



ddPLEX *EGFR/KRAS/BRAF* Mutation Detection Kit

Detect NSCLC Mutations with High Sensitivity and Precision

The ddPLEX *EGFR/KRAS/BRAF* Mutation Detection Kit delivers highly sensitive and accurate results, so you can confidently detect 37 key *EGFR*, *KRAS*, and *BRAF* mutations in non-small cell lung cancer (NSCLC) down to 0.1% variant allele frequency or lower, as well as key *EGFR* resistance mutations such as *T790M* and *C797S*. The kit supports both tissue-based workflows and less-invasive liquid biopsy methods, enabling analysis of circulating tumor DNA (ctDNA) from plasma.

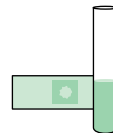
The kit offers:

High Analytical Sensitivity



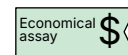
Sensitivity
as low as 0.025%

Plasma or FFPE Tissue



Liquid biopsy
and tissue-based
analysis options

Cost-Effective Mutation Detection



Multimarker analysis
at a fraction of the
cost of NGS workflows

Same Day Results



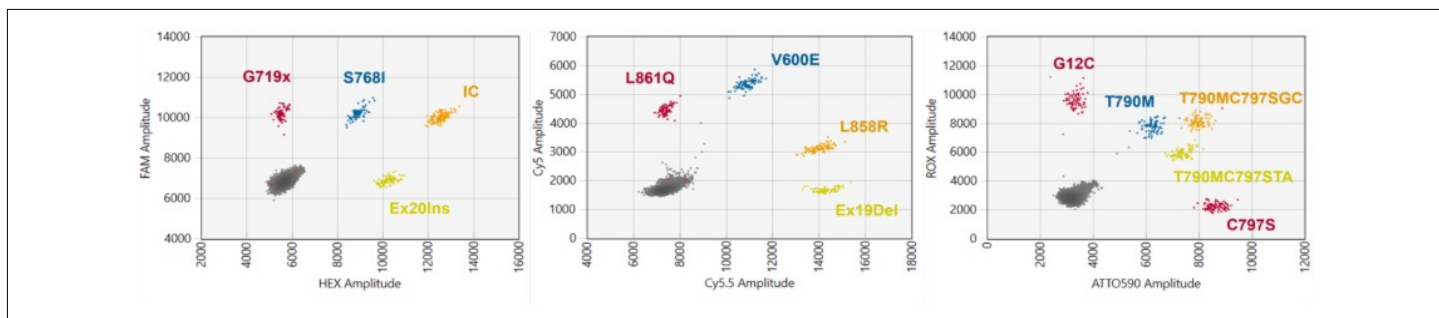
Sample input
to answer in a
single work shift



Scan the QR code or visit
bio-rad.com/ddPLEX-NSCLC
to learn more.

Multiplexed Detection: Visualizing NSCLC Variants Across Three Channels

The kit provides the ability to detect key NSCLC-relevant variants. In this example, separated clusters are shown across six channels (FAM, HEX, Cy5, Cy5.5, ROX, and ATTO590) on the QX600™ Droplet Digital™ PCR System.



Targeted NSCLC-Relevant Variants

Gene	Exon 18	Exon 19	Exon 20	Exon 21	Exon 2	Exon 15
EGFR	G719A, G719S, G719C	Deletions	Insertions, S768I, T790M, C797S (cis/trans)	L861Q		
KRAS					G12C	
BRAF						V600E

Detect Variants with High Sensitivity and Accuracy

The kit exhibits a low limit of detection (LoD) of 0.025% to 0.1% variant allele frequency (VAF). LoD was tested with mutant samples at 0.2, 0.1, 0.05, 0.025, and 0.01% in the background of wild-type DNA. LoD was determined as the lowest variant allele frequency at which all replicates were detected.

Gene	Target	LoD
EGFR	G719x (n=3)	0.025–0.05%
	S768I	0.025%
	Ex20Ins (n=7)	0.025–0.05%
	L861Q	0.025%
	Ex19Del (n=7)	0.05%
	G12C	0.025%
	T790M	0.10%
	T790MC797SGC	0.025%
	T790MC797STA	0.025%
KRAS	C797S	0.025%
	G12C	0.025%
BRAF	V600E	0.025%

Ordering Information

ddPLEX EGFR/KRAS/BRAF Mutation Detection Kit Contents

Catalog #	Component (also available separately)
17010491	ddPLEX EGFR/KRAS/BRAF Mutation Detection Kit
12023484	ddPLEX EGFR/KRAS/BRAF Mutation Detection Assay <ul style="list-style-type: none"> • ddPLEX Mutant Assay, 60 reactions • ddPLEX Total Quant Assay, 60 reactions • ddPLEX Enhancer, 1 vial x 90 µL • ddPLEX Internal Control, 1 vial x 500 µL
12023522	ddPLEX EGFR/KRAS/BRAF Positive Control, 1 vial x 100 µL
12005909*	ddPCR Multiplex Supermix, 2 vials x 600 µL each
12012171	Dithiothreitol (DTT), 2 vials x 1000 µL each

*Other ddPCR Multiplex Supermix packaging options are also available (catalog #12005910 and 12005911).

Visit bio-rad.com/ddPLEX-NSCLC for more information.

For Research Use Only. Not for use in diagnostic procedures.

Bio-Rad, QX600, and Droplet Digital PCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

© 2026 Bio-Rad Laboratories, Inc.

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 800 361 1808 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23