



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



**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 (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