



# VARIANT™ II TURBO

Operator Training Program

**BIO-RAD**

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



**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](http://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