



...brought to light

The PhD™ Ix System builds on the tradition of the original PhD™ System with a focus on bringing simple, yet powerful, solutions to light. Empower your laboratory with streamlined workflows, versatile performance, and robust security. Enjoy the efficiency and quality that comes from trusted automation.



Streamlined
Save "Hands-On" Time



Versatile
Assay Programming



Secure
Traceable IFA/EIA Processing

AUTOIMMUNE TESTING

PhD™ Ix IFA/EIA System

Specifications

Dimensions (WxDxH) 36x27x30 in. (91x68x76 cm)
 Power Input Requirements 100 - 240 VAC at 50/60 Hz
 Power Consumption 220 W max

Workstation Capacity

Sample Capacity EIA: 192 or IFA: 96
 Microplate/Slide Capacity 2 microplates/20 slides
 Number of Different Assays on Board EIA: 8 or IFA: 4
 EIA Walkaway Time 1 full microplate with 3x30 incubations: 2 hr 15 min
 IFA Walkaway Time 96 slide wells with 2x30 incubations: 1 hr 50 min

Workstation Fluidics

Volumetric Pump Proprietary Valveless Ceramic design
 Sample Volume Delivery 1-100µL
 Maximum Reagent Delivery 1000µL
 EIA Incubation Temperature Range up to 40°C
 Precision CV <6% for all volumes
 EIA Plate Washing 3 modes: normal, settable residual wash, well-by-well wash/reagent dispense
 IFA Slide Washing 2 modes: cycle wash & hyperwash with simultaneous dispense/aspiration

Peripheral Specifications

Network Multiple Workstations Yes, with a single computer, microplate reader, and LIS connection
 Bar Code Reading All typical symbologies, including Codabar, Code 39, Code 128, and Interleaved 2 of 5
 Data Reduction (Interpolation) 4PL, 5PL, point-to-point, linear regression
 ASTM LIS Connectivity Bi-Directional with Broadcast Download
 LIS Connection Type RS-232 or TCP/IP
 Microplate Reader Filter Wavelengths 405, 450, 550, 620 nm

Ordering Information

Catalog No.	Description	
426-0400	PhD™ Ix System (110/220V) (includes PhD™ Ix workstation with IFA and EIA hardware, computer, monitor, printer, microplate reader, bar code scanner, surge protector, and Ethernet switch).	1 unit
426-0401	PhD™ Ix Workstation with IFA and EIA hardware (to add to existing PhD™ Ix Systems).	1 unit

Consumables

Catalog No.	Description	
426-0138	PhD™ IFA Conjugate Reagent Tubes and Caps (4/pk).	1 package
426-0139	PhD™ IFA Sample Cups and Caps (20/pk)	1 package
426-0219	Uncoated microplate	1 unit
426-0220	PhD™ Clear EIA Reagent Bottle (30mL, 24/pk)	1 package
426-0221	PhD™ Amber EIA Reagent Bottle (30mL, 8/pk)	1 package



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/diagnostics

Clinical Diagnostics Group

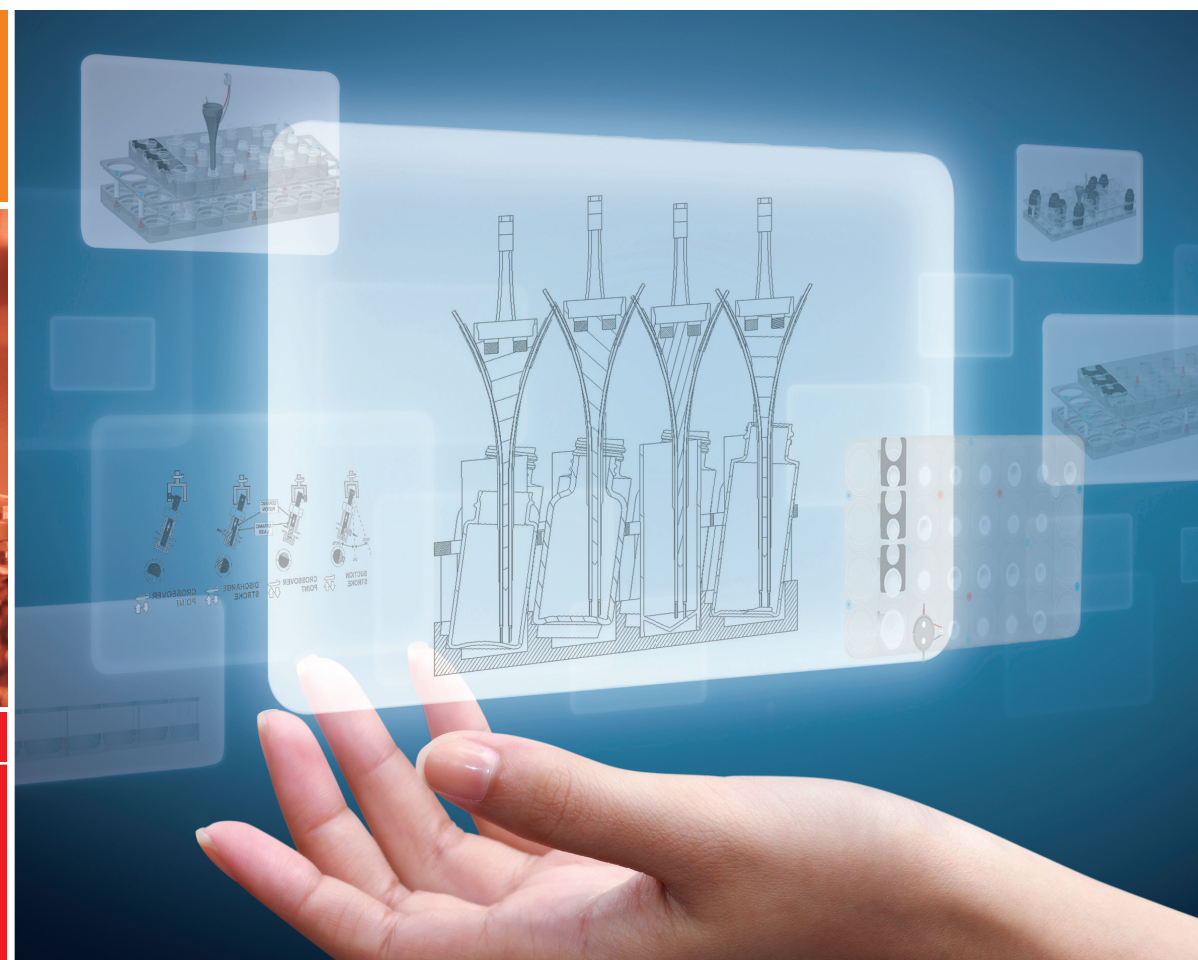
Website www.bio-rad.com/diagnostics U.S. 1-800-224-6723 Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3237-9400 Canada 1-514-334-4372 China 86-21-61698500 Czech Republic 420-241-430-532 Denmark +45-4452-1000 Finland 358-9-804-22-00 France 33-1-47-95-60-00 Germany +49-089-318-840 Greece 30-210-7774396 Hong Kong 852-2789-3300 Hungary +36-1-459-6100 India 1800-180-1224 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico +52 (55) 5488-7670 The Netherlands +31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-3319999 Portugal 351-21-472-7700 Russia 7-495-721-14-04 Singapore 65-6415-3170 South Africa 27-11-442-65-08 Spain 34-91-590-5200 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Taiwan 886-2-2578-7169 Thailand 662-651-8311 United Kingdom +44-020-8328-2000



PhD™ Ix System

IFA and EIA Instrumentation





Enlightened Engineering

Bio-Rad engineers implemented precise magnetic rack positioning on the PhD™ Ix System. What can it do for you? Fewer probe alignments and quicker clean ups, to start. It adds security, too, with alternated polarities so racks can only go in the right way.

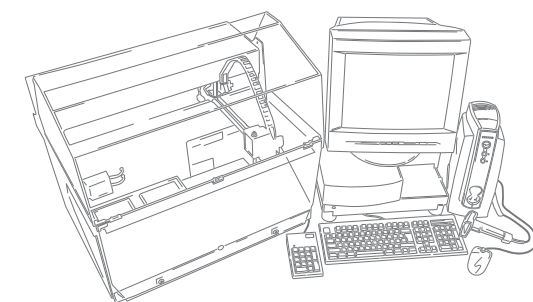


Brighter than ever

Enjoy all the benefits of automation and the confidence of proven reliability. With streamlined workflows, versatile assay programming options and enhanced security features, the PhD™ Ix System provides enlightened solutions for the modern laboratory.

Trusted technology

Laboratories have always loved the reliable performance that comes from the simple, yet powerful PhD™ System. From the proprietary volumetric pump technology to the simple, robust design, the strengths of the original PhD™ System shine through in the PhD™ Ix System.



The Original PhD™ System

Simple, yet powerful

- EIA & IFA processing in one platform
- Network workstations for higher throughput
- Bi-Directional LIS
- Flexible custom assay programming options

Streamlined workflow

Beyond automation, the PhD™ Ix System saves “hands-on” time and reagents, too.

- Instantly create worklists from LIS test orders with broadcast download capabilities
- Align the probe weekly with magnetically positioned racks
- Start runs faster with automatic air bubble removal
- See cleaner IFA backgrounds with dual-probe hyperwashing
- Eliminate dilution plates with in-well dilutions down to 1µL of sample

Versatile

When adding automation, the flexibility to adapt new assays and to bring in send-out testing is a big plus. The PhD™ Ix System has the versatility to fit your needs.

- Automate a wide variety of assays with the flexible Assay Programming Wizard
- Protect sensitive assays from light, dust, and drafts with the dark cover
- Load kit vials directly into the versatile reagent rack
- Easily send IFA & EIA worklists to your laboratory's middleware application

Secure IFA/EIA Processing

Automating routine tasks brings a level of confidence and reproducibility to IFA/EIA processing. The PhD™ Ix System backs up that confidence with robust security features.

- Apply digital signatures with supervisor-level authority
- Track every action with Audit Trails
- Organize reagent caps with the Cap Valet to prevent contamination

