

# **Custom PCR Reagents**

Innovative Ideas Require Unique Solutions

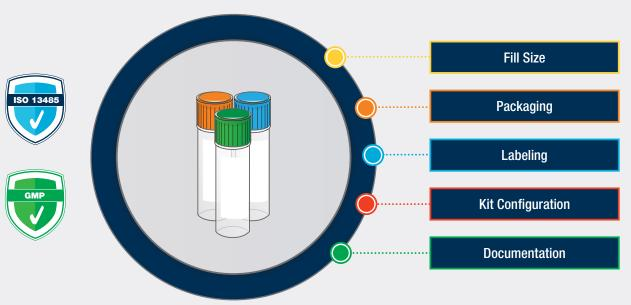
## **Imagine Our Advanced Enzyme Technology Tailored to Your Needs**

#### **High-Performance Reagents**

Do you need a commercial supply partner for high-quality, robust, and reliable PCR, quantitative PCR (qPCR), and reverse transcription reagents? From concept to commercialization, Bio-Rad Laboratories, Inc. offers a range of custom packaging and kitting options for our existing product components, including our advanced enzyme technology and buffer formulations.

With more than 70 years of experience developing and manufacturing products for life sciences and clinical diagnostics customers, Bio-Rad provides the expertise and technology to support your original equipment manufacturer (OEM) amplification needs. Bio-Rad custom PCR reagents are manufactured under good manufacturing process (GMP) and International Quality Standards (ISO 13485:2016) and are quality controlled to guarantee consistent products you can trust.

### **Custom Options**





#### **Custom Options for PCR Reagents**

**Fill Size** — Choose smaller volumes that can be part of a kit or larger volumes for dispensing at your facility.

**Packaging** — Select from various bags, pouches, and boxes to create a finished look.

**Labeling** — Customize labels with your product information and logo to promote your trusted brand.

**Kit Configuration** — Create custom kits using our existing components for efficiency and convenience.

**Documentation** — Request additional quality control testing for your assay or application.

#### **Meet the Bio-Rad PCR Reagent Family**

As your custom manufacturing partner, we are committed to delivering reproducible, reliable, and high-quality products on demand and in your specified format. Choose from our existing portfolio of products.

#### Real-Time PCR Supermixes and Kit

For superior qPCR performance in any condition, with any detection chemistry, and on any instrument:

- SsoAdvanced Universal Supermixes
- iTaq Universal Supermixes

#### Complementary DNA (cDNA) Synthesis Supermixes and Kits

For easy setup, high reproducibility, and bias-free cDNA synthesis:

- Reliance Select cDNA Synthesis Kit
- iScript cDNA Synthesis Kit

## One-Step Reverse Transcription qPCR (RT-qPCR) Reagents and Kits

For robust and reliable amplification of RNA targets on all qPCR platforms, including both carboxy-X-rhodamine (ROX)-dependent and ROX-independent systems:

- Reliance One-Step Multiplex Supermix
- iTaq Universal One-Step RT-qPCR Kit

#### High-Fidelity and Standard PCR Reagents and Kits

For convenience and improved yields in standard and high-fidelity reverse transcription PCR:

- iProof High-Fidelity DNA Polymerase
- iTaq DNA Polymerase

#### PCR Loading and Reference Dyes

For setup and normalization of real-time PCR experiments



Did you know? You can also customize our Hard-Shell<sup>TM</sup> Plates with branding, barcodes, packaging, and configurations to meet your specific needs.

Visit bio-rad.com/CustomPlastics to learn more.

View our Amplification Reagents and Plastics brochure to learn more about our PCR Reagent family.

For research use only. Not for use in diagnostic procedures.

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. Copyright © 2023 by Bio-Rad Laboratories, Inc. All rights reserved.

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 Gramany 00 800 00 24 67 23 Gramany 00 800 00 24 67 23 Gramany 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

Bulletin 3550 Ver A US/EG 23-0570 0923 Sig 0123

