



VARIANT™ nbs

Operator Training Program



VARIANT™ nbs Customer 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
- Users designated to train other operators on-site
- New users

Course information

Participants:

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

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



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