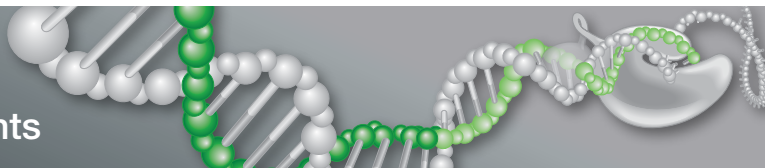


PrimePCR™ Pricing and Bulk Discounts



Assay Bulk Discounts

25% off 5 or more assays

Droplet Digital™ PCR (ddPCR™) Probe Assay Pricing

Product	Catalog #	Description	List Price (USD)
Predesigned ddPCR Copy Number Variation Probe Assays	10031240	ddPCR CNV FAM assay, 200 reactions	\$ 210
	10031241	ddPCR CNV FAM assay, 1,000 reactions	\$ 825
	10031242	ddPCR CNV FAM assay, 2,500 reactions	\$ 1,125
	10031243	ddPCR CNV HEX assay, 200 reactions	\$ 210
	10031244	ddPCR CNV HEX assay, 1,000 reactions	\$ 825
	10031245	ddPCR CNV HEX assay, 2,500 reactions	\$ 1,125
Predesigned ddPCR Mutation Detection Probe Assays	10031246	ddPCR MUT FAM assay, 200 reactions	\$ 350
	10031247	ddPCR MUT FAM assay, 1,000 reactions	\$ 1,375
	10031248	ddPCR MUT FAM assay, 2,500 reactions	\$ 1,875
	10031249	ddPCR MUT HEX assay, 200 reactions	\$ 350
	10031250	ddPCR MUT HEX assay, 1,000 reactions	\$ 1,375
	10031251	ddPCR MUT HEX assay, 2,500 reactions	\$ 1,875
Custom ddPCR Probe Assays	10031276	Custom ddPCR FAM assay, 500 reactions	\$ 175
	10031277	Custom ddPCR FAM assay, 1,000 reactions	\$ 225
	10031278	Custom ddPCR FAM assay, 2,500 reactions	\$ 350
	10031279	Custom ddPCR HEX assay, 500 reactions	\$ 200
	10031280	Custom ddPCR HEX assay, 1,000 reactions	\$ 300
	10031281	Custom ddPCR HEX assay, 2,500 reactions	\$ 400

Discount applies to bulk orders of the same catalog number and is taken off the list price.
Pricing as of November 2013. Prices may vary.

FAM is a trademark of Applera Corporation.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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