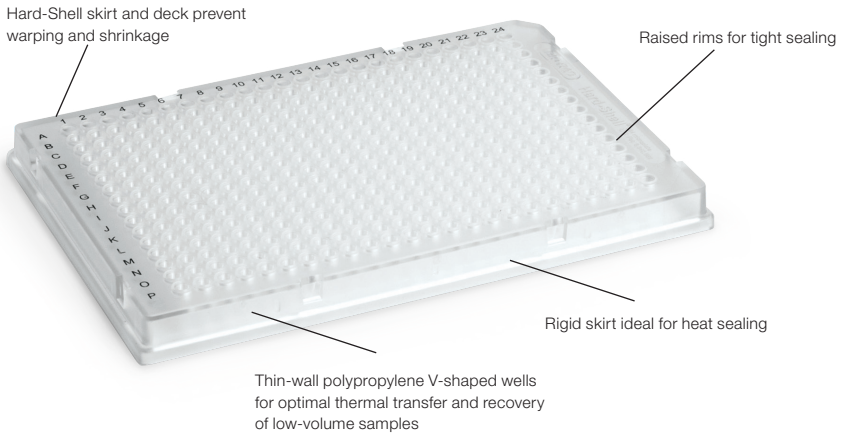


Hard-Shell® 384-Well Standard PCR Plates

Hard-Shell PCR plates represent an innovative concept in PCR plate design. The skirt and deck are molded from a rigid thermostable polymer, preventing robotic handling problems due to plate warping. This hard shell prevents the distortion and shrinkage that may occur when regular single-component polypropylene PCR plates are exposed to the high temperatures of thermal cycling or heat sealing. The thin-wall wells are molded of virgin polypropylene selected for low DNA binding. The raised rims around each well allow tight sealing using a variety of procedures, including pressure-, adhesive-, and heat-sealing methods.



BIO-RAD

Hard-Shell 384-Well Standard PCR Plates

16- x 24-well array, package of 50 plates

HSP-3805	White well, clear shell
HSP-3905	White well, clear shell, bar-coded on 2 sides (Row P and Column 24)
HSP-3865	White well, black shell
HSP-3801	Clear well, clear shell
HSP-3901	Clear well, clear shell, bar-coded on 2 sides (Row P and Column 24)
HSP-3811	Clear well, red shell
HSP-3821	Clear well, yellow shell
HSP-3831	Clear well, blue shell
HSP-3841	Clear well, green shell
HSP-3851	Clear well, white shell
HSP-3866	Black well, black shell

Seals and Accessories

181-4000	PX1™ PCR Plate Sealer , semiautomated heat sealing instrument
181-4030	Optically Clear Heat Seal , for use with PX1 PCR plate sealer, 100 seals
MSA-5001	Microseal® 'A' Film , 50 seals
MSB-1001	Microseal 'B' Adhesive Seals , optically clear, 100 seals
MSC-1001	Microseal 'C' Optical Seals , 100 adhesive seals
MSF-1001	Microseal 'F' Foil , 100 adhesive seals
MSR-0001	Sealing Roller , for adhesive film seals
CHO-1401	Chill-out™ Liquid Wax , red, 100 ml
CHO-1411	Chill-out Liquid Wax , clear, optical grade, 100 ml

Please request bulletin 5496 for more information about Hard-Shell PCR plates, including dimension details, and bulletin 6090 for more information about seals and accessories.

Hard-Shell PCR Plate Compatibility Chart

384-Well Standard PCR Plates



Catalog number	HSP-3xxx
Maximum well volume	50 μ l
ANSI/SBS standard dimensions	Yes
Autoclave specification	121°C, 20 min
Centrifuge specification	Up to 2,250 x g in a swing-out rotor
Application temperature	-80 to 121°C
Storage	Room temperature; avoid sun and UV light
Cleanroom production	Free of DNase, RNase, and human DNA
Thermal Cycler Compatibility*	
Bio-Rad® C1000™, C1000 Touch™, S1000™	•
Bio-Rad® DNA Engine®, DNA Engine Tetrad® and Tetrad 2, DNA Engine Dyad®, Dyad Disciple™	•
Applied Biosystems cyclers (9700, Veriti, etc.)	•
Eppendorf Mastercycler series	•
Real-Time PCR Detection System Compatibility*	
Bio-Rad® CFX384™, CFX384 Touch™	•
Applied Biosystems systems (7900HT, ViiA 7)	•
Other Instrument Compatibility*	
Applied Biosystems capillary DNA sequencers (3700, 3730, etc.)	•
Idaho Technology LightScanner	•

* 384-well instruments.

Note: The Hard-Shell 384-well 480 PCR plates have been specifically designed for the Roche LightCycler 480. Please request bulletin 5496 for more information about these plates.

Eppendorf and Mastercycler are trademarks of Eppendorf AG. LightCycler is a trademark of Roche Diagnostics GmbH. LightScanner is a trademark of Idaho Technology Inc. Veriti is a trademark of Applied Biosystems Corporation. ViiA is a trademark of Life Technologies Corporation.

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.

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 U.S. Patent Number 5,475,610 (Claims 1, 44, 158, 160–163, and 167 only), or corresponding claims in its non-U.S. counterpart, owned by Applied Biosystems Corporation. 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' U.S. Patent Number 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 Numbers 6,767,512 and 7,074,367.



**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site www.bio-rad.com **USA** 800 424 6723
Australia 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11
Brazil 55 11 5044 5699 **Canada** 905 364 3435 **China** 86 21 6169 8500
Czech Republic 420 241 430 532 **Denmark** 44 52 10 00
Finland 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0
Greece 30 210 9532 220 **Hong Kong** 852 2789 3300
Hungary 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460
Mexico 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30
Poland 48 22 331 99 99 **Portugal** 351 21 472 7700
Russia 7 495 721 14 04 **Singapore** 65 6415 3188
South Africa 27 861 246 723 **Spain** 34 91 590 5200
Sweden 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000