

Do You Have All You Need?

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

Amplification Plastics Selection Guide for the T100™ Thermal Cycler



Individual Tubes

TFI-0201 **PCR Tubes with Flat Caps (0.2 ml)**, clear, 1,000
TWI-0201 **PCR Tubes with Domed Caps (0.2 ml)**, clear, 1,000



Tube Strips

TBC-0802 **8-Tube Strips and Domed Cap Strips (0.2 ml)**, clear, 20 bags of 12 x 8-tube strips and 12 x 8-cap strips (1,920 PCR tubes and 1,920 caps)
TBC-1202 **12-Tube Strips and Domed Cap Strips (0.2 ml)**, clear, 20 bags of 8 x 12-tube strips and 8 x 12-cap strips (1,920 PCR tubes and 1,920 caps)
TBS-0201 **8-Tube Strips without Caps (0.2 ml)**, clear, 120 strips (960 PCR tubes)
TBS-1201 **12-Tube Strips without Caps (0.2 ml)**, clear, 100 strips (1,200 PCR tubes)

96-Well Plates

MLP-9601 **Multiplate™ High-Profile 96-Well Unskirted PCR Plates**, clear, 25 plates
HSS-9601 **Hard-Shell® High-Profile 96-Well Semi-Skirted PCR Plates**, clear shell, clear well, 25
HSS-9901 **Hard-Shell High-Profile 96-Well Semi-Skirted PCR Plates**, clear shell, clear well, barcoded, 25
HSS-9641 **Hard-Shell High-Profile 96-Well Semi-Skirted PCR Plates**, green shell, clear well, 25
223-9441 **iQ™ 96-Well PCR Plates**, 25

Sealers

MSA-5001 **Microseal® 'A' Film**, package of 50 seals
MSB-1001 **Microseal 'B' Adhesive Seals**, optically clear, 100
MSF-1001 **Microseal 'F' Foil**, package of 100 seals
TCS-0803 **Optical Flat 8-Cap Strips**, for 0.2 ml PCR tubes and plates, ultraclear, 120
TCS-0801 **Domed 8-Cap Strips**, for 0.2 ml PCR tubes and plates, clear, 120
TCS-1201 **Domed 12-Cap Strips**, for 0.2 ml PCR tubes and plates, clear, 200
223-9442 **96-Well PCR Plate Sealing Mats**, 5

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**, package of 100 optically clear seals
181-4035 **Permanent Clear Heat Seal**, package of 100 nonpeelable, nonpierceable clear seals
181-4040 **Pierceable Foil Heat Seal**, package of 100 foil seals
181-4045 **Peelable Foil Heat Seal**, package of 100 laminate/foil peelable seals

For a complete list of all plastics, visit www.bio-rad.com/web/T100PCRplastics.

Bio-Rad's 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 and iTaq polymerases 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 for the T100™ Thermal Cycler

Application	Reagent	Catalog #	
High-Fidelity, Long, or Fast PCR	iProof™ High-Fidelity DNA Polymerase, 2 U/μl, 100 U	172-5301	
	iProof High-Fidelity DNA Polymerase, 2 U/μl, 500 U	172-5302	
	iProof HF Master Mix, 100 x 50 μl reactions	172-5310	
	iProof GC Master Mix, 100 x 50 μl reactions	172-5320	
	iProof High-Fidelity PCR Kit, 2 U/μl, 50 U, includes 5x reaction buffers, MgCl ₂ solution, DMSO, dNTPs, λ DNA, 1.3 and 10 kb primers, DNA standard	172-5330	
	iProof High-Fidelity PCR Kit, 2 U/μl, 200 U	172-5331	
PCR and Reverse Transcription PCR	iTaq™ DNA Polymerase, 5 U/μl, 250 U	170-8870	
	iTaq DNA Polymerase, 5 U/μl, 5,000 U	170-8875	
	dNTP Mix, 200 μl	170-8874	
	cDNA synthesis using a blend of random and oligo(dT) primers	iScript™ Advanced cDNA Synthesis Kit for RT-qPCR, 50 x 20 μl reactions	170-8842
		iScript Advanced cDNA Synthesis Kit for RT-qPCR, 250 x 20 μl reactions	170-8843
		iScript Reverse Transcription Supermix for RT-qPCR, 25 x 20 μl reactions	170-8840
		iScript Reverse Transcription Supermix for RT-qPCR, 100 x 20 μl reactions	170-8841
	cDNA synthesis using random, oligo(dT), or gene-specific primers	iScript cDNA Synthesis Kit, 25 x 20 μl reactions	170-8890
		iScript cDNA Synthesis Kit, 100 x 20 μl reactions	170-8891
		iScript cDNA Synthesis Kit, 500 x 20 μl reactions	170-8891BUN
iScript Select cDNA Synthesis Kit, 25 x 20 μl reactions		170-8896	
iScript Select cDNA Synthesis Kit, 100 x 20 μl reactions		170-8897	
iScript Select cDNA Synthesis Kit, 500 x 20 μl reactions		170-8897BUN	

To learn more about our products and to choose the product size that's just right for you, visit www.bio-rad.com/web/T100PCRreagents.

