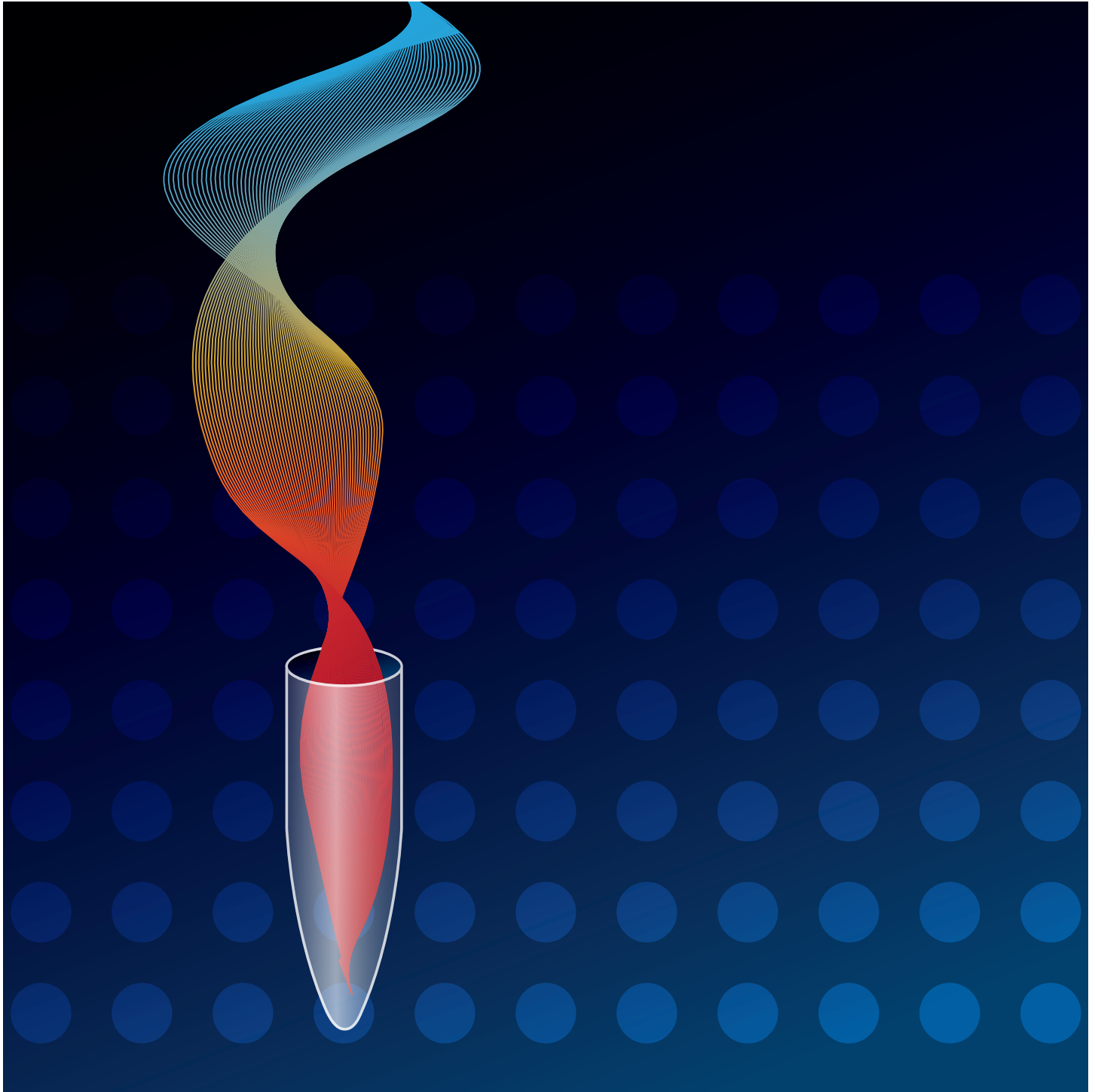
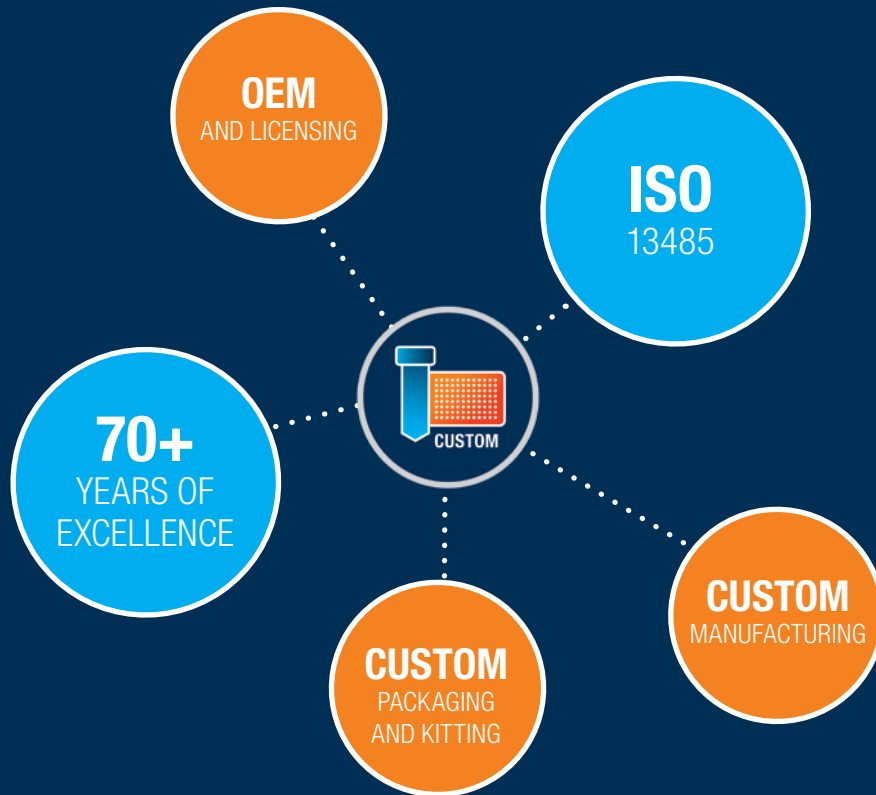


Custom PCR Plastics and Reagents

Partner with the PCR Experts



Committed to Quality



OEM, original equipment manufacturer; ISO, International Organization for Standardization.

A History of Excellence

Make Bio-Rad your trusted partner for tailored, reliable, and high-quality PCR solutions.

With a focus on excellence and customer service since 1952, Bio-Rad has become a global leader in developing and manufacturing innovative products for life science research and clinical diagnostics, helping to advance drug discovery and improve healthcare. We provide a full spectrum of products that complement the areas of cell biology, gene expression, protein purification, protein quantitation, drug discovery and manufacturing, food safety, and science education.

Our products are manufactured under international quality standards (ISO 13485). Stringent manufacturing quality controls guarantee the consistency, reproducibility, and quality you have come to expect from Bio-Rad.

As your custom manufacturing partner, we are committed to delivering exemplary products made to your unique specifications.

Our Commitment

Innovative ideas require unique solutions. We offer a suite of customizable products and services that ensure efficiency, reproducibility, and reliability at each step of discovery, development, and manufacturing. Our products are manufactured in ISO 13485–certified production facilities using the highest quality materials and technologies. An experienced global support team is available to partner with you to provide documentation and technical support.



Custom Amplification Reagents

Reach Your Goals Faster

Bio-Rad is a global leader in developing, manufacturing, and marketing a broad range of innovative products for life science research and clinical diagnostic markets. Finding a partner for your project that delivers the sensitive, reproducible, and robust detection you need in a PCR reagent is key to meeting your goals. Our PCR reagents and kits feature advanced proprietary enzyme technology and buffer formulations to deliver the performance your project requires. We offer customers access to our innovative products supplied in flexible fill sizes and with custom labels and packaging to meet individual requirements.

Our PCR reagents are manufactured to ISO 13485 standards and are quality controlled to ensure consistent production of products you can trust.

Custom Options for PCR Reagents

- **Fill size** — select from a variety of vials and containers
- **Packaging** — choose from various bags, pouches, and boxes
- **Kit configuration** — create a custom kit using existing components
- **Documentation** — request additional QC testing

Original Equipment Manufacturer (OEM) and Licensing

Our extensive expertise in molecular biology has led to the development of a broad product portfolio of reagents. Partner with us for OEM solutions or licensing of Bio-Rad technology to ensure success in your project. We welcome collaborations throughout all stages, from discovery to commercialization.

Trusted Performance

Advanced Buffer Formulation

Refined PCR reagent buffers deliver top-notch performance. Use them confidently with a broad range of complex sample types and target regions, as well as under varying reaction parameters (e.g., primer concentrations, annealing temperature, salt concentration, and cycling products).

Universal Reference Dyes

A proprietary blend of reference dyes in our supermixes means you won't need different system-specific dye concentrations. Instead, reaction setup is easy and data become less variable. Our reference dyes are compatible with high-ROX, low-ROX, and ROX-independent systems.



Robust Reliance Reverse Transcriptase

Reliance Reverse Transcriptase is genetically engineered for high thermostability and efficient, sensitive cDNA synthesis — even with challenging samples and targets.

Powerful Sso7d Fusion DNA Polymerase

Patented Sso7d Fusion DNA Polymerase* is designed to succeed where others fail. It features improved processivity for effective amplification through regions of strong secondary structures, GC-rich regions, and longer products — even in the presence of PCR inhibitors.

* U.S. patent 7,560,260.

PCR Reagent Selection

We offer a wide range of reagents for reverse transcription, PCR, and real-time PCR, which are optimized to generate accurate and reproducible data. Our PCR reagent supermixes feature a blend of passive reference dyes for universal real-time instrument compatibility. Customizable PCR reagents include the following:

<p>Real-Time PCR Supermixes</p>  <p>Visit bio-rad.com/Supermixes</p>	<p>SsoAdvanced Universal Supermixes</p> <ul style="list-style-type: none"> Highly tolerant to common PCR inhibitors Based on highly processive, proprietary Sso7d fusion technology for unparalleled performance 	<p>iTaq Universal Supermixes</p> <ul style="list-style-type: none"> Supplied as 2x supermixes Ideal for reliable and reproducible performance Available in both SYBR® and probe chemistries
<p>cDNA Synthesis Kits</p>  <p>Visit bio-rad.com/RTReagents</p>	<p>Reliance Select cDNA Synthesis Kit</p> <ul style="list-style-type: none"> Formulated with a novel reverse transcriptase engineered to perform well with challenging samples and targets 	<p>iScript cDNA Synthesis Kits</p> <ul style="list-style-type: none"> Versatile; family includes multiple kits designed for various applications Built around an RNase H+ MMLV reverse transcriptase plus an RNase inhibitor for reliable yields
<p>One-Step Reverse Transcription Quantitative PCR (RT-qPCR) Reagents and Kits</p>  <p>Visit bio-rad.com/OneStep</p>	<p>Reliance One-Step Multiplex RT-qPCR Supermix</p> <ul style="list-style-type: none"> Supplied as a 4x supermix Based on a proprietary reverse transcriptase Designed for highly sensitive and reliable detection of up to 5 targets 	<p>iTaq Universal One-Step Kits</p> <ul style="list-style-type: none"> Easy 2-tube system for simple reaction setup Compatible with SYBR® and probe-based RT-qPCR
<p>PCR Reagents</p>  <p>Visit bio-rad.com/StandardPCRReagents</p>	<p>iTaq DNA Polymerase</p> <ul style="list-style-type: none"> Antibody-mediated, hot-start Highly specific, sensitive, and easy to use 	<p>iProof High-Fidelity DNA Polymerase</p> <ul style="list-style-type: none"> Unique thermostable fusion polymerase Accurately amplifies long products from a variety of DNA templates
<p>Loading and Reference Dyes</p>  <p>Visit bio-rad.com/PCRDyes</p>	<p>Precision Blue Real-Time PCR Dye</p> <ul style="list-style-type: none"> Supplied as a concentrated, ready-to-use reagent Useful for increasing the visibility of real-time PCR reaction mixtures 	<p>ROX Passive Reference Dye</p> <ul style="list-style-type: none"> For use with PCR reagents without ROX reference dye for ROX-dependent PCR systems <p>Fluorescein Calibration Dye</p> <ul style="list-style-type: none"> Useful for normalization of well-to-well variations that may result from pipetting errors or instrument limitations

Our portfolio of patented and proprietary chemistries offer unique and unrivaled performance. Need help choosing the right PCR reagent? Visit bio-rad.com/ReagentsSelector to use our interactive selection tool to find the right products for your needs.



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

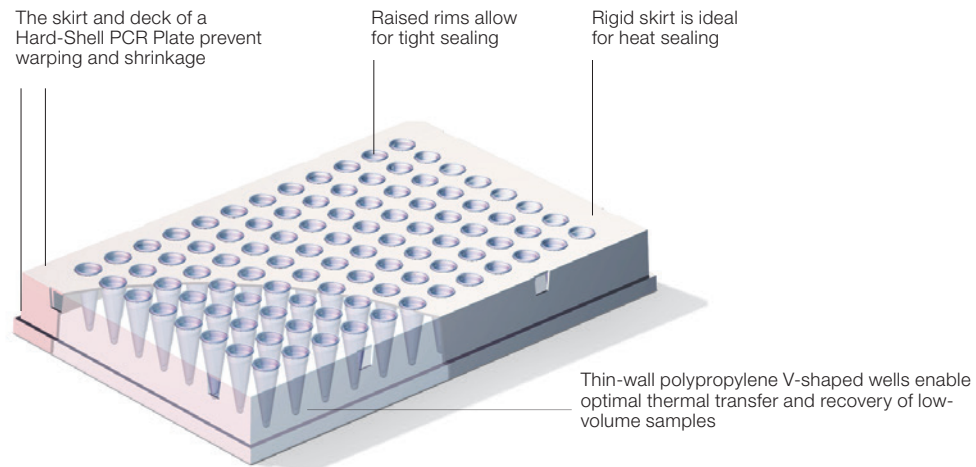
Custom PCR Plastics: Hard-Shell PCR Plates

Precisely Manufactured to Meet Your Needs

We are committed to manufacturing high quality Hard-Shell PCR Plates you can trust. We customize Hard-Shell PCR Plates to your specifications.

Hard-Shell PCR Plates

Customizing our Hard-Shell PCR Plates means taking advantage of our two-component design, delivering superior uniformity and precise handling in any situation. In addition to our broad selection of Hard-Shell PCR Plates, we offer a range of custom options to best fit your specific needs.



Features of Hard-Shell PCR Plates.

Hard-Shell PCR Plates offer:

- Automation compatibility
- Superior well-to-well uniformity for reliable and reproducible real-time PCR results
- Warp resistance and superior stability for durability during robotic handling, high-speed centrifugation, and storage
- Optional user-readable barcode with alphanumeric labels for easy well identification in high-throughput settings
- Footprint and well spacing designed to match American National Standards Institute (ANSI) and Society for Biomolecular Screening (SBS) standard dimensions
- Plastics certified free from DNase, RNase, and human DNA
- Choice of 96- or 384-well format

Customization Process for Plates and Reagents

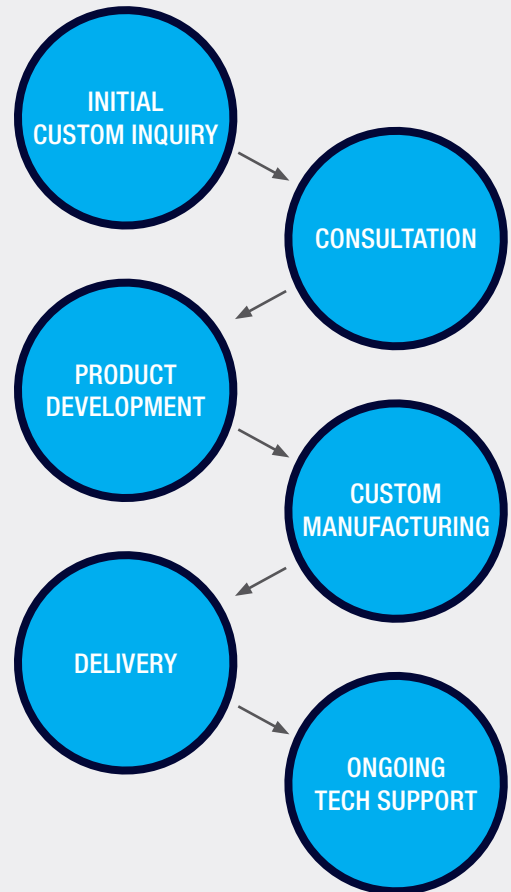
Custom Options for Hard-Shell PCR Plates

- **Barcoding** — specify placement and sequence of 1-D or 2-D barcodes
- **Packaging** — select from various box sizes and packaging formats
- **Color** — choose from a range of shell and well color combinations
- **Branding** — include your company name and logo on PCR plates and packaging
- **Labeling** — customize to your product specifications

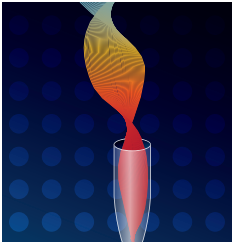
Description	Clear Well	White Well	Black Well
Hard-Shell Low-Profile 96-Well Skirted PCR Plates, 50 plates			
White shell	HSP9601	HSP9655	—
Red shell	HSP9611	—	—
Yellow shell	HSP9621	—	—
Blue shell	HSP9631	HSP9635	—
Green shell	HSP9641	HSP9645	—
Black shell	HSP9661	HSP9665	HSP9666
White shell, barcoded (Row H and Column 12)	HSP9901	HSP9955	—
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	—	—
Black shell	—	HSP3865	HSP3866
Clear shell, barcoded (Row P and Column 24)	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	—	—

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

No custom manufacturing process is the same. Our knowledgeable staff will be your guide through each step of the process. From the first inquiry and consultation, through the first article inspection and production scale-up, and continuing throughout the partnership, our team of experts is dedicated to ensuring your custom product meets your specific needs on your timeline.



Visit bio-rad.com/CustomPCR to learn more about our custom program.



Our Promise

To take pride in being your global partner and use a customer-first approach at every step of the journey.

Visit bio-rad.com/CustomPCR for more information.

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. 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.

BIO-RAD

**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** 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 2 3473 4460 **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

