



VARIANT™ II

Operator Training Program



VARIANT™ II Customer Training Program

Description

The course is design 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
- Users designated to train other operators on-site
- New users

Course information

Participants:

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

Fee:

\$3,000 including lodging and meals. Other expenses, including gratuities, are the responsibility of each individual. Airfare not included.



Course Outline		
<p>DAY ONE</p> <ol style="list-style-type: none"> 1 Introduction <ul style="list-style-type: none"> • Disease States 2 System Overview <ul style="list-style-type: none"> • Hardware • Software • Chemistry 3 Sample Preparation 4 Initiating Assay Run 5 Review and Questions 	<p>DAY TWO</p> <ol style="list-style-type: none"> 1 Result Review 2 Data Management <ul style="list-style-type: none"> • Query • Manual Printouts • Manual LIS Transmission 3 Maintenance <ul style="list-style-type: none"> • Daily • Monthly • Periodic 4 Troubleshooting <ul style="list-style-type: none"> • Pressure Issues • Hardware • Chromatography 5 Review and Questions 	<p>DAY THREE</p> <ol style="list-style-type: none"> 1 Operational Exercises 2 Troubleshooting Review and Exercise 3 Clinical System Network™ Application <ul style="list-style-type: none"> • Overview • Navigation • Operation • Remote Connectivity 4 Course Exam



**Bio-Rad
Laboratories**

For further information and registration, please contact your local Bio-Rad representative at 1-800-224-6723.

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3461-5202 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-64260808 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Mexico 5255-5534-2552 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 7-095-721-14-00 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 44-208-328-2000