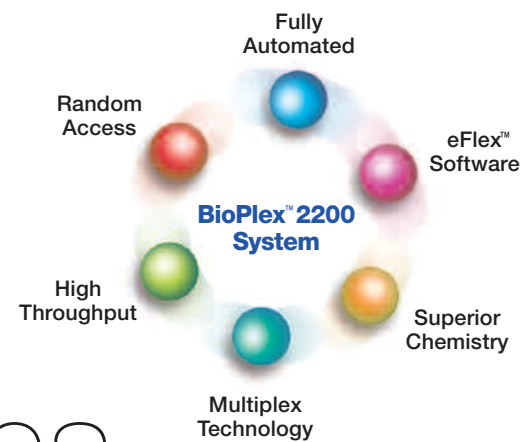
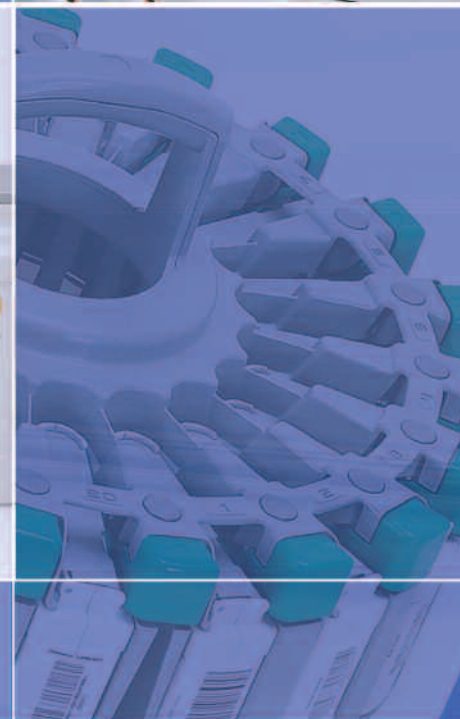




BioPlex™ 2200 System

Like no other. The first and only fully-automated, random access, multiplex testing platform

BIO-RAD



The BioPlex 2200 combines the diagnostic power of proprietary multiplex chemistry, a state-of-the-art software package, and the benefits of full automation and random access capability...all brought together to bring a new level of flexibility and operational benefits to your laboratory.

System Benefits

Fully-Automated/Random Access

- Improved workflow efficiency
- Decreased hands-on time
- Improved turnaround time for test results
- Increased testing capacity
- Fewer errors from specimen and reagent handling

State-of-the-Art eFlex™ Software

- Intuitive, single window, touchscreen interface
- “Add-on” new test orders for previously run patient samples
- LIS connectivity
- Remote diagnostics for faster service response and maximum uptime
- Algorithm-driven interpretive software modules (assay dependent)

Innovative Multiplex Chemistry

- Quality assurance with integrated QC technologies in every test
- Proprietary magnetic bead chemistry for superior performance
- Liquid, ready-to-use reagents, calibrators and controls

Forget limitations.

Leveraging the power of multiplex testing, the BioPlex 2200 brings limitless workflow improvement to your laboratory. You've heard about the promise and future potential of multiplex technology...now Bio-Rad has made that promise a reality.

Disease Focused Multiplex Panel Menu

Currently, multiplex panels are in development for the following disease-state areas:

- Autoimmune
- Serology/Infectious Diseases
- Cardiac
- Diabetes
- Toxicology

Leveraging the Power of Fully-Automated, Random Access, Multiplex Testing

Laboratory and Operator Benefits

- Dramatically reduces labor and turn-around time for specialized testing areas of the lab
- Reduction of errors through automation of manual methods and direct connection with the LIS
- Lower total cost due to vendor and platform consolidation

Physician and Patient Benefits

- Enhanced confidence in results through error reduction and built-in, multiplex, QC checks
- Faster turn-around time for results leads to earlier treatment decisions
- Valuable test "add-on" feature allows for follow-up test orders without a patient redraw

Patented Multiplex Detection Magnetic Bead Technology

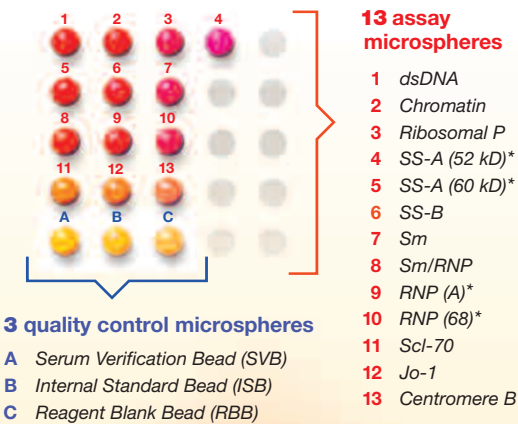
BioPlex™ 2200 immunoassays use heterogeneous sets of magnetic 8 µm beads. The beads are infused with varying ratios of fluorescent dyes, creating unique bead sets. Beads within each set are coated with a ligand (i.e., antigen, antibody, analyte, etc.) specific to a particular assay. Bead sets are then mixed together in a single reagent pack allowing for simultaneous detection of multiple analytes from a single sample.

Each microsphere is dyed with a unique combination of 2 colors



Example: ANA Screen Kit

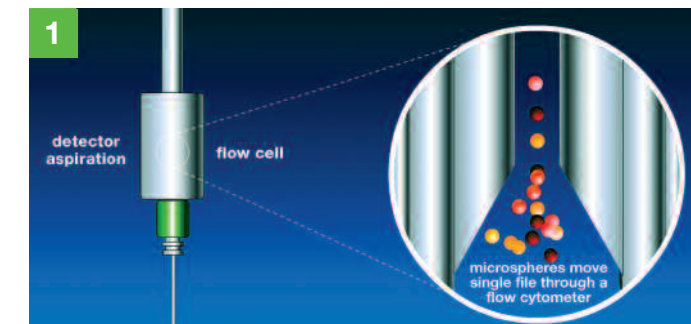
The ANA Screen Reagent Pack uses 16 microspheres: 13 assay beads and 3 quality control checks



*Individual SS-A and RNP results not available in the U.S. Reported as SS-A and RNP for a total of 11 individual results.

Detection Method

The dual-laser, multiplex flow detection method processes a minimum of 150 individual bead results for each analyte, representing more than 150 assay readings. This enhances the overall quality of each result, when compared to standard immunoassays which rely on a single reading.



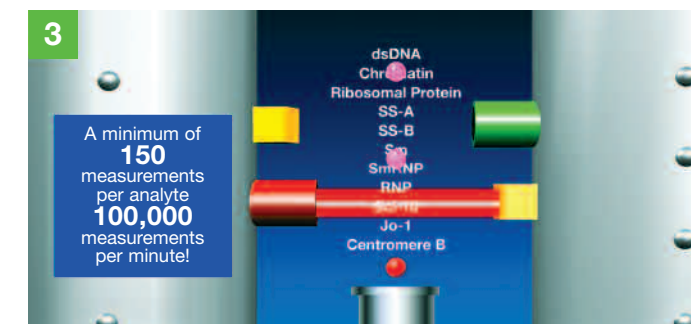
Detection 1

A sample of the final reaction mixture is aspirated to the detector.



Detection 2

Bead mixture passes through the detector where one laser identifies the bead analyte and another measures the analyte concentration.



Detection 3

The system detector performs a minimum of 150 measurements per analyte at the rate of 100,000 measurements per minute.



Quality Assurance with IQ Bead™ Technology

Every test conducted with the BioPlex™ 2200 also features up to three remarkable internal quality control technologies that dramatically improve the quality and accuracy of assay results:

- **Serum Verification Bead (SVB)**
Confirms the presence of serum/plasma
- **Internal Standard Bead (ISB)**
Appears in every tube to standardize detector performance
- **Reagent Blank Bead (RBB)**
Identifies sample problems arising from non-specific binding

BioPlex™ 2200 System

System Specifications

Physical Dimensions

Height 4'-5" (1.3 m)
 Width 4'-3" (1.3 m)
 Depth 2'-10" (0.9 m)
 Weight Approximately 1032 lbs. (468 kg)

Performance Characteristics

Maximum Throughput 22 results per sample at 100 samples per hour, yielding up to 2200 results per hour
 Time to First Result 20-45 minutes (depending on panel)
 Sample Tubes Any CLSI primary sample tubes
 Sample Racks BioPlex™ 2200 racks or Hitachi™ racks
 Sample Volume 5 - 50 µL (depending on panel)
 STAT Capability Continuous Access
 Walkaway Capacity 8 hours of continuous use
 Operational Mode Random and continuous access, priority (STAT) loading
 Operator Interfaces Touch screen monitor, keyboard, trackball mouse
 Aspiration Capability Robotic precision with clot detection
 Quality Control Levey-Jennings, Westgard Rules
 Sample Barcode Acceptable Symbolologies Code 128, Codabar, Code 39, Interleaved 2 of 5
 Sample Types Serum or Plasma For more information, refer to BioPlex™ 2200 assay insert
 Sample Dilutions 1:4, 1:10, and 1:100 (depending on panel)
 Access During Use Continuous sample input/output, consumables replenishment, and waste disposal without interruption of sample processing
 LIS Connectivity Bi-directional: ASTM Standard Class 2 Laser Product

Power Requirements

Input Voltage 100-240 VAC at 50-60 Hz, 1000 Watt
 UPS Optional

Water Requirements

Water Quality CLSI Type II or better
 Direct Plumbing Maximum pressure of 50 kPa

Installation Requirements

Crate Clearance Requirements

Rear 5' (1.5 m)
 Left 4' (1.2 m)
 Right 4' (1.2 m)
 Front 10' (3.0 m)
 Vertical 6' (1.8 m)

Instrument Operational Clearance

Rear 6" (0.2 m)
 Left 3' (1.0 m)
 Right 6" (0.2 m)
 Front 3' (1.0 m)
 Vertical 6' (1.8 m)



Ordering Information

Catalog No. Description

660-0000	BioPlex™ 2200	1 instrument
660-0800	Sample Racks (30 per box)	1 box
660-0801	Sample Tray	1 tray
660-0817	Instrument Sheath Fluid (2 per box)	1 box
660-0818	Instrument Wash Buffer	1 box
660-2003	Reaction Vessels (1000 per bag)	1 bag
666-0001	Instrument Detector Calibration Pack	1 pack
666-0002	Instrument Detector Clean Pack	1 pack
666-0003	Instrument Probe Cleaning Solution Bottles (4 per box)	1 box



**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.** 1-800-2BIO-RAD **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** 86-21-64260808 **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-(0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** 36-1-455-8800 **India** 91-124-4029300 **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** 65-6415-3188 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200
Sweden 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000