



## Droplet Digital PCR (ddPCR) Multiplex Supermix

Bio-Rad's ddPCR Multiplex Supermix is a ready-to-use 4x reaction mix that contains all of the components required for probe-based Droplet Digital PCR except for primers, probe(s), and template. The ddPCR Multiplex Supermix has been optimized to deliver maximum PCR efficiency, specificity, and sensitivity when used with the QX200 Droplet Digital PCR System.

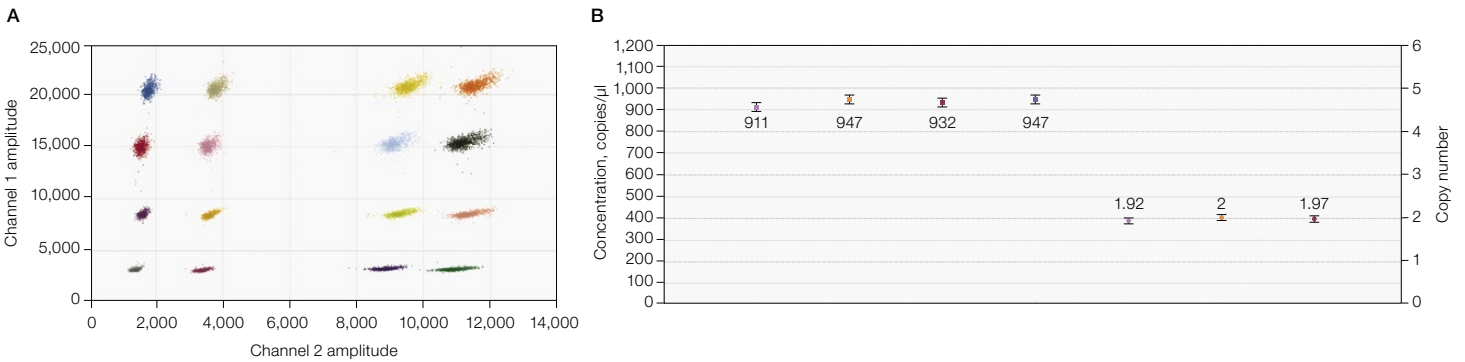
Combined with Bio-Rad's QX200 ddPCR System, the ddPCR Multiplex Supermix allows you to:

- Amplify and detect multiple targets using commercially available probe-based assays
- Input larger amount of samples
- Obtain precise and sensitive absolute measurements for a variety of applications, such as copy number analysis and mutation detection

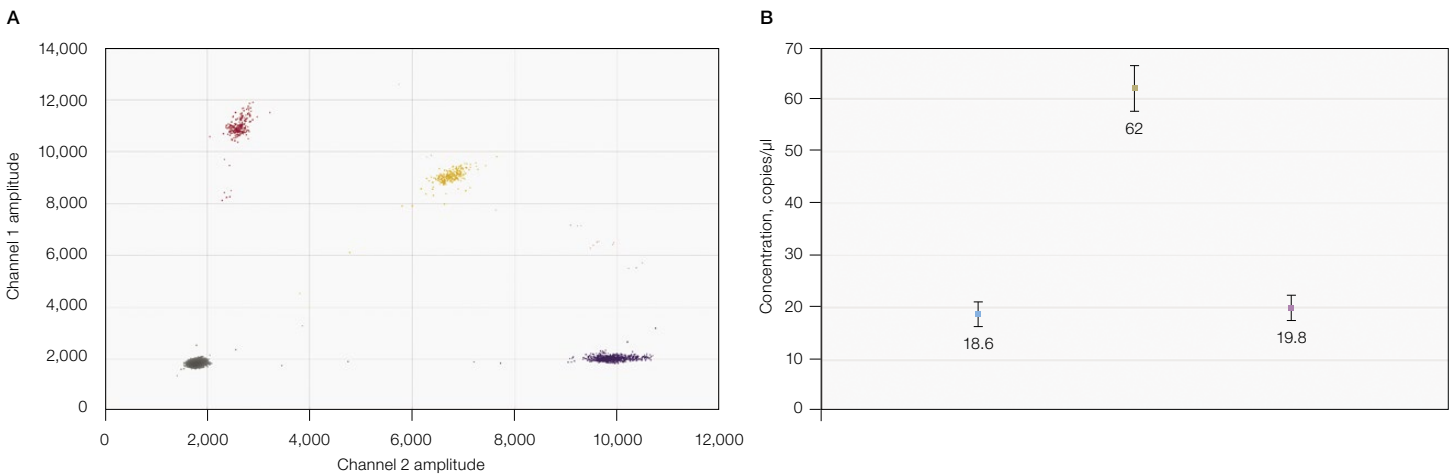
Visit [bio-rad.com/ddPCRMultiplexsmx](https://www.bio-rad.com/ddPCRMultiplexsmx) for more information.



## ddPCR Multiplex Supermix Sample Data



**Determination of copy number of four different targets in a single well.** **A**, 2-D fluorescence amplitude plot shows amplitude multiplexing of *ERBB2*, *PIK3CA*, *BRAF V600*, and *RPP30* using Channel 1 (FAM) and Channel 2 (HEX) of the QX200 ddPCR System; **B**, concentration plot shows copies/ $\mu$ l of *ERBB2* (■), *PIK3CA* (■), *BRAF V600* (■), and *RPP30* (■) in a single well. Copy number shows data points of *ERBB2* (■), *PIK3CA* (■), *BRAF V600* (■). Error bars represent a 95% confidence interval.



**Radial multiplexing of three targets (two mutants and one wild type).** **A**, 2-D fluorescence amplitude plot shows radial multiplexing of *KRAS G12C*, *KRAS G12D*, and wild type using Channel 1 (FAM) and Channel 2 (HEX) of the QX200 ddPCR System. *KRAS G12C* is detected by a FAM probe, *KRAS G12D* by a 50% FAM and 50% HEX probe, and *KRAS G12* wild type by a HEX probe. A cluster in gray is double-negative for the targets, red is *KRAS G12C*, orange is *KRAS G12D*, and purple is *KRAS G12* wild type; **B**, concentration plot shows copies/ $\mu$ l of *KRAS G12C* (■), *KRAS G12D* (■), and *KRAS G12* wild type (■). Error bars represent a 95% confidence interval.

## Ordering Information

Catalog #	Description
12005909	<b>ddPCR Multiplex Supermix</b> , 1 ml (2 x 0.5 ml), 200 x 20 $\mu$ l reactions, 4x supermix
12005910	<b>ddPCR Multiplex Supermix</b> , 2.5 ml (5 x 0.5 ml), 500 x 20 $\mu$ l reactions, 4x supermix
12005911	<b>ddPCR Multiplex Supermix</b> , 12.5 ml (5 x 2.5 ml), 2,500 x 20 $\mu$ l reactions, 4x supermix

Bio-Rad, Droplet Digital PCR, and ddPCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

The QX200 Droplet Digital PCR System, and the consumables and reagents designed to work with this 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. See [bio-rad.com/en-us/trademarks](http://bio-rad.com/en-us/trademarks) for details. 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.**

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600  
**Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30  
**Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580  
**Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300  
**United Kingdom** 44 01923 47 1301

