



PCR Plastics That Perform: Every Run, Every Workflow

Everything You Need for PCR and Real-Time PCR

Bio-Rad's complete line of PCR plates, seals, tubes, and accessories support a wide range of molecular applications by delivering:

- **Precision you can count on:** high-quality injection molding delivers consistent fit and reliable performance
- **Built-in traceability for confidence:** lot-to-lot batch management supports quality control for regulated or diagnostic use
- **Automation readiness:** seamless transition from manual to automated workflows without compromising PCR performance

Use the table below to identify compatible consumables for your Bio-Rad PCR and real-time PCR systems.

Table 1. Instrument compatibility for supported Bio-Rad PCR and real-time PCR systems.

Hard-Shell™ Plates and Multiplates	96- and 48-Well Plates						384-Well Plates
	Hard-Shell Semi-Skirted High-Profile	Hard-Shell Skirted Low-Profile	Hard-Shell Semi-Skirted Low-Profile	Multiplate Unskirted High-Profile	Multiplate Unskirted Low-Profile	iQ Semi- Skirted High-Profile	Hard-Shell Skirted Standard
System #	HSS9xxx	HSP9xxx	HSL9xxx	MLPxxx	MLLxxx	2239441	HSP3xxx
PTC Harmony and PTC Harmony Deepwell Thermal Cyclers	✓	✓	✓	✓	✓	✓	
PTC Tempo and PTC Tempo Deepwell Thermal Cyclers	✓	✓	✓	✓	✓	✓	✓
T100 Thermal Cycler	✓			†		†	
C1000, C1000 Touch, S1000 Thermal Cyclers	†	✓	†	†	†	†	✓
CFX Opus, CFX Touch, CFX Connect and CFX Duet Real-Time PCR Systems	†	✓	†	†	✓	†	✓
CFX Opus and CFX Touch Deepwell Real-Time PCR System	✓	✓	†	✓	✓	†	

✓ Recommended: full validation passed. † Compatible: fit/visual inspection only.



Table 1. Instrument compatibility for supported Bio-Rad PCR and real-time PCR systems (continued).

Tubes					Cap Strips			
PCR Tubes and Caps	Individual High Profile Tube With Flat Cap	Individual High Profile Tube With Doomed Cap	Strips, High Profile	Strips, Low-Profile	Doomed 8-Strips	Doomed 12-Strips	Flat 8-Cap Strips, Optical	
System #	TFI0201	TWI0201	TBSxxxx, TBCxxxx 8-strips, 12-Strips	TLSxxxx	TCS0801	TCS1201	TCS0803	
PTC Harmony and PTC Harmony Deepwell Thermal Cyclers	✓	✓	‡	✓	‡	‡	✓	
PTC Tempo and PTC Tempo Deepwell Thermal Cyclers	✓	✓	‡	✓	‡	‡	✓	
T100 Thermal Cycler	✓	✓	✓		‡	‡	‡	
C1000, C1000 Touch, S1000 Thermal Cyclers	✓	✓	‡	✓	‡	‡	‡	
CFX Opus, CFX Touch, CFX Connect and CFX Duet Real-Time PCR Systems				✓			✓	
CFX Opus and CFX Touch Deepwell Real-Time PCR System	✓		✓	✓			✓	
MicroSeal™ PCR Sealing Film				PCR Plate Heat Seal				
Sealing Films and Heat Seals	Microseal A	Microseal B	Microseal C	Microseal F	Optically Clear	Clear, Permanent	Pierceable Foil	Peelable Foil
System #	MSA5001	MSB1001B	MSC1001	MSF1001B	1814030	1814035	1814040	1814045
T100, C1000, C1000 Touch, S1000 PTC Tempo	‡	✓	✓	✓	✓	✓	✓	✓
PTC Tempo Deepwell, PTC Harmony, and PTC Harmony Deepwell Thermal Cyclers	✓	✓	✓	✓	✓	✓	✓	✓
CFX Opus and Deepwell, CFX Touch and Deepwell, CFX Connect and CFX Duet Real-Time PCR Systems		✓	✓		✓	✓		

✓ Recommended: full validation passed. ‡ Compatible: fit/visual inspection only.

Bio-Rad, Hard-Shell, and Microseal are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.
© 2026 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 55 11 3065 7550 **Canada** 1 800 361 1808 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Greece** 30 210 7774396 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 000 800 00 24 67 23 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 55 5488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Russian Federation** 7 495 721 14 04 **Singapore** 65 6415-3170 **South Africa** 27 21 531 7504 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 662 651 8311 **United Arab Emirates** 971 4 818 7300 **United Kingdom** 00 800 00 24 67 23