



Droplet Digital PCR (ddPCR) Expert Design Assays

ddPCR Expert Design Assays have been designed by Bio-Rad Droplet Digital PCR experts for collaborations or high-value targets outside of the current core application assay menu (mutation detection, copy number variation, gene edit detection, and gene expression).

ddPCR Expert Design Assays can be ordered at bio-rad.com/digital-assays using their Unique Assay ID. They are not wet-lab validated; it is recommended that the assays be validated prior to use by the end user. Information for each assay is available to download online.



Available Assays

For research use only.

Unique Assay ID	Assay Name
dEXD12170542	AAV-ITR2 (FAM)
dEXD76440315	AAV-ITR2 (HEX)
dEXD24979861	eGFP (HEX)
dHsaEXD29194342	BCR-ABL p190/ABL1 mRNA Fusion
dHsaEXD26637642	BCR-ABL p190/GusB mRNA Fusion
dHsaEXD16080615	BCR-ABL p210 mRNA Fusion
dHsaEXD39032815	ABL1 T315I Mutation
dHsaEXD64763042	BRAF V600E TG>AA Mutation
dHsaEXD91642442	EGFR T790M/C797S Cis/Trans
dHsaEXD42003661	EGFR C797S Multiplex Tube 1
dHsaEXD93464115	EGFR T790M/L858R Multiplex Tube 2
dHsaEXD40054642	EGFR G719X/S768I/L861Q Multiplex
dHsaEXD41810842	EGFR T790M/L858R/C797S Multiplex (QX ONE)
dHsaEXD61571942	IDH1 132 Multiplex
dHsaEXD35841715	IDH2 140 Multiplex
dHsaEXD10111488	IDH2 172 Multiplex
dHsaEXD86850342	EML4-ALK mRNA Fusions
dHsaEXD81378442	RET mRNA Fusions
dHsaEXD73338942	ROS1 mRNA Fusions
dHsaEXD64200342	KRAS G12C/D/V Semi Discriminatory Multiplex
dHsaEXD59551642	KRAS G12C/D/V Semi Discriminatory Multiplex (GMP + QC)
dHsaMDXE91450042	ESR1 Multiplex 1
dHsaMDXE65719815	ESR1 Multiplex 2
dEXD63637642	Library Quantification Assay for Ion Torrent
dHsaEXD20945488	TERT C228T_88
dHsaEXD72405942	TERT C228T_113 (recommended assay)
dHsaEXD59485034	TERT C228_163
dHsaEXD85215261	TERT C250T_88
dHsaEXD46675715	TERT C250T_113 (recommended assay)

Unique Assay ID	Assay Name
dHsaEXD33754807	TERT C250T_163
dHsaEXD29156242	HLA-DRB1*01 (FAM)
dHsaEXD61594642	HLA-DRB1*01 (HEX)
dHsaEXD93426015	HLA-DRB1*03 (FAM)
dHsaEXD35864415	HLA-DRB1*03 (HEX)
dHsaEXD67695788	HLA-DRB1*04 (FAM)
dHsaEXD10134188	HLA-DRB1*04 (HEX)
dHsaEXD41965561	HLA-DRB1*07 (FAM)
dHsaEXD74403961	HLA-DRB1*07 (HEX)
dHsaEXD16235334	HLA-DRB1*08 (FAM)
dHsaEXD48673734	HLA-DRB1*08 (HEX)
dHsaEXD80505107	HLA-DRB1*11 (FAM)
dHsaEXD22943507	HLA-DRB1*11 (HEX)
dHsaEXD54774880	HLA-DRB1*13 (FAM)
dHsaEXD87213280	HLA-DRB1*13 (HEX)
dHsaEXD29044653	HLA-DRB1*15 (FAM)
dHsaEXD61483053	HLA-DRB1*15 (HEX)
dHsaEXD93314426	HLA-DRB1 Region 1 (Pre-Amp)
dHsaEXD67584199	HLA-DRB1 Region 2 (Pre-Amp)
dEXD90907415	ZIKA NS5
dEXD65177188	EBOLA VP24
dHsaEXD54774880	HLA-DRB1*13
dHsaEXD32377342	Pre-AMP EGFR ex19
dHsaEXD45186661	Pre-AMP BRAF V600
dHsaEXD70916888	Pre-AMP EGFR L858
dHsaEXD82171642	Pre-AMP EBK Mix 2 Tube 2
dHsaEXD83726207	Pre-AMP EBK Mix 2 Tube 1
dHsaEXD96647115	Pre-AMP EGFR T79X
dEXD88164642	CD-19 CAR-T Assay

Unique Assay ID	Assay Name	Unique Assay ID	Assay Name
dEXD15274642	AAV ITR-2 (FAM)	dEXD76433053	SARS-CoV-2 Influenza Multiplex
dEXD23004642	AAV ITR-2 (HEX)	dEXD14924415	<i>Enterococcus</i> spp.
dEXD19233256	CMV enhancer (HEX)	dEXD27733734	Norovirus
dEXD44963483	CMV enhancer (FAM)	dEXD53463961	<i>Legionella</i>
dEXD19344845	SV40 poly A (FAM)	dEXD79194188	HF183
dEXD83614618	SV40 polyA (HEX)	dEXD92003507	Adenovirus
dEXD55784415	SV40 Poly A (Cy5)	dEXD77482599	RSV A & RSV B
dEXD32154164	BGH poly A (HEX)	dEXD13324415	WPRE (HEX)
dEXD57884391	BGH poly A (FAM)	dEXD14812826	HivPsi (FAM)
dEXD70693710	CMV promoter (HEX)	dEXD26133734	NeoR/KanR (FAM)
dEXD96423937	CMV promoter (FAM)	dEXD38943053	PuroR (HEX)
dEXD30054188	CMV promoter (CY5.5)	dEXD39054642	WPRE (FAM)
dEXD72134642	hGFP (FAM)	dEXD40543053	HivLtr (HEX)
dEXD96535526	hGFP (HEX)	dEXD51863961	KanR (HEX)
dEXD97544642	HGH poly A (FAM)	dEXD64673280	PuroR (FAM)
dEXD71814415	HGH poly A (HEX)	dEXD66273280	HivLtr (FAM)
dEXD22434642	eGFP (HEX)	dEXD77594188	KanR (FAM)
dEXD45075072	eGFP (FAM)	dEXD79082599	HivPsi (HEX)
dEXD84844642	Influenza A/B Triplex Assay	dEXD90403507	NeoR/KanR (HEX)
dEXD15704188	SARS-CoV-2 N1/N2 Assay	dEXD15841942	Variant SARS-CoV-2 Multiplex
dEXD28513507	SARS-CoV-2 N2 (FAM/HEX)	dHsaEXD25084188	ddPCR ABL1_E255K cDNA
dEXD41434415	SARS-CoV-2 N1/N2 BovineRSV Assay	dHsaEXD89353961	ddPCR ABL1_F317L cDNA
dEXD54243734	SARS-CoV-2 N1 (FAM)	dHsaEXD63623734	ddPCR ABL1_Y253H cDNA
dEXD79973961	Bovine RSV (HEX)	dHsaEXD37893507	ddPCR ABL1_V299L cDNA
dEXD40654642	SARS-CoV-2 Influenza (RT Primers)	dHsaEXD12163280	ddPCR ABL1_F359V cDNA

BIO-RAD, DDPCR, DROPLET DIGITAL PCR, and QX ONE are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

Purchase of Digital PCR products from Bio-Rad Laboratories includes a limited, non-transferable right under our intellectual property for you to use the products for your internal research purposes only. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the products or to use the products for commercial applications, such as contract services or fee for services. No rights are granted for diagnostic uses. 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 QX ONE 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. See bio-rad.com/en-us/trademarks for details.



**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 1 905 364 3435 **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 2 3473 4460 **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

