



BioPlex[®] 2200 System Profile



Sample Processing Module

- Processes reaction vessels with patient sample and multiplex reagents
- Inner ring for 37°C incubations
- Outer ring for magnetic wash and separations
- Throughput capacity of up to 100 multiplexed patient samples per hour

Detector Module

- Multiplex flow detection
- Minimum of 200 measurements per analyte

Sample Input Area

- Holds up to 100 primary samples
- Continuous, random access sample loading

STAT Sample Entry

- On demand STAT capability

On-board Bulk Reagent Storage

- On-the-fly loading and unloading of bulk reagents and waste
- Bulk reagent and waste capacity to process up to 800 primary samples



Reagent Carousel

- On-board refrigerated storage
- Store up to 20 reagent packs giving the capacity to process up to 2,000 samples
- Real time inventory management

eFlex™ Software

- Intuitive touch screen operation
- Flexible worklist, testing, and results reporting set-up
- Ability to “Add on” additional test orders for previously run samples
- Bi-directional LIS interface
- Medical Decision Support Software (MDSS)

Sample Output Area

- Holds up to 100 primary samples
- Continuous access sample unloading

Central Processing Computer

- CPU for instrument and software communication

Reaction Vessel Storage/Waste

- Above – hopper for storage of up to 800 reaction vessels
- Below – waste bin storage for used reaction vessels

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	Maximum 45 minutes (depending on panel)
Sample Tubes	Any CLSI primary sample tubes
Sample Racks	BioPlex® 2200 racks or Hitachi™ racks
Sample Volume	Requirements Approx. 5 µL (depending on sample tube)
STAT Capability	1 rack of 5 sample tubes
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 Symbologies	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, 1:100
Access During Use	Continuous sample input/output, consumables replenishment, and waste disposal without interruption of sample processing
LIS Connectivity	Bi-directional: ASTM Standard

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)



BIO-RAD

**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