

# Do You Have All You Need?

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



## Amplification Plastics Selection Guide

Catalog #	Description	C1000 Touch or S1000 Thermal Cycler	CFX Opus 96, CFX96 Touch, CFX96 Touch Deep Well, or CFX Connect Real-Time PCR System	CFX Opus 384 or CFX384 Touch Real-Time PCR System
<b>Tubes and Caps</b>				
TLS0801	<b>Low-Profile 8-Tube Strips without Caps (0.2 ml)</b> , clear, 120 strips (960 PCR tubes)	•	•	
TLS0851	<b>Low-Profile 8-Tube Strips without Caps (0.2 ml)</b> , white, 120 strips (960 PCR tubes)		•	
TCS0803	<b>Optical Flat 8-Cap Strips</b> , for 0.2 ml PCR tubes and plates, ultraclear, 120 strips	•	•	
<b>48-Well Plates</b>				
MLL4801	<b>Multiplate Low-Profile 48-Well Unskirted PCR Plates</b> , clear, 50	•		
<b>96-Well Plates</b>				
HSP9601	<b>Hard-Shell Low-Profile 96-Well Skirted PCR Plates</b> , white shell, clear well, 50	•	•	
HSP9655	<b>Hard-Shell Low-Profile 96-Well Skirted PCR Plates</b> , white shell, white well, 50	•	•	
<b>384-Well Plates</b>				
HSP3801	<b>Hard-Shell 384-Well Skirted PCR Plates</b> , clear shell, clear well, 50	•		
HSP3901	<b>Hard-Shell 384-Well Skirted PCR Plates</b> , clear shell, clear well, barcoded, 50	•		
HSP3805	<b>Hard-Shell 384-Well Skirted PCR Plates</b> , clear shell, white well, 50			•
HSP3905	<b>Hard-Shell 384-Well Skirted PCR Plates</b> , clear shell, white well, barcoded, 50			•
<b>Adhesive Plate Seals</b>				
MSB1001	<b>Microseal 'B' PCR Plate Sealing Film</b> , optically clear seals, 100	•	•	•
MSC1001	<b>Microseal 'C' PCR Plate Sealing Film</b> , optically clear seals, 100	•	•	•
<b>PCR Plate Sealer and Heat Seals</b>				
1814000	<b>PX1 PCR Plate Sealer</b>	•	•	•
1814030	<b>Optically Clear Heat Seal</b> , 100	•	•	•
1814040	<b>Pierceable Foil Heat Seal</b> , 100	•		

Visit [bio-rad.com/AmplificationPlastics](http://bio-rad.com/AmplificationPlastics) for a complete list of all plastics.

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

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.

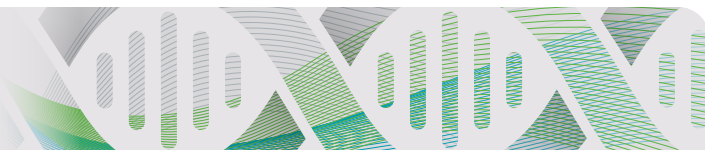
SYBR is a trademark of Thermo Fisher Scientific Inc. Bio-Rad Laboratories, Inc. is licensed by Thermo Fisher Scientific Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

All trademarks used herein are the property of their respective owner.



# Do You Have All You Need?

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



## Amplification Reagents Selection Guide

Catalog #	Description	Catalog #	Description
<b>Two-Step Reverse Transcription (RT) Reagents</b>		<b>Cell Lysis RT-qPCR Kits</b>	
1725037	iScript Advanced cDNA Synthesis Kit for RT-qPCR, 25 x 20 µl reactions	1725080	SingleShot Cell Lysis Kit, 100 x 50 µl reactions
1725038	iScript Advanced cDNA Synthesis Kit for RT-qPCR, 100 x 20 µl reactions	1725085	SingleShot SYBR® Green Kit for Cell Lysis and RT-qPCR, 100 x 50 µl reactions
1708840	iScript Reverse Transcription Supermix for RT-qPCR, 25 x 20 µl reactions	1725090	SingleShot Probes Kit for Cell Lysis and RT-qPCR, 100 x 50 µl reactions
1708841	iScript Reverse Transcription Supermix for RT-qPCR, 100 x 20 µl reactions	1725095	SingleShot SYBR® Green One-Step Kit for Cell Lysis and RT-qPCR, 100 x 50 µl reactions
1708890	iScript cDNA Synthesis Kit, 25 x 20 µl reactions	1725070	SingleShot Probes One-Step Kit for Cell Lysis and RT-qPCR, 100 x 50 µl reactions
1708891	iScript cDNA Synthesis Kit, 100 x 20 µl reactions		
1708896	iScript Select cDNA Synthesis Kit, 25 x 20 µl reactions		
1708897	iScript Select cDNA Synthesis Kit, 100 x 20 µl reactions		
<b>Quantitative PCR (qPCR) Supermixes with SYBR® Green</b>		<b>Multiplex Supermixes</b>	
1725270	SsoAdvanced Universal SYBR® Green Supermix, 200 x 20 µl reactions	1725848	iQ Multiplex Powermix, 50 x 50 µl reactions
1725271	SsoAdvanced Universal SYBR® Green Supermix, 500 x 20 µl reactions	1725849	iQ Multiplex Powermix, 200 x 50 µl reactions
1725016	SsoAdvanced Universal Inhibitor-Tolerant SYBR® Green Supermix, 200 x 20 µl reactions	12010176	Reliance One-Step Multiplex Supermix, 200 x 20 µl reactions
1725017	SsoAdvanced Universal Inhibitor-Tolerant SYBR® Green Supermix, 500 x 20 µl reactions	12010220	Reliance One-Step Multiplex Supermix, 1,000 x 20 µl reactions
1725120	iTaq Universal SYBR® Green Supermix, 200 x 20 µl reactions	<b>Preamplification Reagents</b>	
1725121	iTaq Universal SYBR® Green Supermix, 500 x 20 µl reactions	1725160	SsoAdvanced PreAmp Supermix, 50 x 50 µl reactions
<b>Probe-Based qPCR Supermixes</b>		12004856	iScript Explore One-Step RT and PreAmp Kit, 50 x 50 µl reactions
1725280	SsoAdvanced Universal Probes Supermix, 200 x 20 µl reactions	<b>High Resolution Melt Analysis Supermixes</b>	
1725281	SsoAdvanced Universal Probes Supermix, 500 x 20 µl reactions	1725110	Precision Melt Supermix, 200 x 20 µl reactions
1725130	iTaq Universal Probes Supermix, 200 x 20 µl reactions	1725112	Precision Melt Supermix, 1,000 x 20 µl reactions
1725131	iTaq Universal Probes Supermix, 500 x 20 µl reactions	<b>Standard PCR Reagents and Kits</b>	
<b>One-Step Reverse Transcription qPCR (RT-qPCR) Reagents</b>		1725310	iProof HF Master Mix, 100 x 50 µl reactions
12010176	Reliance One-Step Multiplex Supermix, 200 x 20 µl reactions	1725301	iProof High-Fidelity DNA Polymerase, 2 U/µl, 100 U
1725150	iTaq Universal SYBR® Green One-Step Kit, 100 x 20 µl reactions	1725330	iProof High-Fidelity PCR Kit, 2 U/µl, 50 U
1725140	iTaq Universal Probes One-Step Kit, 100 x 20 µl reactions	1708870	iTaq DNA Polymerase, 5 U/µl, 250 U
		1708874	dNTP Solution Mix, 200 µl

Visit [bio-rad.com/AmplificationReagents](http://bio-rad.com/AmplificationReagents) for more product size options, or to find the best reagents for your application.



Bio-Rad's PCR reagents and analytical instruments are manufactured under an ISO 13485:2016 certified Quality Management System and are quality control tested to ensure consistent product performance you can trust.

**BIO-RAD**