

An underwater scene featuring a vibrant coral reef on the right side, populated with numerous small, orange-colored fish. The water is a deep blue, with some light rays visible in the upper left corner. The overall composition is dynamic and colorful.

# Bio-Plex™

System and Technology

A World of Information  
From a Single Drop

**BIO-RAD**

## Introduction to the Bio-Plex System

The Bio-Plex suspension array system is a flexible, easy-to-use bioassay system for the simultaneous detection and quantitation of up to 100 different analytes in a single microplate well. Multiplexing with the Bio-Plex system yields data that are linked within a system so that complex relationships and pathways of biomolecules can be revealed.

### Multiplex With a Single Drop

- Use as little as 12 µl of serum or other biological sample per multiplex assay
- Dramatically increase the amount of useful data per sample
- Increase sample throughput
- Directly correlate data from multiple analytes to reveal complex relationships and pathways of biomolecules
- Perform immunoassays, enzyme assays, receptor-ligand assays, DNA hybridization assays, and more

# Unsurpassed Multianalyte Detection

## System Integration for Optimal Results

The Bio-Plex system is designed, manufactured, and tested as a complete, fully integrated system combining hardware, Bio-Plex Manager™ software, system validation and calibration tools, assays, and beads. The fully integrated and validated system ensures accurate and reproducible assay results. It also gives you access to xMap technology and all these other benefits:

- Standardized single-source system components for error-free operation
- Simplified operation with automated startup, shutdown, validation, and calibration functions
- Enhanced dynamic range and recovery with immunoassays
- Powerful dedicated data reduction software
- Easy-to-use reports
- Installation and operation qualification (IQ/OQ) certification

## Support That Maintains Your Productivity

The integrated design of the Bio-Plex system means that a single source provides all the support you need.

- Worldwide technical support and on-site field service
- Service agreements for scheduled preventive maintenance



### Bio-Plex Assays and Reagents

Ready-to-run Bio-Plex assay kits are available in premixed multiplex and singleplex configurations.

#### Assay kits

- Cytokines
- Phosphoproteins

#### Assay reagents

- Amine coupling kit
- COOH beads

### Bio-Plex System Software

The only integrated system control and data analysis package for xMAP technology.

#### Unique features

- StatLIA 4PL and 5PL curve fitting
- Automated functions
- Percentage recovery calculations

### High-Throughput Fluidics (HTF) System

The Bio-Plex HTF system allows the processing of microplates without refilling system fluid containers after each run.

# and Quantitation

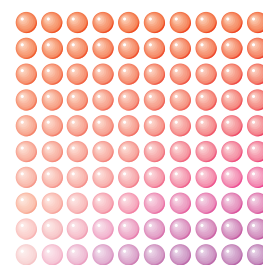
## The Technology

The Bio-Plex array system utilizes xMAP technology to permit the multiplexing of up to 100 different assays within a single sample. The system uses a liquid suspension array of 100 sets of 5.5 µm beads, each internally dyed with different ratios of two spectrally distinct fluorophores to assign it a unique spectral address (Figure 1). Each set of beads can be conjugated with a different capture molecule; the conjugated beads can then be mixed and incubated with sample in a microplate well to react with specific analytes. Capture molecules can include:

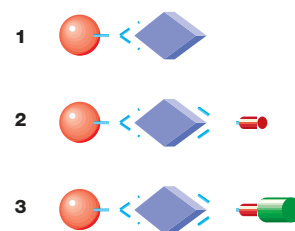
- Enzyme substrates
- DNA
- Receptors
- Antigens
- Antibodies



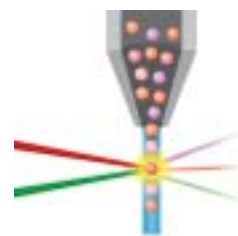
To detect and quantitate each captured analyte, a fluorescently labeled reporter molecule that specifically binds the analyte is added. Figure 2 illustrates use of a capture molecule for a Bio-Plex sandwich immunoassay. Following incubation, the contents of each microplate well are drawn into the Bio-Plex array reader, and precision fluidics align the beads in single file through a flow cell where two lasers excite the beads individually (Figure 3). The red classification laser excites the dyes in each bead, identifying its spectral address. The green reporter laser excites the reporter molecule associated with the bead, which allows quantitation of the captured analyte. High-speed digital signal processors and Bio-Plex Manager software record the fluorescent signals simultaneously for each bead, translating the signals into data for each bead-based assay.



**Fig. 1.** Bio-Plex assays are designed using microscopic beads, each with a different color code, or spectral address, to permit discrimination among multiplex assays.



**Fig. 2.** In a typical Bio-Plex assay, a capture molecule conjugated to a color-coded bead binds to a target analyte (1) followed by binding with biotinylated detection antibody (2) and a reporter molecule, streptavidin-PE (3).



**Fig. 3.** In the Bio-Plex array reader, a red classification and a green reporter laser illuminate individual beads to identify each bead's spectral address and associated reporter signal.

### Bio-Plex Array Reader and Microplate Platform

The flow-based, dual-laser array reader classifies each bead and its associated assay, and quantitates the amount of analyte captured in each assay. The microplate platform permits automated processing of samples from 96-well microplates in ~30 min.

### Bio-Plex Validation and Calibration Tools

To ensure accurate and reproducible assay results, the Bio-Plex system includes tools for validation of system fluidics, the reporter and classification channels, laser alignment, and daily system calibration.

## Bio-Plex System Specifications

### Bio-Plex Array Reader

Operating temperature	15–30°C (59–86°F)
Humidity	10–90%, noncondensing
Input voltage range	100–240 V, 1.5 A, 47–63 Hz
<b>Lasers</b>	
Reporter laser	532 nm, >10 mW frequency-doubled diode; 30 x 60 µm elliptical beam
Classification laser	635 nm, 10 mW diode; 30 x 60 µm elliptical beam

### Fluidics

Sheath flow rate	90 µl/sec
Cuvette pathlength	200 µm
Sample injection rate	60 µl/min

### Electronics

Reporter channel detection	Photomultiplier tube, A/D resolution 14 bits
Classification and doublet discriminator channel detection	Avalanche photodiodes with temperature compensation, A/D resolution 12 bits
Communications interface	RS-232

### Signal processing

Measurement resolution	15 bits effective
Processor modes	Linear, with logarithmic or linear display option

### Dynamic range

Dimensions (W x D x H)	43 x 51 x 23 cm (17 x 20 x 9")
Weight	23 kg (60 lb)

### Bio-Plex Microplate Platform

Operating temperature	15–30°C (59–86°F)
Humidity	10–90%, noncondensing
Input voltage range	100–240 V, 0.4 A, 47–63 Hz
Communication interface	RS-232
Plate capacity	One 96-well microplate ≤2 cm (0.75") thick
Dimensions (W x D x H)	44 x 61 x 8 cm (17.3 x 24 x 3")
Weight	14.4 kg (32 lb)

### Bio-Plex High-Throughput Fluidics (HTF) System

Operating temperature	15–30°C (59–86°F)
Humidity	20–80%, noncondensing
Input voltage range	100–240 V, 0.4 A, 47–63 Hz
Dimensions (W x D x H)	20 x 30 x 25 cm (8 x 12 x 10")
Weight	9 kg (20 lb)

## Ordering Information

Catalog #	Description
<b>Bio-Plex System*</b>	
171-000001	<b>Bio-Plex Suspension Array System</b> , 100–240 V
171-000005	<b>Bio-Plex Suspension Array System</b> , 100–240 V, with high-throughput fluidics (HTF) system

\* Bio-Plex systems include array reader, microplate platform, PC, monitor, Bio-Plex Manager software, calibration kit, validation kit, maintenance, calibration, and validation (MCV) plate II, 20 L sheath fluid, instructions

### Bio-Plex System Accessories

171-000050	<b>High-Throughput Fluidics (HTF) System</b> , 100–240 V
171-000055	<b>Sheath Fluid</b> , 20 L
171-002001	<b>Communication Cable</b> , 5 ft
171-002010	<b>Sheath Fluid Bottle</b> , 1 L
171-002012	<b>Sheath Waste Bottle</b> , 2 L
171-002020	<b>Sample Needle</b> , long (11.6 cm/4.6")
171-002023	<b>Needle Guide</b>
171-002026	<b>Needle Adjustment Wrench</b>
171-002030	<b>Protective Shield for Sample Needle</b>

### Bio-Plex Validation and Calibration Accessories

171-203000	<b>Bio-Plex Validation Kit 3.0</b> , includes optics validation, reporter validation, classify validation, and fluidics validation bead sets for approximately 50 validation routines
171-203031	<b>Bio-Plex MCV (Maintenance, Calibration, and Validation) Plate II</b>
171-203060	<b>Bio-Plex Calibration Kit</b> , includes CAL1 and CAL2 calibration beads for approximately 50 daily calibration routines

StatLIA is a trademark of Brendan Scientific Corp. xMAP is a trademark of Luminex Corp.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191  
Canada (905) 712-2771 Czech Republic + 420 2 41 43 05 32 China (86-21) 63052255 Denmark 44 52 10 00 Finland 09 804 22 00  
France 01 47 95 69 65 Germany 089 318 84-177 Hong Kong 852-2789-3300 India (91-124)-6398112/113/114, 6450092/93 Israel 03 951 4127  
Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 52 5 534 2552 to 54  
The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 8126 672 Portugal 351-21-472-7700  
Russia 7 095 721 1404 Singapore 65-62729877 South Africa 00 27 11 4428508 Spain 34 91 590 5200 Sweden 08 555 12700  
Switzerland 061 717-9555 Taiwan (8862) 2578-7189/2578-7241 United Kingdom 020 8328 2000