

Do You Have All You Need?

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

Amplification Plastics Selection Guide

		C1000 Touch™ or S1000™ Thermal Cycler	CFX96 Touch™, CFX96 Touch Deep Well, or CFX Connect™ Real-Time PCR Detection System	CFX384 Touch™ Real-Time PCR Detection System
Tubes				
TFI-0201	PCR Tubes with Flat Caps (0.2 ml), clear, 1,000	•		
TLS-0801	Low-Profile 8-Tube Strips without Caps (0.2 ml), clear, 120 strips (960 PCR tubes)	•	•	
TLS-0851	Low-Profile 8-Tube Strips without Caps (0.2 ml), white, 120 strips (960 PCR tubes)		•	
48-Well Plates				
MLL-4801	Multiplate™ Low-Profile 48-Well Unskirted PCR Plates, clear, 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, clear, 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, barcoded, 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, barcoded, 50			•
Sealers				
MSB-1001	Microseal® 'B' Adhesive Seals, optically clear, 100	•	•	•
MSC-1001	Microseal 'C' Optical Seals, 100	•	•	•
TCS-0803	Optical Flat 8-Cap Strips, for 0.2 ml PCR tubes and plates, ultraclear, 120 strips	•	•	
PCR Plate Sealer and Heat Seals				
181-4000	PX1™ PCR Plate Sealer, includes heat sealing instrument, plate support block that holds 96-well and 384-well plates, sealing frame, power cord			
181-4030	Optically Clear Heat Seal, 100	•	•	•
181-4040	Pierceable Foil Heat Seal, 100	•		

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

SYBR is a trademark 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.

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 Numbers 6,767,512 and 7,074,367.

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of iProof, iQ, iTag, and SsoAdvanced 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.

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 Numbers 7,347,977; 6,340,589; and 6,528,302.

BIO-RAD

Do You Have All You Need?

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

Amplification Reagents Selection Guide

Catalog #	Description	Catalog #	Description
Two-Step Reverse Transcription Reagents		iTaq Universal Probes Supermixes	
170-8842	iScript™ Advanced cDNA Synthesis Kit for RT-qPCR, 50 x 20 µl reactions	172-5130	iTaq Universal Probes Supermix, 200 x 20 µl reactions
170-8843	iScript Advanced cDNA Synthesis Kit for RT-qPCR, 250 x 20 µl reactions	172-5131	iTaq Universal Probes Supermix, 500 x 20 µl reactions
170-8840	iScript Reverse Transcription Supermix for RT-qPCR, 25 x 20 µl reactions	172-5132	iTaq Universal Probes Supermix, 1,000 x 20 µl reactions
170-8841	iScript Reverse Transcription Supermix for RT-qPCR, 100 x 20 µl reactions	172-5134	iTaq Universal Probes Supermix, 2,500 x 20 µl reactions
170-8890	iScript cDNA Synthesis Kit, 25 x 20 µl reactions	172-5135	iTaq Universal Probes Supermix, 5,000 x 20 µl reactions
170-8891	iScript cDNA Synthesis Kit, 100 x 20 µl reactions		
170-8896	iScript Select cDNA Synthesis Kit, 25 x 20 µl reactions	High Resolution Melt Analysis Supermixes	
170-8897	iScript Select cDNA Synthesis Kit, 100 x 20 µl reactions	172-5110	Precision Melt Supermix, 200 x 20 µl reactions
		172-5112	Precision Melt Supermix, 1,000 x 20 µl reactions
SsoAdvanced™ Universal SYBR® Green Supermixes		Multiplex Supermixes	
172-5270	SsoAdvanced™ Universal SYBR® Green Supermix, 200 x 20 µl reactions	172-5848	iQ™ Multiplex Powermix, 50 x 50 µl reactions
172-5271	SsoAdvanced™ Universal SYBR® Green Supermix, 500 x 20 µl reactions	172-5849	iQ Multiplex Powermix, 200 x 50 µl reactions
172-5272	SsoAdvanced™ Universal SYBR® Green Supermix, 1,000 x 20 µl reactions		
172-5274	SsoAdvanced™ Universal SYBR® Green Supermix, 2,500 x 20 µl reactions	Standard PCR Supermixes*	
172-5275	SsoAdvanced™ Universal SYBR® Green Supermix, 5,000 x 20 µl reactions	172-5301	iProof™ High-Fidelity DNA Polymerase, 2 U/µl, 100 U
		172-5302	iProof High-Fidelity DNA Polymerase, 2 U/µl, 500 U
SsoAdvanced Universal Probes Supermixes		170-8870	iTaq DNA Polymerase, 5 U/µl, 250 U
172-5280	SsoAdvanced Universal Probes Supermix, 200 x 20 µl reactions	170-8875	iTaq DNA Polymerase, 5 U/µl, 5,000 U
172-5281	SsoAdvanced Universal Probes Supermix, 500 x 20 µl reactions	170-8874	dNTP Mix, 200 µl
172-5282	SsoAdvanced Universal Probes Supermix, 1,000 x 20 µl reactions	172-5858	ROX Passive Reference Dye, 0.5 ml
172-5284	SsoAdvanced Universal Probes Supermix, 2,500 x 20 µl reactions		
172-5285	SsoAdvanced Universal Probes Supermix, 5,000 x 20 µl reactions		
iTaq™ Universal SYBR® Green Supermixes			
172-5120	iTaq™ Universal SYBR® Green Supermix, 200 x 20 µl reactions		
172-5121	iTaq™ Universal SYBR® Green Supermix, 500 x 20 µl reactions		
172-5122	iTaq™ Universal SYBR® Green Supermix, 1,000 x 20 µl reactions		
172-5124	iTaq™ Universal SYBR® Green Supermix, 2,500 x 20 µl reactions		
172-5125	iTaq™ Universal SYBR® Green Supermix, 5,000 x 20 µl reactions		

* For more product size options, or to learn more about our PCR reagents, visit www.bio-rad.com/web/AmplificationReagents.