



EZ-Check™ Real-Time PCR Solution

Simple Workflow, Reliable Results from the Brand You Trust

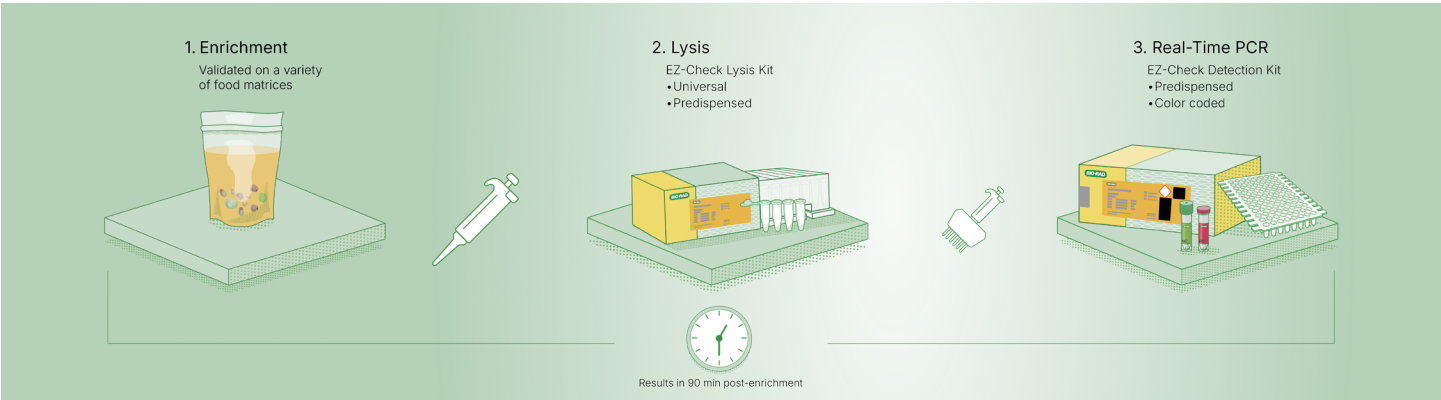
The EZ-Check Real-Time PCR Solution from Bio-Rad is the next generation in rapid and accurate detection of food pathogens. Our EZ-Check Solution allows the qualitative detection of target bacteria in a broad range of food products, animal feed, and environmental samples.

- **Simple workflow:** only 2 pipetting steps
- **Fast time to results:** ~90 minutes from enrichment to results
- **Built-in control:** an internal amplification control in every well to monitor false negatives
- **Validated instruments:** compatible with multiple Bio-Rad thermal cyclers for maximum flexibility
- **Automation:** fully integrated with the iQ-Check™ Prep System and Free DNA Removal Solution



EZ-Check Streamlined Workflow

Only Two Pipetting Steps Needed



Choose the EZ-Check Solution for Efficiency and Confidence

The EZ-Check Solution has a streamlined design that reduces labor, handling time, and training, boosting operational efficiency and lowering user error. The solution meets both AOAC and ISO standards, ensuring regulatory compliance, and the workflow is scalable for labs of all sizes. What's more, the EZ-Check Solution seamlessly integrates with the current Bio-Rad portfolio of food pathogen detection methods and uses nontoxic reagents with minimal waste, generating a safe and sustainable lab environment.

Ordering Information

Catalog #	Description	Catalog #	Description
Supplies		Accessories	
12018075	EZ-Check Lysis Kit	12019781	EZ-Check Lysis Rack
12018082	EZ-Check <i>Salmonella</i> spp.	12019771	EZ-Check PCR Frame
17010251	EZ-Check <i>Salmonella</i> spp. with Lysis Kit, includes 12018082 and 12018075	12022466	MHR13 EZ-Check Block
12018844	EZ-Check <i>Listeria</i> spp. Kit	12022580	EZ-Check Lysis Storage Caps, 600
17010252	EZ-Check <i>Listeria</i> spp. with Lysis Kit, includes 12018844 and 12018075	Equipment and Software	
12018858	EZ-Check <i>Listeria monocytogenes</i> Kit	17010702	Micropipet 8-Channel, 20–200 µL
17010253	EZ-Check <i>L. monocytogenes</i> with Lysis Kit, includes 12018858 and 12018075	17010254	MHR13 with EZ-Check Block
3594970	Free DNA Removal Solution	17007992	CFX Opus 96 Real-Time PCR System
17010688	Xcluda Aerosol Barrier Pipet Tips (Style D), sterile, 20–200 µL	17007991	CFX Opus Deepwell Real-Time PCR System
		17010275	CFX Duet Real-Time PCR System for Food Applications
		3593893	CFX Maestro Software, Industrial Diagnostic Edition

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

Bio-Rad and EZ-Check are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. IQ-Check is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner. ©2026 Bio-Rad Laboratories, Inc.



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 55 11 3065 7550 Canada 1 800 361 1808 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 Greece 30 210 7774396 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 000 800 00 24 67 23 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 55 5488 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 7 495 721 14 04 Singapore 65 6415-3170 South Africa 27 21 531 7504 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 662 651 8311 United Arab Emirates 971 4 818 7300 United Kingdom 00 800 00 24 67 23									