

## ddPCR™ *KRAS* Screening Multiplex Kit



The ddPCR *KRAS* Screening Multiplex Kit is designed to screen for seven *KRAS* mutations. The kit contains a 20x ddPCR *KRAS* Screening Multiplex Assay and ddPCR Supermix for Probes (No dUTP).

Combined with Bio-Rad's QX100™, QX200™, or QX200™ AutoDG™ Droplet Digital PCR (ddPCR) System, the ddPCR *KRAS* Screening Multiplex Kit allows you to:

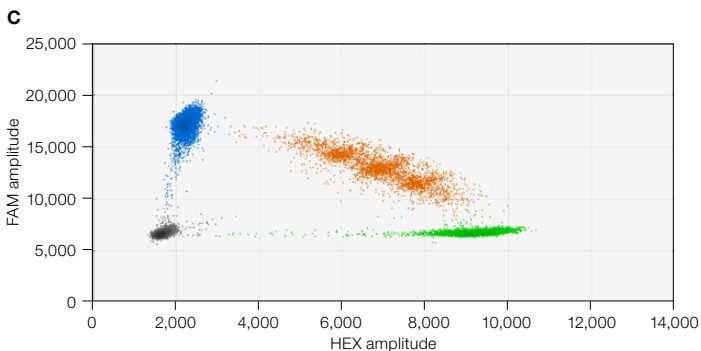
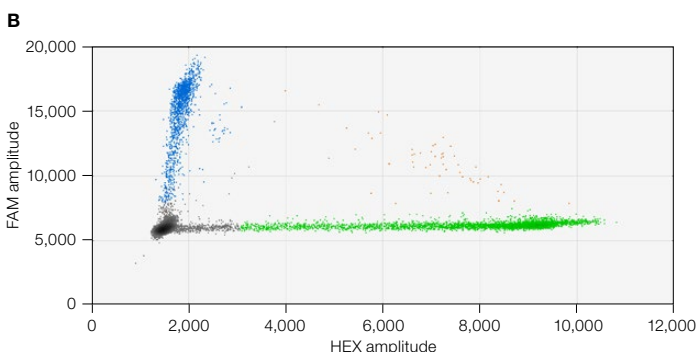
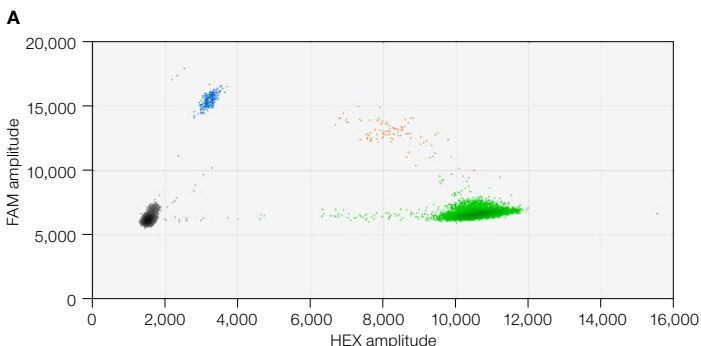
- Quantify and screen for multiple *KRAS* mutations in a single well
- Obtain sensitive and precise detection down to 0.2% mutant levels
- Screen multiple samples in a rapid and cost-effective manner

Visit [bio-rad.com/web/KRASscrmxkit](http://bio-rad.com/web/KRASscrmxkit) for more information.

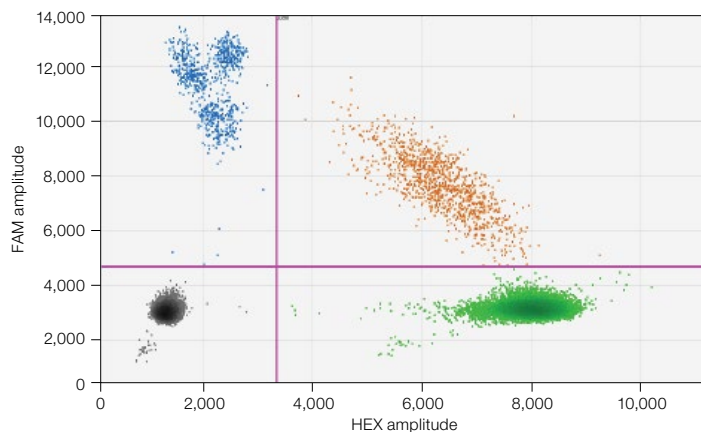


## Quantification and Screening of KRAS Mutations

The ddPCR KRAS Screening Multiplex Kit contains seven KRAS mutant assays (G12A, G12C, G12D, G12V, G12R, G12S, G13D) and one wild-type assay. The kit allows for rapid identification of multiple KRAS mutations down to 0.2% mutant in a single well. No preamplification step is required for use of this kit.



**ddPCR KRAS Screening Multiplex Kit used to screen various sample types.** 2-D fluorescence amplitude plots show the wells of cell-line (A), formalin-fixed, paraffin-embedded (B), and cell-free DNA samples (C). The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for wild-type DNA only, the blue cluster represents the droplets that are positive for mutant DNA only, and the orange cluster represents the droplets that are positive for both mutant and wild-type DNA.



**Multiplexed single-well detection of seven KRAS mutations.** A 2-D fluorescence amplitude plot generated by QuantaSoft™ Software shows the wells of a mixed mutant and wild-type sample. The black cluster on the plot represents the negative droplets, the green cluster represents the droplets that are positive for wild-type DNA only, the blue cluster represents the droplets that are positive for mutant DNA only, and the orange cluster represents the droplets that are positive for both mutant and wild-type DNA.

## Ordering Information

Catalog #	Description
186-3506	ddPCR KRAS Screening Multiplex Kit, 200 x 20 µl reactions, includes 20x ddPCR KRAS Screening Multiplex Assay (1 x 200 µl vial), ddPCR Supermix for Probes (No dUTP) (2 x 1 ml vials)

FAM is a trademark of Applied Biosystems.

The QX100 or 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**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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 3065 7550  
**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** 81 3 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** 1800 88 22 88 **United Kingdom** 020 8328 2000