




Do You Have All You Need?

Bio-Rad has the right plastics for all your PCR needs

Amplification Plastics Selection Guide

		C1000™ or S1000™ Thermal Cycler	CFX96™ or CFX384™ Real-Time PCR Detection System
Tubes			
	TFI-0201	0.2 ml Tubes With Flat Caps, natural, 1,000	•
	TLS-0801	Low-Profile 0.2 ml 8-Tube Strips Without Caps, natural, 120 strips (960 tubes)	•
	TLS-0851	Low-Profile 0.2 ml 8-Tube Strips Without Caps, white, 120 strips (960 tubes)	•
48-Well Plates			
	MLL-4801	Multiplate™ Low-Profile 48-Well Unskirted PCR Plates, natural, 50	•
96-Well Plates			
	HSP-9601	Hard-Shell® Low-Profile 96-Well Skirted PCR Plates, white shell, clear well, 50	•
	HSP-9655	Hard-Shell Low-Profile 96-Well Skirted PCR Plates, white shell, white well, 50	•
	MLL-9601	Multiplate Low-Profile 96-Well Unskirted PCR Plates, natural, 25	•
	MLL-9651	Multiplate Low-Profile 96-Well Unskirted PCR Plates, white, 25	•
384-Well Plates			
	HSP-3801	Hard-Shell 384-Well Skirted PCR Plates, clear shell, clear well, 50	•
	HSP-3901	Hard-Shell 384-Well Skirted PCR Plates, clear shell, clear well, bar-coded, 50	•
	HSP-3805	Hard-Shell 384-Well Skirted PCR Plates, clear shell, white well, 50	•
	HSP-3905	Hard-Shell 384-Well Skirted PCR Plates, clear shell, white well, bar-coded, 50	•
	Sealers		
	MSB-1001	Microseal® 'B' Adhesive Seals, optically clear, 100	•
	TCS-0803	Optical Flat 8-Cap Strips, for 0.2 ml tubes and plates, ultraclear, 120 strips	•

For a complete list of all plastics, please visit www.bio-rad.com/pcrplastics.

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. SYBR is a trademark of Invitrogen Corporation. Bio-Rad Laboratories, Inc. is licensed by Invitrogen Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Hard-Shell plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 7,347,977, 6,340,589, and 6,528,302.

Notice regarding Bio-Rad thermal cyclers and real-time systems: Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under one or more of U.S. Patents 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160-163 and 167 only), and 6,703,236 (claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied Biosystems' United States Patent 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of these products includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.



Do You Have All You Need?

Bio-Rad has the right reagents for all your PCR needs

Amplification Reagent Selection Guide

Application	Reagents	Catalog #
High-Fidelity, Long, or Fast PCR	Hot-start iProof™ DNA polymerase	172-5351
	iProof™ high-fidelity DNA polymerase	172-5301
	iTaq™ DNA polymerase; dNTP mix	170-8870; 170-8874
Standard or Quantitative PCR		
Quantitative PCR		
Dye-based detection	SsoFast™ EvaGreen® supermix	172-5201
	iQ™ SYBR® Green supermix	170-8882
	iTaq™ SYBR® Green supermix with ROX	172-5851
	iTaq™ fast SYBR® Green supermix with ROX	172-5101
Probe-based detection	iQ supermix; iQ multiplex powermix	170-8862; 172-5849
	iTaq supermix with ROX	172-5855
	iTaq fast supermix with ROX	172-5106
One-Step Quantitative RT-PCR		
Dye-based detection	iScript™ one-step RT-PCR kit with SYBR® Green	170-8892
Probe-based detection	iScript one-step RT-PCR kit for probes	170-8894
Two-Step RT-PCR		
cDNA synthesis using a blend of random and oligo(dT) primers	iScript cDNA synthesis kit	170-8890
	Hot-start iProof DNA polymerase	172-5351
	iProof high-fidelity DNA polymerase	172-5301
	iTaq DNA polymerase; dNTP mix	170-8870; 170-8874
cDNA synthesis using random, oligo(dT), or gene-specific primers	iScript Select cDNA synthesis kit	170-8896
	Hot-start iProof DNA polymerase	172-5351
	iProof high-fidelity DNA polymerase	172-5301
	iTaq DNA polymerase; dNTP mix	170-8870; 170-8874
Two-Step Quantitative RT-PCR		
ROX-independent thermal cyclers	iScript Select cDNA synthesis kit	170-8896
	iScript cDNA synthesis kit	170-8890
	iQ supermix; iQ multiplex powermix	170-8862; 172-5849
	iQ™ SYBR® Green supermix	170-8882
ROX-dependent thermal cyclers	iScript Select cDNA synthesis kit	170-8896
	iScript cDNA synthesis kit	170-8890
	iTaq supermix with ROX	172-5855
	iTaq™ SYBR® Green supermix with ROX	172-5851
	iTaq fast supermix with ROX	172-5106
	iTaq™ fast SYBR® Green supermix with ROX	172-5101
	iQ multiplex powermix	172-5849
	ROX passive reference dye	172-5858

To learn more about our products and to choose the product size that's just right for you, please visit us on the Web at www.bio-rad.com/pcrreagents.