

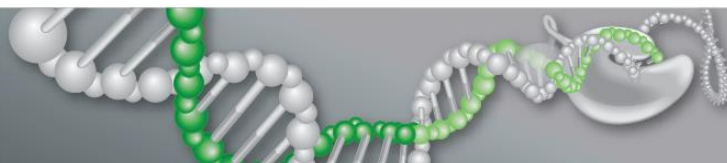
Droplet Digital™ PCR (ddPCR™) Assay Pricing

Product	Catalog #	Description	List Price (JPY)
ddPCR Mutation Detection Assays (validated)	10031246	ddPCR MUT FAM assay, 200 reactions, validated	¥ 38,000
	10031247	ddPCR MUT FAM assay, 1,000 reactions, validated	¥ 114,000
	10031248	ddPCR MUT FAM assay, 2,500 reactions, validated	¥ 190,000
	10031249	ddPCR MUT HEX assay, 200 reactions, validated	¥ 42,000
	10031250	ddPCR MUT HEX assay, 1,000 reactions, validated	¥ 126,000
	10031251	ddPCR MUT HEX assay, 2,500 reactions, validated	¥ 210,000
	10049550	ddPCR MUT FAM and HEX assay, 200 reactions, validated	¥ 72,000
	10049551	ddPCR MUT FAM and HEX assay, 1,000 reactions, validated	¥ 216,000
	10049552	ddPCR MUT FAM and HEX assay, 2,500 reactions, validated	¥ 360,000
	12005576	ddPCR MUT Cy5 and Cy5.5 assay, 200 reactions, validated	¥ 94,000
	12005577	ddPCR MUT Cy5 and Cy5.5 assay 1,000 reactions, validated	¥ 238,000
	12005578	ddPCR MUT Cy5 and Cy5.5 assay 2,500 reactions, validated	¥ 390,000
ddPCR Mutation Detection Assays	10042964	ddPCR MUT FAM assay, 200 reactions	¥ 36,000
	10042965	ddPCR MUT FAM assay, 1,000 reactions	¥ 106,000
	10042966	ddPCR MUT FAM assay, 2,500 reactions	¥ 176,000
	10042967	ddPCR MUT HEX assay, 200 reactions	¥ 40,000
	10042968	ddPCR MUT HEX assay, 1,000 reactions	¥ 118,000
	10042969	ddPCR MUT HEX assay, 2,500 reactions	¥ 196,000
	12008220	ddPCR MUT Cy5 assay, 200 reactions	¥ 42,000
	12008221	ddPCR MUT Cy5 assay, 1,000 reactions	¥ 160,000
	12008222	ddPCR MUT Cy5 assay, 2,500 reactions	¥ 220,000
	12008223	ddPCR MUT Cy5.5 assay, 200 reactions	¥ 42,000
	12008224	ddPCR MUT Cy5.5 assay, 1,000 reactions	¥ 160,000
	12008225	ddPCR MUT Cy5.5 assay, 2,500 reactions	¥ 220,000
	10049047	ddPCR MUT FAM and HEX assay, 200 reactions	¥ 66,000
	10049048	ddPCR MUT FAM and HEX assay, 1,000 reactions	¥ 198,000
	10049049	ddPCR MUT FAM and HEX assay, 2,500 reactions	¥ 330,000
	12005579	ddPCR MUT Cy5 and Cy5.5 assay, 200 reactions	¥ 86,000
	12005580	ddPCR MUT Cy5 and Cy5.5 assay, 1,000 reactions	¥ 220,000
	12005581	ddPCR MUT Cy5 and Cy5.5 assay, 2,500 reactions	¥ 360,000
ddPCR Copy Number Variation Assays (validated)	10031240	ddPCR CNV FAM assay, 200 reactions, validated	¥ 34,000
	10031241	ddPCR CNV FAM assay, 1,000 reactions, validated	¥ 84,000
	10031242	ddPCR CNV FAM assay, 2,500 reactions, validated	¥ 132,000
	10031243	ddPCR CNV HEX assay, 200 reactions, validated	¥ 38,000
	10031244	ddPCR CNV HEX assay, 1,000 reactions, validated	¥ 94,000
	10031245	ddPCR CNV HEX assay, 2,500 reactions, validated	¥ 148,000
ddPCR Copy Number Variation Assays	10042958	ddPCR CNV FAM assay, 200 reactions	¥ 32,000
	10042959	ddPCR CNV FAM assay, 1,000 reactions	¥ 78,000
	10042960	ddPCR CNV FAM assay, 2,500 reactions	¥ 124,000
	10042961	ddPCR CNV HEX assay, 200 reactions	¥ 36,000
	10042962	ddPCR CNV HEX assay, 1,000 reactions	¥ 88,000
	10042963	ddPCR CNV HEX assay, 2,500 reactions	¥ 140,000
	12005520	ddPCR CNV Cy5 assay, 200 reactions	¥ 44,000
	12005571	ddPCR CNV Cy5 assay 1,000 reactions	¥ 122,000
	12005572	ddPCR CNV Cy5 assay, 2,500 reactions	¥ 186,000
	12005573	ddPCR CNV Cy5.5 assay, 200 reactions	¥ 44,000
	12005574	ddPCR CNV Cy5.5 assay, 1,000 reactions	¥ 122,000
	12005575	ddPCR CNV Cy5.5 assay, 2,500 reactions	¥ 186,000

continued

注1: 本製品のプライマーおよびプローブ配列は非開示です。コントロール配列等は製品検索ページにありますMIQE Contextをご参照ください。

注2: 本製品は研究用試薬であり、QX100/QX200 ddPCR PCRシステムおよびQX ONE ddPCRシステムにおける専用試薬です。診断および治療目的には使用できません。



Droplet Digital™ PCR (ddPCR™) Assay Pricing (continued)

Product	Catalog #	Description	List Price (JPY)
ddPCR Gene Expression Assays	10031252	ddPCR GEX FAM assay, 200 reactions	¥ 32,000
	10031253	ddPCR GEX FAM assay, 1,000 reactions	¥ 78,000
	10031254	ddPCR GEX FAM assay, 2,500 reactions	¥ 124,000
	10031255	ddPCR GEX HEX assay, 200 reactions	¥ 36,000
	10031256	ddPCR GEX HEX assay, 1,000 reactions	¥ 88,000
	10031257	ddPCR GEX HEX assay, 2,500 reactions	¥ 140,000
	12005582	ddPCR GEX Cy5 assay, 200 reactions	¥ 44,000
	12005583	ddPCR GEX Cy5 assay, 1,000 reactions	¥ 78,000
	12005584	ddPCR GEX Cy5 assay, 2,500 reactions	¥ 152,000
	12005585	ddPCR GEX Cy5.5 assay, 200 reactions	¥ 44,000
	12005586	ddPCR GEX Cy5.5 assay, 1,000 reactions	¥ 78,000
	12005587	ddPCR GEX Cy5.5 assay, 2,500 reactions	¥ 152,000
	10031258	ddPCR™ GEX EvaGreen® assay, 200 reactions	¥ 14,000
	10031259	ddPCR™ GEX EvaGreen® assay, 1,000 reactions	¥ 36,000
	10031260	ddPCR™ GEX EvaGreen® assay, 2,500 reactions	¥ 56,000
Custom ddPCR Assays	10031276	Custom ddPCR FAM assay, 200 reactions	¥ 32,000
	10031277	Custom ddPCR FAM assay, 1,000 reactions	¥ 74,000
	10031278	Custom ddPCR FAM assay, 2,500 reactions	¥ 116,000
	10031279	Custom ddPCR HEX assay, 200 reactions	¥ 36,000
	10031280	Custom ddPCR HEX assay, 1,000 reactions	¥ 84,000
	10031281	Custom ddPCR HEX assay, 2,500 reactions	¥ 132,000
	12005588	Custom ddPCR Cy5 assay, 200 reactions	¥ 44,000
	12005589	Custom ddPCR Cy5 assay, 1000 reactions	¥ 62,000
	12005590	Custom ddPCR Cy5 assay, 2,500 reactions	¥ 86,000
	12005591	Custom ddPCR Cy5.5 assay, 200 reactions	¥ 44,000
	12005592	Custom ddPCR Cy5.5 assay, 1,000 reactions	¥ 62,000
	12005593	Custom ddPCR Cy5.5 assay, 2,500 reactions	¥ 86,000
	10031282	Custom ddPCR™ EvaGreen® assay, 200 reactions	¥ 20,000
	10031283	Custom ddPCR™ EvaGreen® assay, 1,000 reactions	¥ 30,000
	10031284	Custom ddPCR™ EvaGreen® assay, 2,500 reactions	¥ 48,000

注1: 本製品のプライマーおよびプローブ配列は非開示です。コントロール配列等は製品検索ページにありますMIQE Contextをご参照ください。

注2: 本製品は研究用試薬であり、QX100/QX200 ddPCR PCRシステムおよびQX ONE ddPCRシステムにおける専用試薬です。診断および治療目的には使用できません。

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen dye for use in real-time PCR, for research purposes only. FAM and HEX are trademarks of Life Technologies Corporation.

バイオ・ラッド ラボラトリーズ 株式会社
ライフサイエンス

Visit us at <http://www.bio-rad.com>

本 社	〒140-0002	東京都品川区東品川 2-2-24 天王洲ビル7F	TEL:03-6361-7000	FAX:03-5463-8480
大阪営業所	〒532-0025	大阪市淀川区新北野1-14-11 第一生命ビル	TEL:06-6308-6568	FAX:06-6308-3064
		*学術的お問い合わせは	TEL:03-6404-0331	FAX:03-6404-0334

取扱店