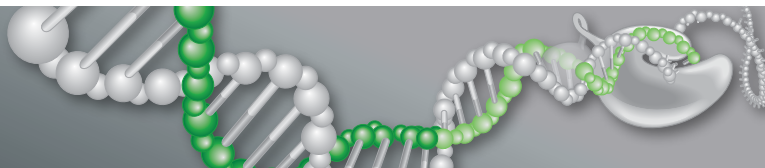


PrimePCR™ Pricing

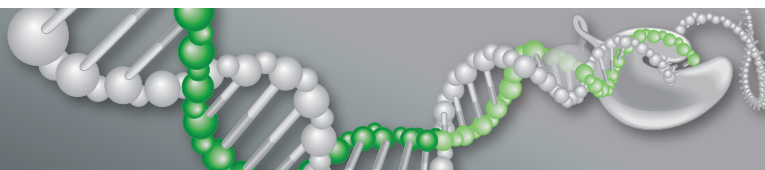


Assays and Controls

Primer Assay Pricing

Product	Catalog #	Description	List Price (JPY)
Primer Assays (desalted) (Wet-lab validated — human, mouse, and rat)	10025636	Primer assay desalted, 200 reactions	¥ 16,000
	10025637	Primer assay desalted, 1,000 reactions	¥ 48,000
	10025638	Primer assay desalted, 2,500 reactions	¥ 80,000
Primer Assays (HPLC) (Wet-lab validated — human, mouse, and rat)	10025639	Primer assay HPLC, 200 reactions	¥ 26,000
	10025640	Primer assay HPLC, 1,000 reactions	¥ 78,000
	10025641	Primer assay HPLC, 2,500 reactions	¥ 130,000
Primer Assays (desalted) (Expertly designed — <i>Arabidopsis</i> , chicken, cow, dog, pig, rabbit, rhesus monkey, yeast, zebrafish)	10042976	Primer assay desalted, 200 reactions	¥ 12,000
	10042977	Primer assay desalted, 1,000 reactions	¥ 36,000
	10042978	Primer assay desalted, 2,500 reactions	¥ 60,000
Primer Assays (HPLC) (Expertly designed — <i>Arabidopsis</i> , chicken, cow, dog, pig, rabbit, rhesus monkey, yeast, zebrafish)	10042979	Primer assay HPLC, 200 reactions	¥ 24,000
	10042980	Primer assay HPLC, 1,000 reactions	¥ 72,000
	10042981	Primer assay HPLC, 2,500 reactions	¥ 120,000
Custom Primer Pairs (desalted)	10025220	Custom primer pair desalted, 200 reactions	¥ 6,000
	10025221	Custom primer pair desalted, 1,000 reactions	¥ 16,000
	10025222	Custom primer pair desalted, 2,500 reactions	¥ 26,000
Custom Primer Pairs (HPLC)	10025223	Custom primer pair HPLC, 200 reactions	¥ 22,000
	10025224	Custom primer pair HPLC, 1,000 reactions	¥ 66,000
	10025225	Custom primer pair HPLC, 2,500 reactions	¥ 110,000
PreAmp Assays	10041594	PreAmp custom assay, 400 reactions	¥ 4,000
	10041595	PreAmp SYBR® Green assay, 400 reactions	¥ 16,000
	10041596	PreAmp probe assay, 400 reactions	¥ 16,000

Pricing as of September 2016. Prices may vary.



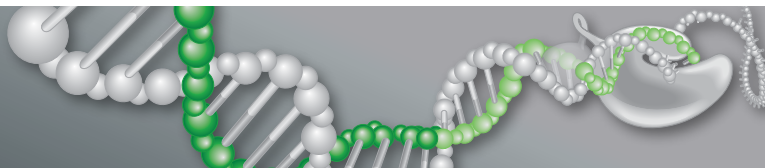
Assays and Controls

Probe Assay Pricing

Product	Catalog #	Description	List Price (JPY)
qPCR Probe Assays (Wet-lab validated — human, mouse, and rat)	12001950	qPCR FAM probe assay, 200 reactions	¥ 24,000
	10031225	qPCR FAM probe assay, 500 reactions	¥ 44,000
	10031226	qPCR FAM probe assay, 1,000 reactions	¥ 62,000
	10031227	qPCR FAM probe assay, 2,500 reactions	¥ 86,000
	10031228	qPCR HEX probe assay, 500 reactions	¥ 48,000
	10031229	qPCR HEX probe assay, 1,000 reactions	¥ 68,000
	10031230	qPCR HEX probe assay, 2,500 reactions	¥ 96,000
	10031234	qPCR TEX 615 probe assay, 500 reactions	¥ 48,000
	10031235	qPCR TEX 615 probe assay, 1,000 reactions	¥ 68,000
	10031236	qPCR TEX 615 probe assay, 2,500 reactions	¥ 96,000
	10031231	qPCR Cy5 probe assay, 500 reactions	¥ 48,000
	10031232	qPCR Cy5 probe assay, 1,000 reactions	¥ 68,000
	10031233	qPCR Cy5 probe assay, 2,500 reactions	¥ 96,000
	10031237	qPCR Cy5.5 probe assay, 500 reactions	¥ 48,000
	10031238	qPCR Cy5.5 probe assay, 1,000 reactions	¥ 68,000
	10031239	qPCR Cy5.5 probe assay, 2,500 reactions	¥ 96,000
	qPCR Probe Assays (Expertly designed — <i>Arabidopsis</i> , chicken, cow, dog, pig, rabbit, rhesus monkey, yeast, zebrafish)	12001961	qPCR FAM probe assay, 200 reactions
10042982		qPCR FAM probe assay, 500 reactions	¥ 34,000
10042983		qPCR FAM probe assay, 1,000 reactions	¥ 48,000
10042984		qPCR FAM probe assay, 2,500 reactions	¥ 68,000
10042985		qPCR HEX probe assay, 500 reactions	¥ 40,000
10042986		qPCR HEX probe assay, 1,000 reactions	¥ 56,000
10042987		qPCR HEX probe assay, 2,500 reactions	¥ 78,000
10042988		qPCR TEX 615 probe assay, 500 reactions	¥ 40,000
10042989		qPCR TEX 615 probe assay, 1,000 reactions	¥ 56,000
10042990		qPCR TEX 615 probe assay, 2,500 reactions	¥ 78,000
10042991		qPCR Cy5 probe assay, 500 reactions	¥ 78,000
10042992		qPCR Cy5 probe assay, 1,000 reactions	¥ 56,000
10042993		qPCR Cy5 probe assay, 2,500 reactions	¥ 78,000
10042994		qPCR Cy5.5 probe assay, 500 reactions	¥ 40,000
10042995		qPCR Cy5.5 probe assay, 1,000 reactions	¥ 56,000
10042996		qPCR Cy5.5 probe assay, 2,500 reactions	¥ 78,000

Pricing as of September 2016. Prices may vary.

continued

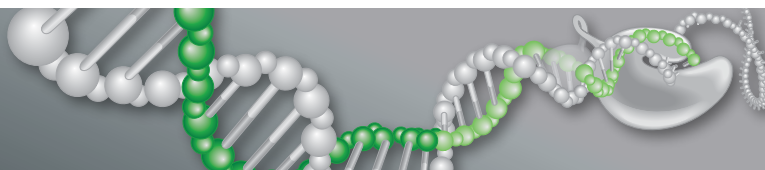


Assay and Controls

Probe Assay Pricing (continued)

Product	Catalog #	Description	List Price (JPY)
Custom qPCR Probe Assays	10031261	Custom qPCR FAM assay, 500 reactions	¥ 22,000
	10031262	Custom qPCR FAM assay, 1,000 reactions	¥ 32,000
	10031263	Custom qPCR FAM assay, 2,500 reactions	¥ 44,000
	10031264	Custom qPCR HEX assay, 500 reactions	¥ 30,000
	10031265	Custom qPCR HEX assay, 1,000 reactions	¥ 38,000
	10031266	Custom qPCR HEX assay, 2,500 reactions	¥ 52,000
	10031270	Custom qPCR TEX 615 assay, 500 reactions	¥ 34,000
	10031271	Custom qPCR TEX 615 assay, 1,000 reactions	¥ 48,000
	10031272	Custom qPCR TEX 615 assay, 2,500 reactions	¥ 68,000
	10031267	Custom qPCR Cy5 assay, 500 reactions	¥ 34,000
	10031268	Custom qPCR Cy5 assay, 1,000 reactions	¥ 48,000
	10031269	Custom qPCR Cy5 assay, 2,500 reactions	¥ 68,000
	10031273	Custom qPCR Cy5.5 assay, 500 reactions	¥ 34,000
	10031274	Custom qPCR Cy5.5 assay, 1,000 reactions	¥ 48,000
	10031275	Custom qPCR Cy5.5 assay, 2,500 reactions	¥ 68,000

Pricing as of September 2016. Prices may vary.



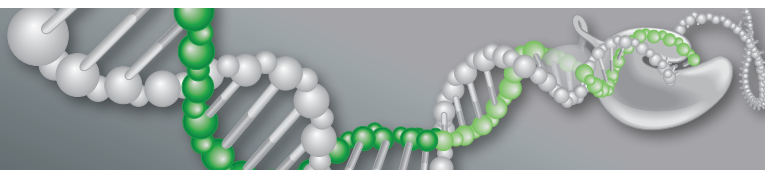
Template and Control Assays

Template and Control Assay Pricing

Product	Catalog #	Description	List Price (JPY)
Positive PCR and Reverse Transcription Control Assays — All Species	10025591	PCR control assay, 200 reactions	¥ 16,000
	10025695	RT control assay, 200 reactions	¥ 16,000
	10031288	PCR control FAM assay, 200 reactions	¥ 20,000
	10031289	PCR control HEX assay, 200 reactions	¥ 20,000
	10031286	RT control FAM assay, 200 reactions	¥ 20,000
Template and Control Assays — Human	10025716	Synthetic DNA control template, human	¥ 16,000
	10031285	Synthetic DNA control template probe, human	¥ 20,000
	10025694	RNA quality control assay, human	¥ 16,000
	10025352	gDNA control assay, human	¥ 16,000
	10031287	RNA quality control probe assay, human	¥ 20,000
	10031290	gDNA control probe assay, human	¥ 20,000
Template and Control Assays — Mouse	10029101	Synthetic DNA control template, mouse	¥ 16,000
	10031285	Synthetic DNA control template probe, mouse	¥ 20,000
	10029103	RNA quality control assay, mouse	¥ 16,000
	10029105	gDNA control assay, mouse	¥ 16,000
	10032858	RNA quality control probe assay, mouse	¥ 20,000
	10032859	gDNA control probe assay, mouse	¥ 20,000
Template and Control Assays — Rat	10044155	Synthetic DNA control template, rat	¥ 16,000
	10044156	Synthetic DNA control template probe, rat	¥ 20,000
	10044157	RNA quality control assay, rat	¥ 16,000
	10044158	gDNA control assay, rat	¥ 16,000
	10044159	RNA quality control probe assay, rat	¥ 20,000
	10044160	gDNA control probe assay, rat	¥ 20,000
Template and Control Assays — Zebrafish	10047261	Synthetic DNA control template, zebrafish	¥ 16,000
	10047262	Synthetic DNA control template probe, zebrafish	¥ 20,000
	10047263	RNA quality control assay, zebrafish	¥ 16,000
	10047264	gDNA control assay, zebrafish	¥ 16,000
	10047265	RNA quality control probe assay, zebrafish	¥ 20,000
	10047266	gDNA control probe assay, zebrafish	¥ 20,000
Template and Control Assays — Yeast	10047267	Synthetic DNA control template, yeast	¥ 16,000
	10047268	Synthetic DNA control template probe, yeast	¥ 20,000
	10047269	RNA quality control assay, yeast	¥ 16,000
	10047270	gDNA control assay, yeast	¥ 16,000
	10047271	RNA quality control probe assay, yeast	¥ 20,000
	10047272	gDNA control probe assay, yeast	¥ 20,000
Template and Control Assays — Rhesus Monkey	10047273	Synthetic DNA control template, rhesus monkey	¥ 16,000
	10047274	Synthetic DNA control template probe, rhesus monkey	¥ 20,000
	10047275	RNA quality control assay, rhesus monkey	¥ 16,000
	10047276	gDNA control assay, rhesus monkey	¥ 16,000
	10047277	RNA quality control probe assay, rhesus monkey	¥ 20,000
	10047278	gDNA control probe assay, rhesus monkey	¥ 20,000

Pricing as of September 2016. Prices may vary.

continued

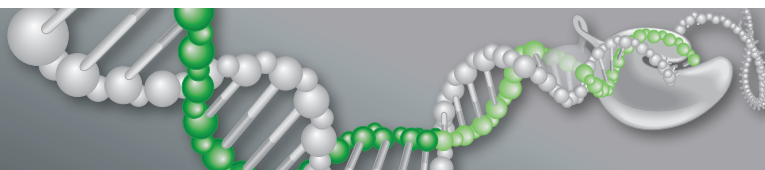


Template and Control Assays

Template and Control Assay Pricing (continued)

Product	Catalog #	Description	List Price (JPY)
Template and Control Assays — Pig	10047279	Synthetic DNA control template, pig	¥ 16,000
	10047280	Synthetic DNA control template probe, pig	¥ 20,000
	10047281	RNA quality control assay, pig	¥ 16,000
	10047282	gDNA control assay, pig	¥ 16,000
	10047283	RNA quality control probe assay, pig	¥ 20,000
	10047284	gDNA control probe assay, pig	¥ 20,000
Template and Control Assays — Cow	10047285	Synthetic DNA control template, cow	¥ 16,000
	10047286	Synthetic DNA control template probe, cow	¥ 20,000
	10047287	RNA quality control assay, cow	¥ 16,000
	10047288	gDNA control assay, cow	¥ 16,000
	10047289	RNA quality control probe assay, cow	¥ 20,000
	10047290	gDNA control probe assay, cow	¥ 20,000
Template and Control Assays — Rabbit	10047315	Synthetic DNA control template, rabbit	¥ 16,000
	10047316	Synthetic DNA control template probe, rabbit	¥ 20,000
	10047317	RNA quality control assay, rabbit	¥ 16,000
	10047318	gDNA control assay, rabbit	¥ 16,000
	10047319	RNA quality control probe assay, rabbit	¥ 20,000
	10047320	gDNA control probe assay, rabbit	¥ 20,000
Template and Control Assays — Dog	10047291	Synthetic DNA control template, dog	¥ 16,000
	10047292	Synthetic DNA control template probe, dog	¥ 20,000
	10047293	RNA quality control assay, dog	¥ 16,000
	10047294	gDNA control assay, dog	¥ 16,000
	10047295	RNA quality control probe assay, dog	¥ 20,000
	10047296	gDNA control probe assay, dog	¥ 20,000
Template and Control Assays — Chicken	10047297	Synthetic DNA control template, chicken	¥ 16,000
	10047298	Synthetic DNA control template probe, chicken	¥ 20,000
	10047299	RNA quality control assay, chicken	¥ 16,000
	10047300	gDNA control assay, chicken	¥ 16,000
	10047301	RNA quality control probe assay, chicken	¥ 20,000
	10047302	gDNA control probe assay, chicken	¥ 20,000
Template and Control Assays — <i>Arabidopsis</i>	10047303	Synthetic DNA control template, <i>Arabidopsis</i>	¥ 16,000
	10047304	Synthetic DNA control template probe, <i>Arabidopsis</i>	¥ 20,000
	10047305	RNA quality control assay, <i>Arabidopsis</i>	¥ 16,000
	10047306	gDNA control assay, <i>Arabidopsis</i>	¥ 16,000
	10047307	RNA quality control probe assay, <i>Arabidopsis</i>	¥ 20,000
	10047308	gDNA control probe assay, <i>Arabidopsis</i>	¥ 20,000

Pricing as of September 2016. Prices may vary.



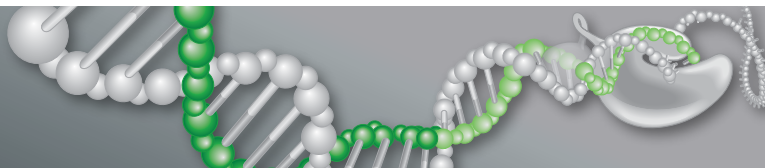
Pathway and Collection Arrays

Array Pricing

Product	Catalog #	Description	Minimum Plate Order Requirement	List Price (JPY)
96-Well PCR Arrays	96-Well Categories			
	Variable	16 assays x 6 samples	1	¥ 26,000
	Variable	24 assays x 4 samples	1	¥ 26,000
	Variable	32 assays x 3 samples	1	¥ 30,000
	Variable	48 assays x 2 samples	1	¥ 32,000
	Variable	96 assays x 1 sample	1	¥ 34,000
384-Well PCR Arrays	384-Well Categories			
	Variable	16 assays x 24 samples	1	¥ 38,000
	Variable	24 assays x 16 samples	1	¥ 40,000
	Variable	32 assays x 12 samples	1	¥ 42,000
	Variable	48 assays x 8 samples	1	¥ 42,000
	Variable	96 assays x 4 samples	1	¥ 44,000
	Variable	128 assays x 3 samples	1	¥ 48,000
	Variable	192 assays x 2 samples	1	¥ 50,000
	Variable	384 assays x 1 sample	1	¥ 54,000

Pricing as of September 2016. Prices may vary.

PrimePCR™ Pricing



Custom Plates

- Custom 96-well and 384-well plates with 1–96 unique genes
- Custom 384-well plates with 97 or more unique genes

Custom Plate Pricing

Product	Catalog #	Description	Minimum Plate Order Requirement	List Price (JPY)
Custom 96-Well Plates		96-Well Categories		
	10025216	1–12 genes	3	¥ 24,000
	10025217	13–24 genes	3	¥ 30,000
	10025218	25–36 genes	3	¥ 34,000
	10025219	37–96 genes	3	¥ 40,000
Custom 384-Well Plates		384-Well Categories		
	10025210	1–12 genes	6	¥ 34,000
	10025211	13–24 genes	6	¥ 40,000
	10025213	25–36 genes	6	¥ 46,000
	10025214	37–96 genes	6	¥ 52,000
	10025215	97–192 genes	12	¥ 56,000
	10025212	193–384 genes	12	¥ 68,000

Pricing as of September 2016. Prices may vary.

Cy is a trademark of GE Healthcare. FAM, HEX, SYBR, and TEX are trademarks of Life Technologies Corporation. Bio-Rad Laboratories, Inc. is licensed by Life Technologies Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

The use of PrimePCR PreAmp Assays is covered by one or more of the following U.S. patents and corresponding patent claims outside the U.S.: 5,804,375; 5,994,056; and 6,171,785. The purchase of these products includes a limited, nontransferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. These products are for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.



**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** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

