



# 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

### 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).

### 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 Outline				
<p><b>DAY ONE</b></p> <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>	<p><b>DAY TWO</b></p> <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> </ul> </li> <li>– Initiate calibration and QC assay run</li> </ul> </li> <li><b>3 Review and Questions</b></li> </ol>	<p><b>DAY THREE</b></p> <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> </ul> </li> <li>– Releasing results</li> </ul> </li> <li><b>4 Review and Questions</b></li> </ol>	<p><b>DAY FOUR</b></p> <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>	<p><b>DAY FIVE</b></p> <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