



# BioPlex<sup>®</sup> 2200 System

U.S. Operator Training Program



## U.S. Operator Training Program

### Description

The course is designed to:

- Provide the skills required to operate the instrument on a daily basis
- Setup and initiate assay run
- Present procedures and techniques to properly maintain the instrument
- Learn techniques to interpret results
- Learn skills applicable to troubleshooting

### Who should take this course?

The course is designed:

- For users designated as the primary operators who are not familiar with the platform
- For users designated to train other operators on-site
- As a pre-installation course

### Course Information

#### Participants:

A minimum of two participants is required to conduct a class. Enrollment is limited to six participants to allow for maximum hands-on instructions.

**Fee:** (Training fee is included with new instrument installations)

\$5,000 including lodging and meals. Other expenses, including gratuities, are the responsibility of each individual. Airfare not included (San Francisco).



Course Outline				
DAY ONE	DAY TWO	DAY THREE	DAY FOUR	DAY FIVE
<ol style="list-style-type: none"> <li><b>1 Introduction</b> <ul style="list-style-type: none"> <li>• Multiplex testing</li> </ul> </li> <li><b>2 System Overview</b> <ul style="list-style-type: none"> <li>• Hardware</li> <li>• Software</li> </ul> </li> <li><b>3 Review and Questions</b></li> </ol>	<ol style="list-style-type: none"> <li><b>1 System Overview</b> (continued)                     <ul style="list-style-type: none"> <li>• Software</li> <li>• Chemistry</li> </ul> </li> <li><b>2 Assays</b> <ul style="list-style-type: none"> <li>• Methods available</li> <li>• Assay preparation                             <ul style="list-style-type: none"> <li>– Reagent packs</li> <li>– Calibration</li> <li>– Quality Control (QC)</li> <li>– Initiate calibration and QC assay run</li> </ul> </li> </ul> </li> <li><b>3 Review and Questions</b></li> </ol>	<ol style="list-style-type: none"> <li><b>1 Assays</b> (continued)</li> <li><b>2 Initiate Sample Run</b></li> <li><b>3 Result Review</b> <ul style="list-style-type: none"> <li>• Add-on tests</li> <li>• Review                             <ul style="list-style-type: none"> <li>– Repeat</li> <li>– Dilution</li> <li>– Releasing results</li> </ul> </li> </ul> </li> <li><b>4 Review and Questions</b></li> </ol>	<ol style="list-style-type: none"> <li><b>1 Maintenance</b> <ul style="list-style-type: none"> <li>• Daily</li> <li>• Weekly</li> <li>• Monthly</li> <li>• On demand</li> </ul> </li> <li><b>2 Troubleshooting</b></li> <li><b>3 Review and Questions</b></li> </ol>	<ol style="list-style-type: none"> <li><b>1 Course Exams</b> <ul style="list-style-type: none"> <li>• Written Exam</li> <li>• Wet Lab Exam</li> </ul> </li> <li><b>2 Review and Questions</b></li> </ol>



Bio-Rad Laboratories

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

Clinical Diagnostics Group

Website [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) U.S. 1-800-2BIO-RAD (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 86-21-64260808 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 +36-1-459-6100 India 1-800-180-1224 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 27-11-442-85-08 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