



## Real-Time PCR Solutions

Advanced technologies for data you need, every time



# Trusted Performance

## Advanced Buffer Formulation

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

## Universal Reference Dyes

A proprietary blend of reference dyes 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.

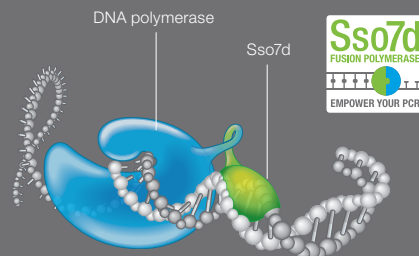


## Robust Reliance Reverse Transcriptase

The 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

Our 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, longer products — even in the presence of PCR inhibitors.





Bio-Rad's PCR Reagents Are:

Accurate  
Efficient Quality  
Reproducible  
Consistent Reliable  
Easy

# qPCR Reagents for All, Data You Can Rely On

Bio-Rad's universal qPCR reagents feature:

- Single tube, 2x ready-to-use supermix
- Advanced buffer formulation for optimal performance under varying conditions
- A proprietary blend of universal reference dyes\* compatible with all qPCR systems
- Flexible cycling compatibility for both fast and standard protocols

## SsoAdvanced Universal Supermixes

Expect unparalleled performance based on the highly processive Sso7d Fusion DNA Polymerase\*\* with maximum sensitivity, efficiency, and reproducibility.

Our SsoAdvanced supermixes are highly tolerant to common PCR inhibitors, making them ideal for compromised samples.



You perform...	Your sample is...	Your target is/has...
<ul style="list-style-type: none"> <li>▪ Gene expression</li> <li>▪ Single-copy detection</li> <li>▪ lncRNA and miRNA expression</li> <li>▪ Low-copy viral/pathogen detection</li> <li>▪ Methylation studies (probe)</li> <li>▪ Copy number variation</li> <li>▪ Mutation detection</li> <li>▪ ChiP</li> <li>▪ Genotyping</li> <li>▪ siRNA knockdown</li> <li>▪ Multiplexing of up to 2 targets</li> </ul>	<ul style="list-style-type: none"> <li>▪ gDNA from various samples, including tissue, cell lines, plant, soil, FFPE, and LCM</li> <li>▪ Cell lysate</li> <li>▪ Crude extraction</li> <li>▪ Blood spot</li> <li>▪ Plant extract</li> <li>▪ cDNA</li> <li>▪ Plasmid</li> </ul>	<ul style="list-style-type: none"> <li>▪ Highly GC rich (&gt;60%)</li> <li>▪ Highly AT rich (&gt;60%)</li> <li>▪ Increased secondary structure</li> <li>▪ &gt;250 bp</li> </ul>

## iTaq Universal Supermixes

These supermixes are ideal for reliable and reproducible performance for various real-time PCR applications.



You perform...	Your sample is...	Your target is/has...
<ul style="list-style-type: none"> <li>▪ Gene expression</li> <li>▪ Genotyping</li> <li>▪ Mutation detection</li> <li>▪ Pathogen detection</li> <li>▪ Multiplexing of up to 2 targets</li> </ul>	<ul style="list-style-type: none"> <li>▪ gDNA mostly from tissue and cell lines</li> <li>▪ cDNA</li> <li>▪ Plasmid</li> </ul>	<ul style="list-style-type: none"> <li>▪ 40–60% GC-rich</li> <li>▪ 50–250 bp</li> </ul>

\* U.S. patent 9,493,824. \*\* U.S. patents 6,627,424; 7,541,170; and 7,560,260.

## Flexible Solutions for Sensitive qPCR Directly from RNA

Bio-Rad's One-Step RT-qPCR Supermix and Kits offer:

- Convenient reaction setup using 1- or 2-tube system
- A proprietary blend of universal reference dyes\* to be compatible with all qPCR systems
- Advanced buffer formulation for optimal performance under varying conditions

### Reliance One-Step Multiplex Supermix

This 4x supermix is expertly formulated for highly sensitive and reliable detection of up to 5 targets directly from RNA. It features genetically engineered ultra-sensitive Reliance Reverse Transcriptase and an optimized blend of powerful DNA polymerases, including Sso7d Fusion DNA Polymerase\*\*, which provide advanced inhibitor tolerance. Assembled reaction stability (24-hour) makes it ideal for automation and high-throughput workflows.



You perform...	Your sample is...	Your target is/has...
<ul style="list-style-type: none"> <li>▪ Multiplexing of up to 5 targets</li> <li>▪ Gene expression</li> <li>▪ Low-copy viral/pathogen detection</li> <li>▪ Food safety studies</li> <li>▪ Diagnostic research</li> <li>▪ High-throughput screening</li> <li>▪ Automated testing</li> </ul>	<ul style="list-style-type: none"> <li>▪ Difficult matrices such as blood, serum, or plasma</li> <li>▪ RNA from various samples, including tissue and cell lines</li> <li>▪ Cell lysate</li> <li>▪ Potentially contaminated with PCR inhibitors</li> </ul>	<ul style="list-style-type: none"> <li>▪ Highly GC rich (&gt;60%)</li> <li>▪ Internal positive control of RNA or DNA</li> <li>▪ 70–250 bp</li> </ul>

### iTaq One-Step Universal Kits

Achieve reliable target detection directly from RNA by setting up a reaction with our one-step kits enabled by a 2x qPCR supermix and reverse transcriptase.



You perform...	Your sample is...	Your target is/has...
<ul style="list-style-type: none"> <li>▪ Gene expression</li> <li>▪ lncRNA and miRNA expression</li> <li>▪ Viral/pathogen detection</li> <li>▪ Mutation detection</li> <li>▪ Pathogen detection</li> <li>▪ siRNA knockdown</li> <li>▪ Multiplexing of up to 2 targets</li> </ul>	<ul style="list-style-type: none"> <li>▪ RNA, mostly from tissue and cell lines</li> <li>▪ Cell lysate</li> </ul>	<ul style="list-style-type: none"> <li>▪ Highly GC rich (&gt;60%)</li> <li>▪ Highly AT rich (&gt;60%)</li> <li>▪ 40–60% GC-rich (SYBR®)</li> <li>▪ 50–250 bp</li> </ul>

cDNA, complementary DNA; Cq, quantification cycle; ChIP, chromatin immunoprecipitation; FFPE, formalin-fixed paraffin embedded; LCM, laser capture microdissection



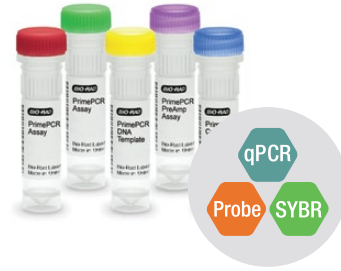
Integrate Bio-Rad's qPCR reagents with the CFX Automation System II to enable walk-away high-throughput qPCR operation.



Visit [bio-rad.com/CFXauto](http://bio-rad.com/CFXauto) for more information.



Simplify your gene expression or lncRNA expression with expertly designed primers and probe assays and predesigned panels.



Visit [bio-rad.com/PrimePCR](http://bio-rad.com/PrimePCR) for more information.



Use high-quality PCR plastics to be confident in your qPCR data. Choose from an extensive range of tube strips, and 48-well, 96-well, and 384-well PCR plates.



Visit [bio-rad.com/PCRPlastics](http://bio-rad.com/PCRPlastics) for more information.



Choose the best cDNA synthesis kit for your application, with formulations optimized for ease, yield, flexibility, fidelity, or gDNA clearance.



Visit [bio-rad.com/iScript](http://bio-rad.com/iScript) for more information.

## Ordering Information

Detection Method	Reaction Size	SsoAdvanced Universal	iTaq Universal	Reliance One-Step Supermix	iTaq Universal One-Step	SsoAdvanced Universal Inhibitor-Tolerant
SYBR® Green	100 x 20 µl reactions	–	–	–	1725150	–
	200 x 20 µl reactions	1725270	1725120	–	–	1725016
	500 x 20 µl reactions	1725271	1725121	–	1725151	1725017
	1,000 x 20 µl reactions	1725272	1725122	–	–	1725018
	2,500 x 20 µl reactions	1725274	1725124	–	–	–
	5,000 x 20 µl reactions	1725275	1725125	–	–	–
Probe	100 x 20 µl reactions	–	–	–	1725140	–
	200 x 20 µl reactions	1725280	1725130	12010176	–	–
	500 x 20 µl reactions	1725281	1725131	–	1725141	–
	1,000 x 20 µl reactions	1725282	1725132	12010220	–	–
	2,000 x 20 µl reactions	–	–	12020221	–	–
	2,500 x 20 µl reactions	1725284	1725134	–	–	–
5,000 x 20 µl reactions	1725285	1725135	–	–	–	

Visit [bio-rad.com/PCRReagents](http://bio-rad.com/PCRReagents) for more information.



Bio-Rad's PCR reagents are manufactured under an ISO 13485:2016 certified Quality Management System and are quality control tested to ensure consistent product performance you can trust.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

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.

All trademarks mentioned herein are the property of their respective owner.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600  
**Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30  
**Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580  
**Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300  
**United Kingdom** 44 01923 47 1301

