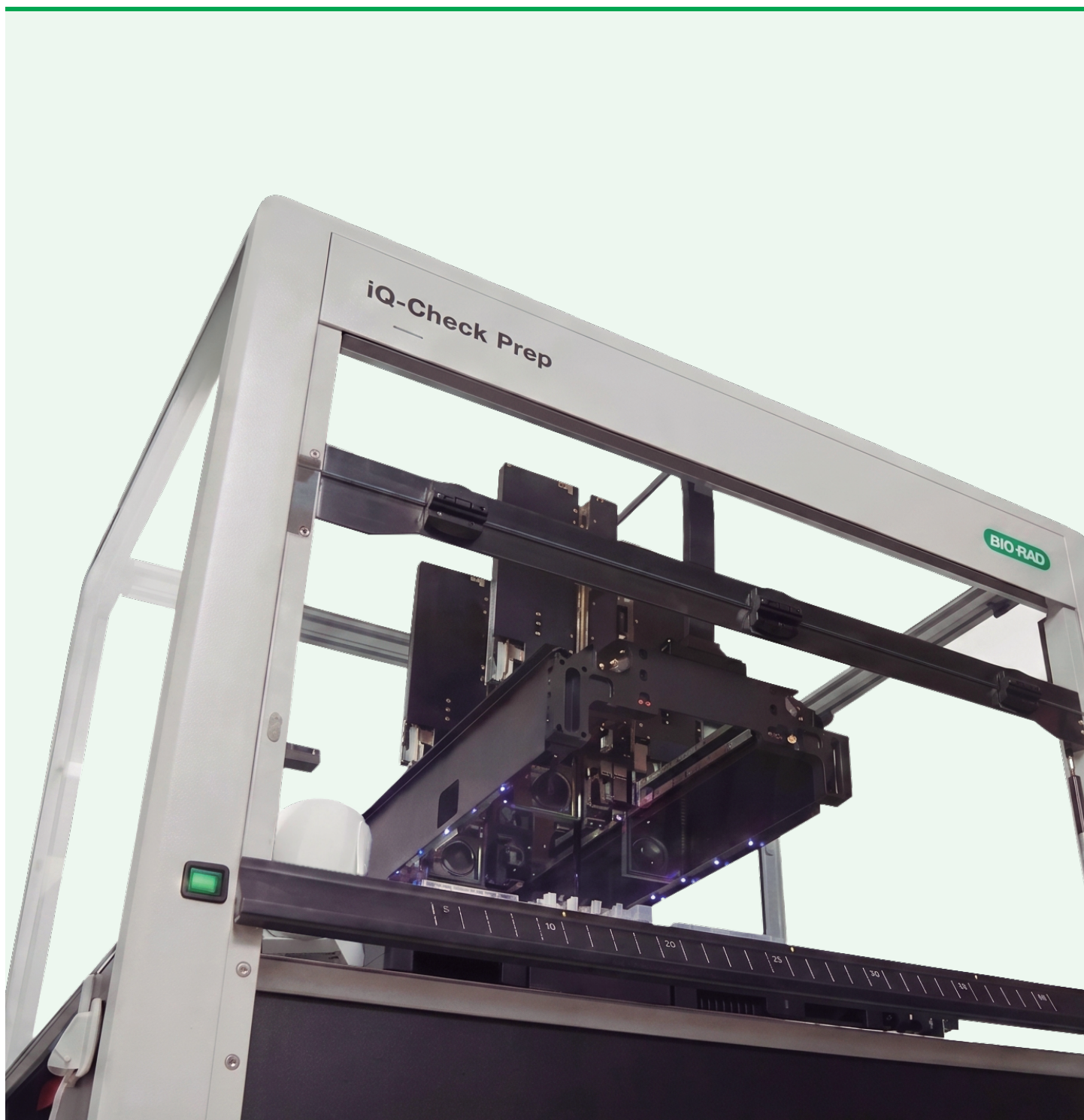


Automating the EZ-Check™ Method

Designed for Simplicity. Built for Reproducibility.

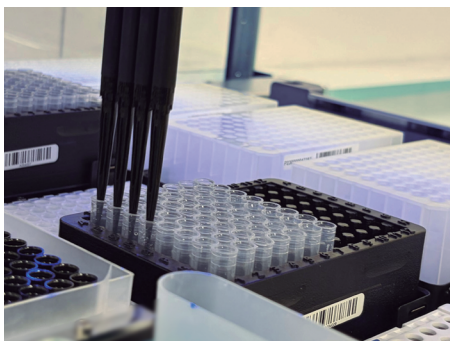


Visit [bio-rad.com/solutions](https://www.bio-rad.com/solutions) for more information.

An Integrated Automation Solution

Designed for Real-World Lab Workflows

When paired with EZ-Check kits, the iQ-Check™ Prep System reinforces consistency, traceability, and reproducibility without adding complexity. Automation standardizes sample preparation and PCR setup, reduces the already few hands-on steps, and supports reliable results across users. These attributes make EZ-Check detection automation an ideal fit for laboratories seeking an easy-to-adopt, scalable solution that balances simplicity with operational control.



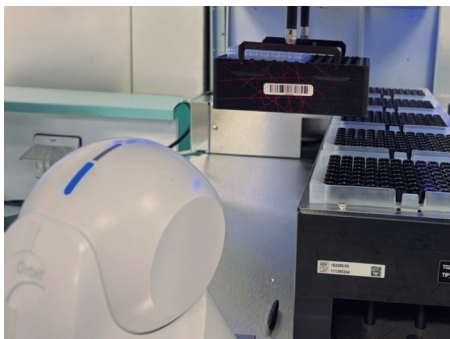
Walk-Away Automation

- Preserve simplicity while improving traceability and reproducibility
- Combine manual and automated workflows to match daily lab needs and sample volumes
- Receive result notifications by email and reallocate laboratory resources to higher-value tasks
- Maintain consistent testing workflows during staffing gaps, weekends, or training periods



Increased Productivity, Flexibility, and Scale

- Prepare two plates at one time with simultaneous processing
- Run multiple targets in parallel within a standardized workflow
- Scale testing from small daily runs to higher volumes without increasing workflow complexity or staffing
- Automate sample dilution preparation and Free DNA Removal Solution protocols

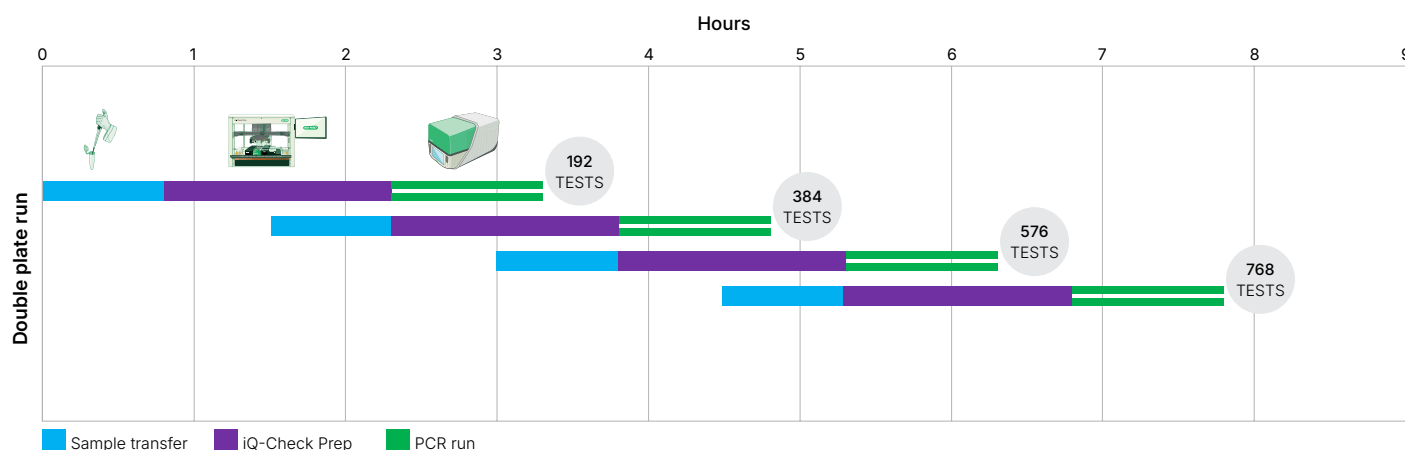
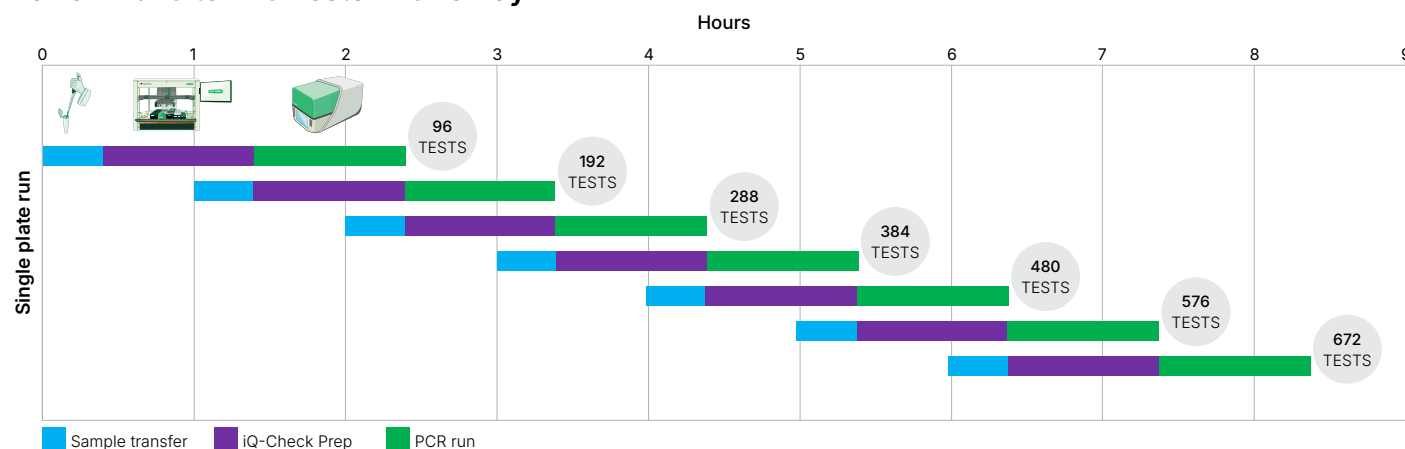


Greater Confidence and Efficiency

- Strengthen traceability with barcodes and laboratory information management system (LIMS) integration
- Rely on internationally validated methods
- Improve reproducibility by reducing user-to-user variability
- Minimize cross-contamination risk through simplified, streamlined automation



Perform 670 to 770 Tests in One Day

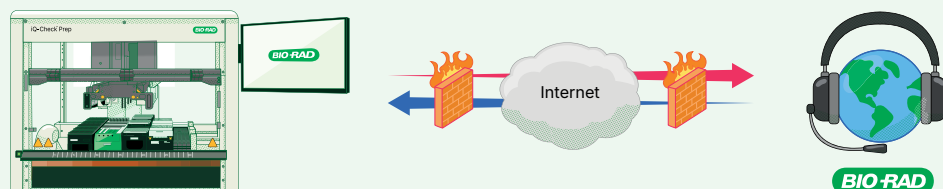


Remote Support via BRiCare

BRiCare is a remote support and monitoring system that gives Bio-Rad the ability to rapidly troubleshoot and solve your system problems.

Benefits

- Minimal interruptions
- Fast guidance and training
- Enhanced lab productivity
- Greater peace of mind



General Specifications

Feature	Specifications
Instrument dimensions (W x D x H)	1,124 x 795 x 903 mm (44.25 x 31.30 x 35.55 in.)
Work area dimensions (W x D x H)	675 x 465 x 195 mm (26.57 x 18.31 x 7.68 in.)
Weight	120 kg (265 lb)
Positional accuracy	X-Y-Z positional accuracy of 0.1 mm
Tip size	Single volume 300 µL
Liquid level detection	Capacitive liquid level detection (cLLD) and pressure (pLLD) on aspiration, cLLD on dispense
Labware	Recommended labware from Bio-Rad

Operating Data

Feature	Specifications
Maximum power consumption	600–1,000 VA (depending on configuration)
Voltage	115/230 VAC
Frequency	50/60 Hz ± 5%
Delayed action fuse	600 VA 115 V: 6.3 A (T6.3AL250) 230 V: 3.15 A (T3.15AL250) 1,000 VA 115 V: 10 A (T10AL250)
Installation category	II
Pollution degree	2
Temperature range	15–35°C
Relative humidity	30–85% (noncondensing, indoors)
Noise level	<65 dBA (per EN27779)
Elevation maximum	2,000 m above sea level

Ordering Information

Catalog #	Description
Instruments	
3594911	iQ-Check Prep System
17007991	CFX Opus Deepwell Real-Time PCR System
17007992	CFX Opus 96 Real-Time PCR System
17010275	CFX Duet Real-Time PCR System for Food Applications

Software	
3593893	CFX Maestro Software, Industrial Diagnostic Edition

Consumables	
3594900	iQ-Check Deep Well Microplates, 96-well plates, 50
12022468	iQCP Tips, 300 µL, case of 5,760
12014476	iQCP Waste Bag, case of 25
12014473	iQCP Reservoir, case of 28
12022581	iQCPv5 Labware Barcodes, Pkg of 1,000

Kits and Reagents	
12018075	EZ-Check Lysis Kit
12018844	EZ-Check <i>Listeria</i> spp. Kit
12018858	EZ-Check <i>Listeria monocytogenes</i> Kit
12018082	EZ-Check <i>Salmonella</i> spp. Kit
3594979	Free DNA Removal Solution, 1 viral

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

Bio-Rad, EZ-Check, and Hard-Shell 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