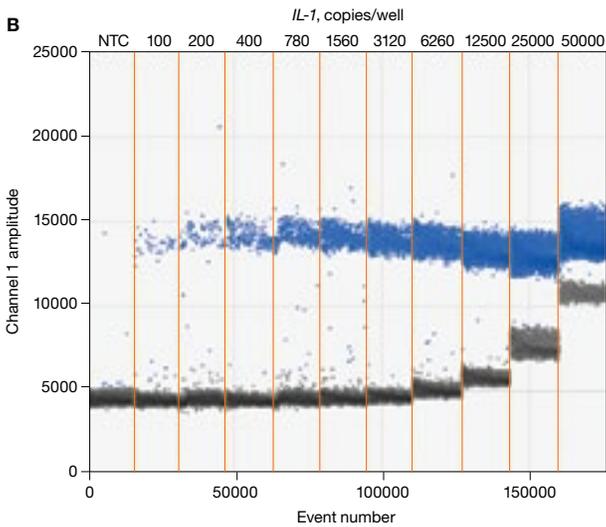
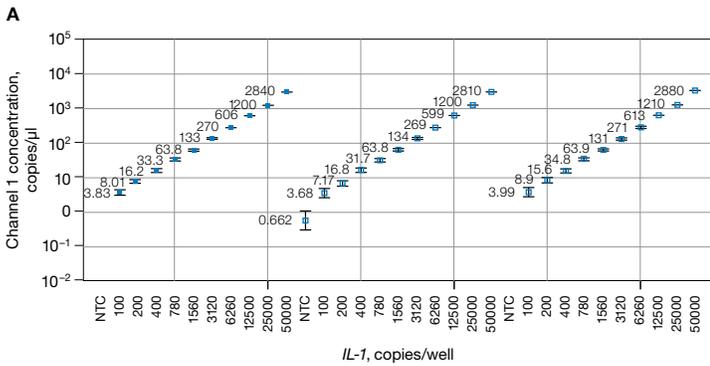
When used with Bio-Rad's QX200 droplet generator, the QX200 ddPCR EvaGreen supermix allows you to:

- Obtain precise and sensitive absolute concentrations for a variety of applications, such as gene expression, copy number detection, and DNA rearrangements
- Simplify your experiment preparation and decrease your running costs by eliminating probes
- Discriminate between spliced variants by size with the use of simple primer pairs

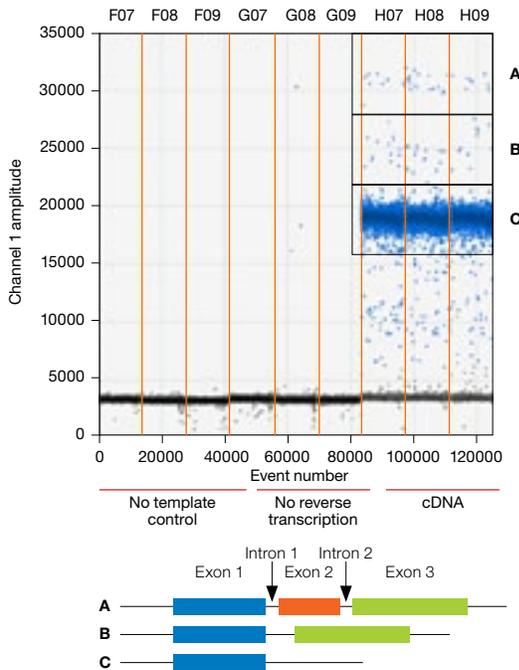
For more information, visit bio-rad.com/web/QX200evagreensmx.

BIO-RAD

Performance Data for QX200 ddPCR EvaGreen Supermix



ddPCR enables precise and reproducible detection calls across a range of sample concentrations. **A**, concentration plot shows data for both merged wells (solid blue markers) and individual replicate wells (open blue markers) across a dilution series. Interleukin 1 (*IL-1*) samples ranging from 0 to 50,000 copies were detected with an EvaGreen assay and plotted on a log scale. There is excellent linearity and reproducibility between replicates. **B**, 1-D fluorescence amplitude plot shows a sample dilution curve detected with an EvaGreen *IL-1* assay in the FAM/EvaGreen channel of the QX200 droplet reader. The data demonstrate good separation between positive (■) and negative (■) data points. All error bars generated by QuantaSoft™ software represent the 95% confidence interval. NTC, no template control.



Splice variant discrimination using the QX200 ddPCR EvaGreen supermix with a single set of primer pairs. **A**, primary transcript containing exons 1, 2, and 3; **B**, transcript missing exon 2; **C**, transcript missing exons 2 and 3. Positive droplets containing longer amplicons (population A) have a higher fluorescence amplitude and can be distinguished from positive droplets containing smaller amplicons (populations B and C).

Ordering Information

Catalog #	Description
186-4033	QX200 ddPCR EvaGreen Supermix , 2 ml (2 x 1 ml), 200 x 20 μl reactions, 2x supermix
186-4034	QX200 ddPCR EvaGreen Supermix , 5 ml (5 x 1 ml), 500 x 20 μl reactions, 2x supermix
186-4035	QX200 ddPCR EvaGreen Supermix , 25 ml (5 x 5 ml), 2,500 x 20 μl reactions, 2x supermix
186-4036	QX200 ddPCR EvaGreen Supermix , 50 ml (10 x 5 ml), 5,000 x 20 μl reactions, 2x supermix

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. FAM is a trademark of Applied Biosystems Corporation.

The QX200 Droplet Digital PCR system and/or its 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 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.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 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** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000