

Customize Hard-Shell™ PCR Plates

Innovative Ideas Require Unique Solutions

Imagine a PCR Plate Manufactured to Meet Your Precise Needs

Do you need a PCR plate with a specific barcode, with your logo engraved on the plate, or packaged in a specific configuration? Bio-Rad Laboratories, Inc. is ready to partner with you to customize our popular Hard-Shell PCR Plates to meet your requirements.

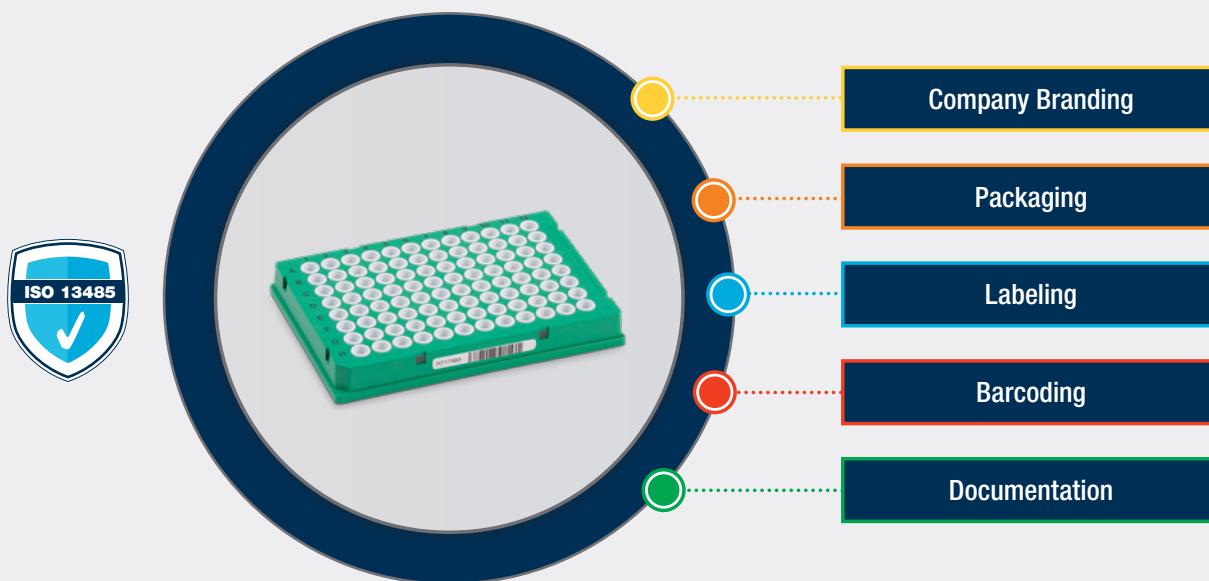
With more than 70 years of experience developing and manufacturing products for life sciences and clinical diagnostics customers, Bio-Rad has the knowledge and expertise to produce your custom PCR plate. As your custom manufacturing partner, we are committed to delivering exemplary products designed to your unique specifications.

Our global manufacturing centers of excellence follow International Quality Standards (ISO 13485:2016) and stringent quality controls to guarantee the consistent, reproducible, and quality Hard-Shell PCR Plates you have come to expect from Bio-Rad.

Custom Process

Our knowledgeable team will be your guide through each step of the process. From the first inquiry and consultation to the first article inspection and production scale-up, and throughout the partnership, our team of experts is dedicated to ensuring your custom PCR plate meets your specific needs on your timeline.

Custom Options



Custom Options for Hard-Shell PCR Plates

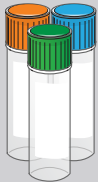
Branding — Include your company name and logo on PCR plates and packaging labels.

Packaging — Select from various box sizes to fit the exact number of plates you require. Determine the packaging format with the number of plates per bag, the number of bags per box, and other specifications.

Labeling — Customize your packaging labels to meet your product specifications, including size, location, and format.

Barcoding — Specify the number, placement, and sequence of one-dimensional or two-dimensional barcodes on your PCR plate.

Documentation — Add product insert and other documentation to your customized product.



Did you know? You can also customize Bio-Rad PCR reagents and kits for reliable amplification, robust reverse transcription, and high-quality real-time PCR data.

Visit www.biorad.com/CustomReagents to learn more.

Shell Description	Clear Well	White Well
Hard-Shell Low-Profile 96-Well Skirted PCR Plates, 50 plates		
White shell	HSP9601	HSP9655
Red shell	HSP9611	—
Yellow shell	HSP9621	—
Blue shell	HSP9631	HSP9635
Green shell	HSP9641	HSP9645
White shell, barcoded	HSP9901	HSP9955

Hard-Shell Low-Profile 96-Well Semi-Skirted PCR Plates, 25 plates		
Clear shell	HSL9601	HSL9605
Green shell	HSL9641	HSL9645
Clear shell, barcoded	HSL9901	HSL9905

Hard-Shell High-Profile 96-Well Semi-Skirted PCR Plates, 25 plates		
Clear shell	HSS9601	—
Green shell	HSS9641	—
Clear shell, barcoded	HSS9901	—

Hard-Shell 384-Well Standard PCR Plates, 50 plates		
Clear shell	HSP3801	HSP3805
Red shell	HSP3811	—
Yellow shell	HSP3821	—
Blue shell	HSP3831	—
Green shell	HSP3841	—
Clear shell, barcoded	HSP3901	HSP3905

Hard-Shell 96-Well 480 PCR Plates, with barcode on Row A side, 25 plates		
Clear shell	—	HSR9905
Clear shell	HSR9901	—

Hard-Shell 384-Well 480 PCR Plates, with barcode on Row A and Column 24 sides, 50 plates		
Clear shell	—	HSR4805
Clear shell	HSR4801	—

High-profile = 20.75 mm
Low-profile = 16.06 mm

Visit www.bio-rad.com/CustomPlastics to initiate the process. A custom production expert will contact you to get your project started.

BIO-RAD and HARD-SHELL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 by Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](http://www.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** 4003 0399 **Canada** 1 905 364 3435 **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 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050 **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** 52 555 488 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** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **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** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

