



ddPCR™ Multiplex Mutation Screening Kits

Droplet Digital™ PCR (ddPCR) Multiplex Mutation Screening Kits are designed for rapid screening of several key cancer mutations in a single reaction with high precision and sensitivity. These kits are ideal for use with formalin-fixed, paraffin-embedded (FFPE) samples, liquid biopsy, fresh/frozen tissue, and samples with low yield or inhibitory substances. The kits contain an optimized, validated ddPCR Mutation Screening Multiplex Assay and ddPCR Supermix for Probes (No dUTP).

- Screen multiple mutations in a single well
- Detect with high sensitivity (limit of detection $\leq 0.5\%$)
- Process numerous samples in a rapid, cost-effective manner
- Work with low amounts of input DNA without preamplification
- Use with Bio-Rad's QX100™, QX200™, or QX200™ AutoDG™ Droplet Digital PCR System

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



Mutations Detected by ddPCR Multiplex Mutation Screening Kits

Catalog Number	ddPCR Multiplex Mutation Screening Kit*	Mutations Detected
12001037	<i>BRAF</i> V600	V600E, V600K, V600R
1863506	<i>KRAS</i> G12/G13	G12A, G12C, G12D, G12R, G12S, G12V, G13D
12001626	<i>KRAS</i> Q61	Q61H (183A>C), Q61H (183A>T), Q61K, Q61L, Q61R
12001094	<i>NRAS</i> G12	G12A, G12C, G12D, G12R, G12S, G12V
12001627	<i>NRAS</i> G12/G13	G12A, G12C, G12D, G12S, G12V, G13D, G13R, G13V
12001006	<i>NRAS</i> Q61	Q61H (183A>C), Q61H (183A>T), Q61K, Q61L, Q61R
12002392	<i>EGFR</i> exon 19 deletions	>15 <i>EGFR</i> exon 19 deletions

* Assays for each of the mutations detected by these kits can also be ordered individually. Each kit contains 20x ddPCR Mutation Screening Multiplex Assay (1 x 200 µl) and 2x ddPCR Supermix for Probes (No dUTP) (2 x 1 ml).

Key Actionable Mutations in Various Cancers

Gene	Lung Cancer	Colorectal Cancer	Ovarian Cancer	Thyroid Cancer	Melanoma	GIST	Glioma	Pancreatic Cancer
<i>BRAF</i>	•	•	•	•	•	•	•	
<i>KRAS</i>	•	•	•	•				•
<i>NRAS</i>	•	•		•	•			
<i>EGFR</i>	•	•			•			

GIST, gastrointestinal stromal tumor.

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 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.

The Droplet Digital PCR (ddPCR) Multiplex Mutation Screening Kits are intended for research use only.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050
Italy 39 02 216091 **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 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 (0) 861 246 723
Spain 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311
United Arab Emirates 971 4 8187300 **United Kingdom** 44 020 8328 2000

