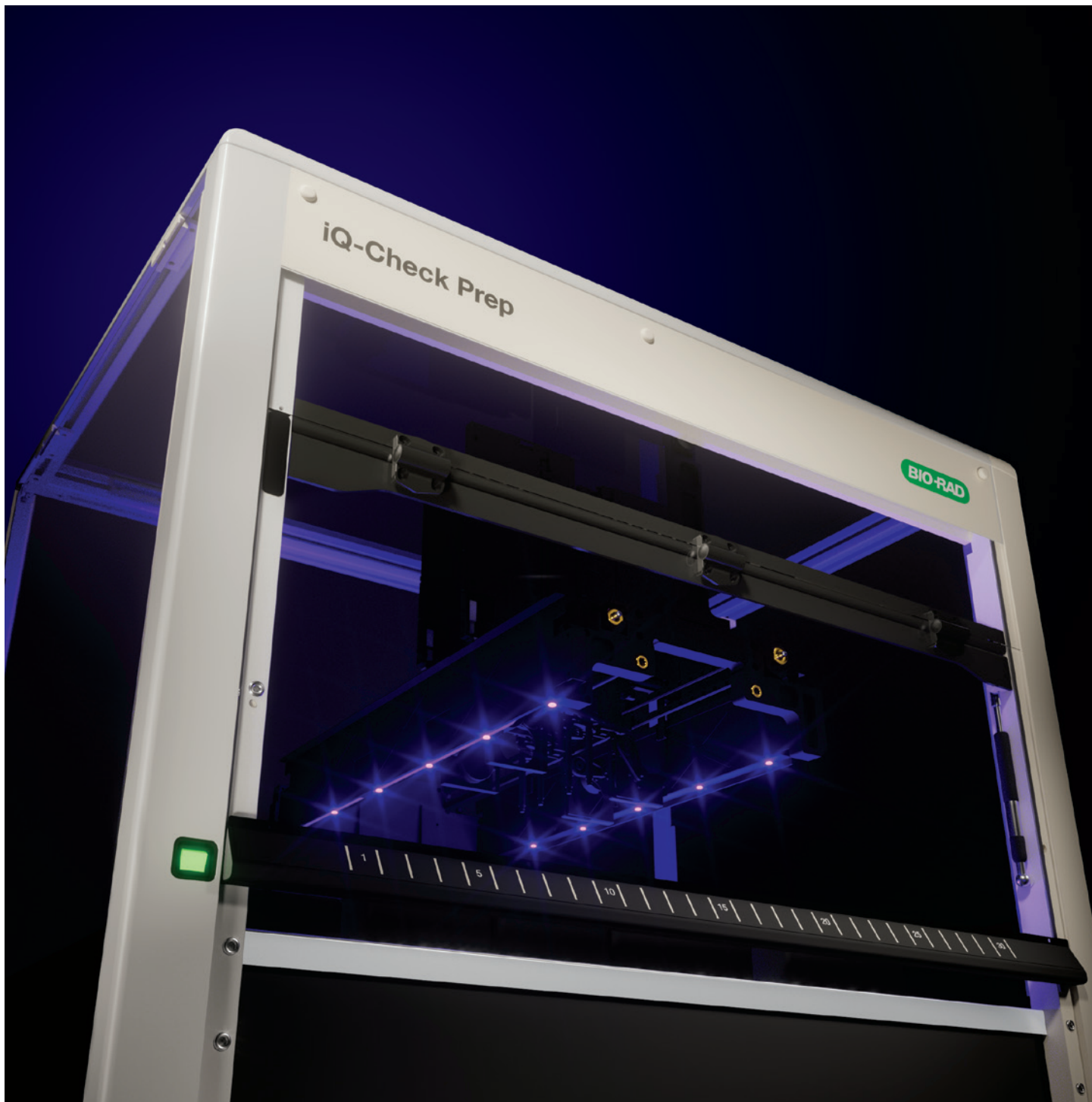


iQ-Check™ Prep Solution

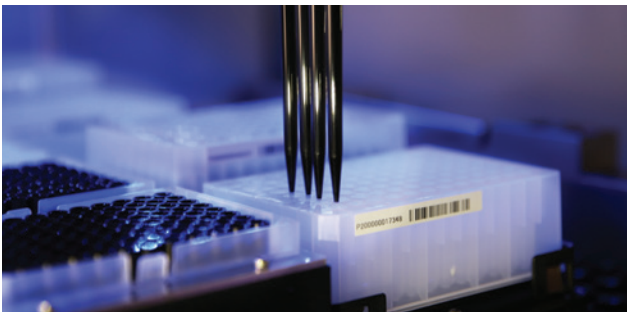
Maximize Lab Performance with the Power of Automation



An Integrated Automation Solution

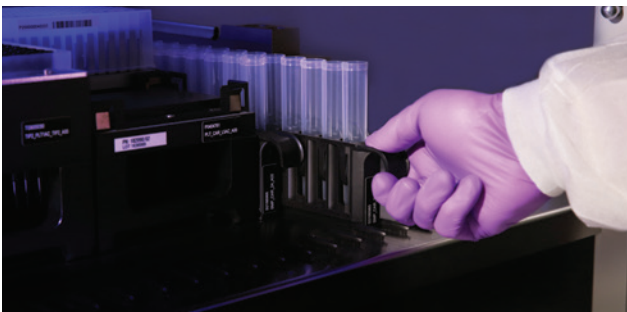
Designed for the Way You Work

The iQ-Check Prep Solution is everything you need to simplify your testing regimen — systems, kits, consumables, software, service, and support. Designed for use with our full range of iQ-Check Food Pathogen Detection Kits together with the CFX96 Touch Deep Well and CFX Opus Real-Time PCR Systems, the iQ-Check Prep Solution streamlines food pathogen and cannabis testing workflows, improves laboratory staff utilization, and significantly reduces technical errors that cause delays and damage credibility. Robustness and flexibility make the iQ-Check Prep Solution an ideal fit for medium- and high-throughput labs.



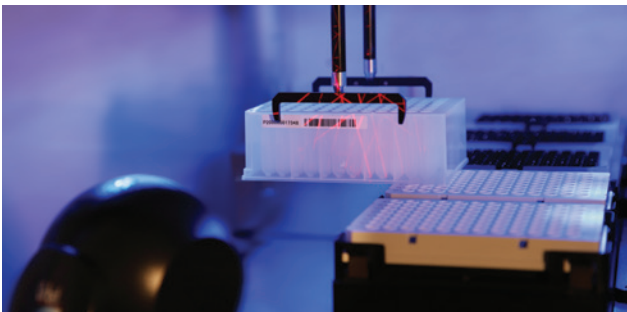
Walk-Away Automation

- Fully automate DNA extraction, PCR setup, detection, and analysis with the most integrated PCR automation solution on the market
- Define automated and manual steps with flexible plate setup
- Receive result notifications by email and reallocate lab resources to perform other tasks
- Maintain throughput during down periods (weekends, vacation)



Increased Productivity, Flexibility, and Scale

- Process samples in individual tubes or deep well plates
- Run up to 4 different iQ-Check Assays simultaneously
- Test a few or as many as 500 tests in an 8 hr shift
- Automate the iQ-Check Free DNA Removal Solution protocol



Greater Confidence and Efficiency

- Improve traceability with barcodes and laboratory information management system (LIMS) integration
- Trust each pipet step with liquid level sensing in real time
- Maximize reproducibility by removing human error
- Avoid cross-contamination



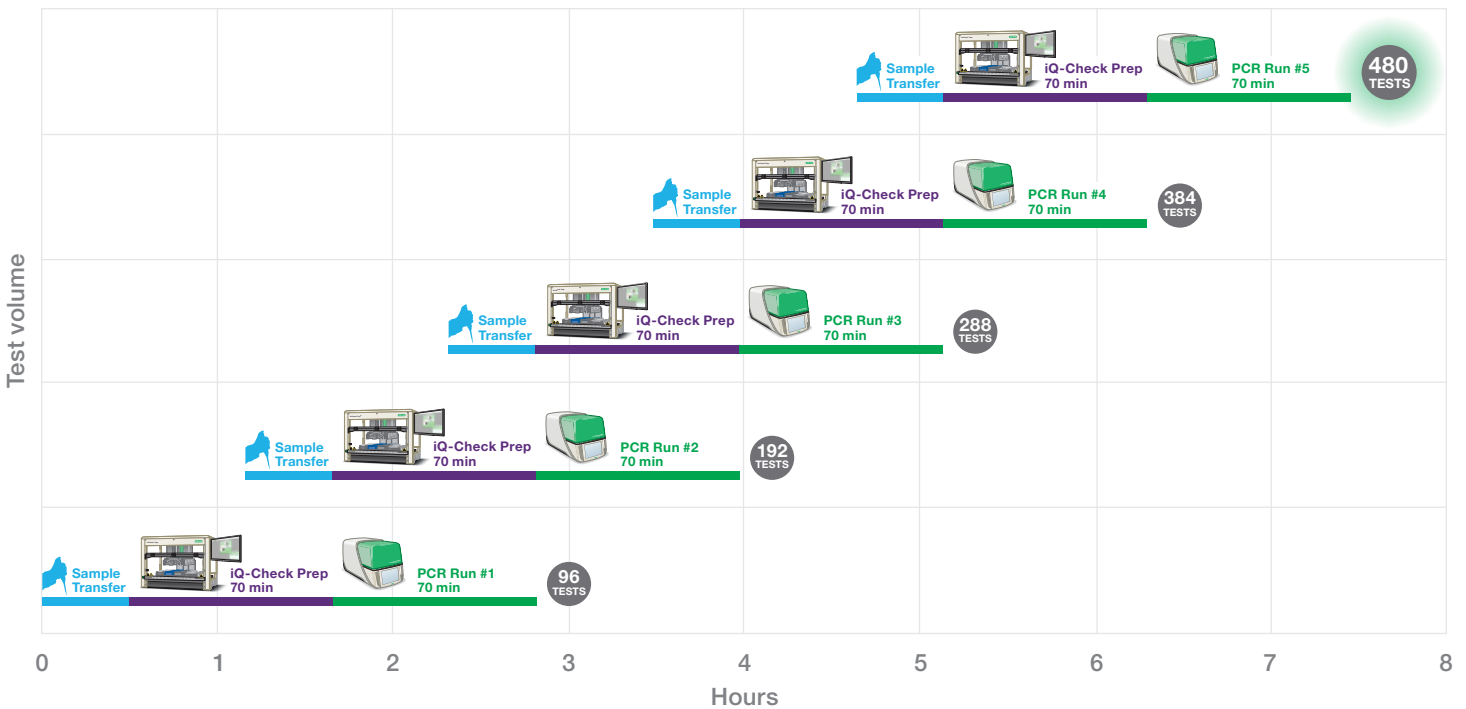
Designed to Make Your Work Easier

- Rely on internationally validated methods
- Reduce hands-on time with ready-to-use, load-and-go reagents
- Free up time with limited maintenance requirements
- Access proactive monitoring and remote support via BRiCare (Bio-Rad™ Instrument Care)

Scan the QR code to view our validations and certifications.



Perform Nearly 500 Tests in One 8 Hour Shift



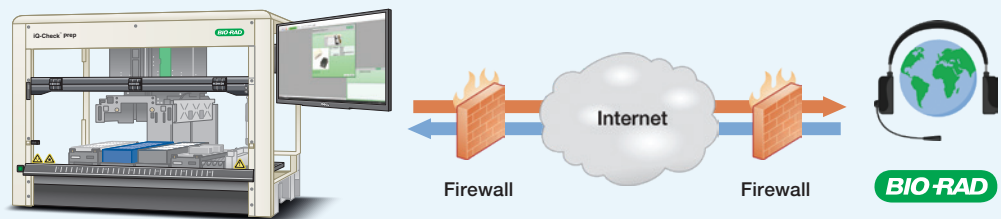
By overlapping sample preparation and PCR steps, you can perform 480 tests in a single 8 hr shift using one iQ-Check Prep System and one real-time PCR system. Several real-time PCR systems can be connected to one computer if required. You can reallocate laboratory resources to perform other tasks during system run time.

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	Specification
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	Low volume: 50 µl; high volume: 1,000 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	Specification
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)
Altitude maximum	2,000 m above sea level

Ordering Information

Catalog # Description

Instruments

3594911 iQ-Check Prep System
3600037 CFX96 Touch Deep Well Real-Time PCR System
17007991 CFX Opus Deepwell Real-Time PCR System
17007992 CFX Opus 96 Real-Time PCR System

Software

3593893 CFX Manager Software, Industrial Diagnostic Edition

Consumables

3594900 iQ-Check Deep Well Microplates, 96-well plates, 50
12016673 Tube for PCR Mix, 5 ml, 25
12014486 iQCP Tips, 50 µl, case of 5,760
12014483 iQCP Tips, 1,000 µl, case of 3,840
12014476 iQCP Waste Bag, case of 25
12014473 iQCP Reservoir, case of 28
HSP9955XTU Hard-Shell™ 96-Well PCR Plates, barcoded

Kits and Reagents

3578135 iQ-Check *Campylobacter* PCR Detection Kit, 96 tests
3578137 iQ-Check *Cronobacter* PCR Detection Kit, 96 tests
3578114 iQ-Check *E. coli* O157:H7 PCR Detection Kit, 96 tests
12003068 iQ-Check *Enterobacteriaceae* PCR Detection Kit, 96 tests
3578124 iQ-Check *Listeria monocytogenes* II PCR Detection Kit, 96 tests
3578113 iQ-Check *Listeria* spp. PCR Detection Kit, 96 tests
3578123 iQ-Check *Salmonella* II PCR Detection Kit, 96 tests
3578142 iQ-Check *S. Enteritidis* PCR Detection Kit, 96 tests
3578139 iQ-Check STEC VirX PCR Detection Kit, 96 tests
12013174 iQ-Check STEC SerO II PCR Detection Kit, 32 tests
12004306 iQ-Check *S. Typhimurium* PCR Detection Kit, 96 tests
12006574 iQ-Check *Vibrio* PCR Detection Kit, 96 tests
3594970 iQ-Check Free DNA Removal Solution, 1 vial

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

BIO-RAD 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.
© 2022 Bio-Rad Laboratories, Inc.

“In a high-throughput environment like a cannabis testing laboratory, having reliable and consistent automation is invaluable. Partnering with Bio-Rad and utilizing the iQ-Check Prep Automated platform has allowed us to virtually eliminate operator error from our qPCR preparation. With the iQ-Check Prep, we are able to train even the newest laboratory technicians to execute qPCR at an expert level over the course of a few days.”

MICHAEL ESPOSITO
MCR LABS

“Speed, simplicity, quality, better work organization, and waste reduction. The Bio-Rad automation system allowed us to greatly cut down the analysis time for food pathogens while maintaining excellent quality of results.”

REMIGIO BERTO
EPTA NORD SRL

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.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

