



# EDUCATION SERVICES TO EMPOWER YOUR LAB

**OneSource**  
Laboratory Services

Customer Classroom Training Program 2019

  
**PerkinElmer**  
For the Better



# YOUR BENEFIT FROM TRAINING

Our classroom training offers an ideal setting for learning how to

get the best possible performance from your instrument. Free from everyday distractions, you'll receive hands-on experience in a laboratory environment complete with the latest equipment, software and accessories. You'll receive guidance from expert faculty with comprehensive instrument knowledge and experience. You'll also have the opportunity to network with other participants to problem solve.

Training courses are offered at PerkinElmer Centers of Excellence, conveniently located throughout the United States and Canada. We maintain a small class size to ensure you receive individual attention and ample hands-on instrument use – everything you need for focused learning that can lead to increased productivity.

Participants are accepted on a first-come, first-served basis, so reserve your spot soon.

## UNABLE TO GET AWAY FROM YOUR LAB?

Contact us to discuss bringing the training directly to you – saving you time and travel expenses. Email [OnsiteTraining@perkinelmer.com](mailto:OnsiteTraining@perkinelmer.com) for more information.

## Table of Contents

2019 Calendar	3
<b>Atomic Spectroscopy</b>	4
Atomic Absorption	4
Inductively Coupled Plasma	5
Inductively Coupled Mass Spectrometry	6
<b>Chromatography</b>	7
Liquid Chromatography	7
Chromatography Data Systems	7
Gas Chromatography	8
Gas Chromatography Mass Spectrometry	9
LC Mass Spectrometry	10
<b>Molecular Spectroscopy</b>	11
Infrared Spectroscopy	11
<b>Thermal/Elemental Analysis</b>	12
Thermal Analysis	12
Locations	13
Enrollment Information	13
Course Agreement	13
Registration Document	14

# CUSTOMER CLASSROOM TRAINING PROGRAM - 2019

## 2019 CALENDAR

ATOMIC SPECTROSCOPY									
Course Name	Course No.	Days	Price	Atlanta, GA	Chicago, IL	San Jose, CA	Shelton, CT	Toronto, CAN	
<b>ATOMIC ABSORPTION (AA)</b>									
Flame Atomic Absorption	N0205001	2	\$1,450.00	FEB 11-12	JUN 24-25		APR 8-9		
Graphite Furnace Atomic Absorption	N0200017	3	\$2,050.00	FEB 13-15	JUN 26-28		APR 10-12		
Advanced Furnace AA	N0200024	2	\$1,450.00	Call for Availability					
<b>INDUCTIVELY COUPLED PLASMA (ICP-OES)</b>									
Inductively Coupled Plasma	N0205010	3	\$2,050.00	JAN 8-10 MAR 4-6 MAY 7-9	JAN 28-30 APR 2-4 JUN 25-27	JAN 9-11 FEB 20-22 APR 23-25 JUN 5-7	JAN 15-18 MAR 12-14 APR 23-25 JUN 17-19	MAR 27-29 JUN 11-13	
Advanced ICP Training	N0200027	2	\$1,450.00	MAR 7-8	JAN 31-FEB 1		JUN 20-21		
<b>ICP MASS SPECTROMETRY (ICP-MS)</b>									
NexION® ICP-MS	N0200193	5	\$3,150.00	JAN 28-FEB 1 MAR 25-29 APR 29-MAY 3 JUN 3-7	FEB 4-8 MAY 20-24 JUN 17-21	JAN 14-18 MAR 18-22 MAY 6-10 JUN 10-14	FEB 4-8 MAR 4-8 MAY 13-17 JUN 24-28	APR 1-5	
LC/ICP-MS Instrumentation-Metal Speciation	N0200522	3	\$2,050.00	Call for Availability					
<b>CHROMATOGRAPHY</b>									
Course Name	Course No.	Days	Price	Atlanta, GA	Chicago, IL	San Jose, CA	Shelton, CT	Toronto, CAN	
<b>GAS CHROMATOGRAPHY (GC)</b>									
Fundamentals of Gas Chromatography	N0200407	3	\$2,050.00	FEB 6-8	MAY 8-10	JUN 5-7	APR 3-5		
<b>GC MASS SPECTROMETRY (GC/MS)</b>									
Clarus® GC/MS Operator	N0200417	4	\$2,700.00	FEB 11-14	MAY 13-16	JUN 10-13	APR 8-11		
Torion® GC/MS Basic Operations	N0239010	3	\$2,050.00	Call for Availability					
Torion® GC/MS Advanced Operations	N0239011	5	\$3,150.00				AUG 13-17		
Torion® GC/MS Advanced Operations and Maintenance for Distributors	N0239012	10	\$6,300.00				Call for Availability		
<b>LIQUID CHROMATOGRAPHY (LC)</b>									
Altus™ LC Training	N0233024	2	\$1,450.00						
Fundamental LC	N0200501	2	\$1,450.00						
<b>LC MASS SPECTROMETRY (LC-MS)</b>									
AxION® 2 DSA	N0233019	3	\$2,050.00						
AxION 2 LC/DSA	N0233020	5	\$3,150.00						
AxION 2 LCMS	N0233022	3	\$2,050.00						
QSight® Triple Quad LC/MS/MS	N0233042	5	\$3,150.00				FEB 25-MAR 1		
<b>CHROMATOGRAPHY DATA SYSTEMS (CDS)</b>									
Empower® Quickstart Software	N0233025	3	\$2,050.00				Call for Availability		
TotalChrom® Principles	N0201043	3	\$2,050.00	MAR 12-14					
<b>MOLECULAR SPECTROSCOPY</b>									
Course Name	Course No.	Days	Price	Atlanta, GA	Chicago, IL	San Jose, CA	Shelton, CT	Toronto, CAN	
<b>INFRARED SPECTROSCOPY (IR)</b>									
AssureID™ Software	N0200188	2	\$1,450.00						
Intro to FT-IR Spectroscopy	N0200126	2	\$1,450.00	APR 9-10		FEB 25-26			
Spectral Interpretation	N0200152	2	\$1,450.00	APR 11-12		FEB 27-28			
<b>THERMAL/ELEMENTAL ANALYSIS</b>									
Course Name	Course No.	Days	Price	Atlanta, GA	Chicago, IL	San Jose, CA	Shelton, CT	Toronto, CAN	
Intro to Differential Scanning Calorimetry	N0200607	2	\$1,450.00						
Intro to Thermogravimetric Analysis	N0200605	2	\$1,450.00				Call for Availability		



## Atomic Absorption

N020-5001 | 2 Days | Tuition \$1,450

### Flame Atomic Absorption

This two-day course provides the analyst with the knowledge and skills needed for optimizing and troubleshooting basic flame atomic absorption. Laboratory exercises include experiments in controlling interferences as well as procedures for developing methods. WinLab and Syngistix™ software will also be covered.

#### Prerequisite

The analyst should have spent at least one month becoming familiar with the instrument and should have a working knowledge of Microsoft® Windows and AA WinLab software before attending this course.

#### Lectures and Labs

- Hardware optimization and calibration
- Setting up the element parameter file
- Maintenance procedures
- Background correction
- Flame emissions analysis
- Interferences

N020-0017 | 3 Days | Tuition \$2,050 (if not included w/instrument purchase)

### Graphite Furnace Atomic Absorption

This three-day course will provide the analyst with the knowledge necessary to set up, run, troubleshoot, and operate the graphite furnace. A considerable portion of this course is devoted to understanding the capabilities of the WinLab and Syngistix software used to operate the system. The course begins with the basics of graphite furnace atomic absorption (GFAA) and concludes with the most recent developments in the technique including simultaneous graphite furnace analysis.

#### Prerequisite

The student should have spent at least one month becoming familiar with the instrument, and should have a working knowledge of Microsoft® Windows and the AA WinLab software before attending this course.

#### Lectures and Labs

- Introduction to GFAA
- Setting up the element parameter file
- Interferences
- Matrix modifiers
- Background correction systems
- Quality control
- Method development
- Optimization of simultaneous analysis parameters (SIMAA)
- Troubleshooting and maintenance

N020-0024 | 2 Days | Tuition \$1,450

### Advanced Furnace AA

This two-day laboratory intensive course, which will focus on how to handle difficult samples. The modifier and its role in solving analytical problems will be discussed in detail. The class will examine complex matrices, focusing on peak plot interpretation as a means to method development. Techniques such as QA/QC as a means of verifying the accuracy of data will be discussed. Class participants will be expected to have a good working knowledge of the software as this will not be addressed in the course. This course is not for the novice. The basic techniques of graphite furnace and an extensive knowledge of the Microsoft® Windows-based software will be expected.

#### Prerequisite

The student should have completed course N020-0017 (basic Graphite Furnace) or have a minimum of six months experience with graphite furnace analysis and the WinLab software before attending this course.

#### Lectures and Labs

- Method optimization
- Quality control for method validation
- Peak plot interpretation
- Solid sampling
- Understanding modifiers

## Inductively Coupled Plasma



N020-5010 | 3 Days | Tuition \$2,050 (if not included w/instrument purchase)

### Inductively Coupled Plasma Series

This three-day comprehensive course comprises a study of basic emission theory, radial versus axial viewing, ICP hardware, and WinLab 32 and Syngistix software. All lab experiments will be conducted on current models of the PerkinElmer ICP series.

#### Prerequisite

Analysts should have spent at least one month becoming familiar with the instrument and should have a working knowledge of WinLab 32 and/or Syngistix software before attending this course.

#### Lectures and Labs

Theory of atomic emission	Method development
Plasma performance	Background correction
Identifying and overcoming interferences	Multicomponent Spectral Fitting (MSF)
ICP component	Interfering Element Correction (IEC)
Data reprocessing	Echelle optical system maintenance
Radial vs. axial viewing	Parameter optimization

N020-0027 | 2 Days | Tuition \$1,450

### Advanced Optima ICP

This two-day course provides analysts with the knowledge needed to develop a method for ICP-OES analysis of complex samples. This lab intensive course covers when and how to use internal standards, inter-element correction factors (IECs) and multi-component spectral fitting (MSF) effectively to obtain accurate results.

#### Prerequisite

The student should have completed course N020-5010 (ICP with ICP WinLab and Syngistix software) or have a minimum of six months experience with the Optima/Avio ICP and the WinLab/Syngistix software before attending this course. This course is not recommended for beginners.

#### Lectures and Labs

- Identifying interferences
- Inter-element correction factors
- Using internal standards
- Multicomponent spectral fitting



## Inductively Coupled Mass Spectrometry

N020-0193 | 5 Days | Tuition \$3,150 (if not included w/instrument purchase)

### NexION ICP-MS Instrumentation

This five-day course will cover the NexION® series of ICP-MS. This course will be a combination of classroom lectures and hands on laboratory experiments and exercises. The items covered will consist of the following:

- ICP-MS theory and method development
- Overview and discussion of the hardware of the
- NexION/Syngistix software training
- Discussion on optimization and setup of the instrument
- Introduction into the optimization and the use of
- Universal Cell Technology (UCT)
- Troubleshooting, and maintenance
- Reporting and data exporting

#### Prerequisite

Students should have spent at least one month becoming familiar with the instrument, and should have a working knowledge of the software before attending this course.

*\*An experiment on how to use the autosampler will be covered in class however ESI Autosampler material will NOT be covered.*

Also as part of this course material, appropriate lab exercises will be conducted that are related to the discussions held. Laboratory experiments will be conducted to show how to set up, optimize and maintain the ICP-MS instrument. In addition, data will be collected by means of various analytical methods available with this technique. All lab experiments will be conducted on the NexION models of ICP-MS instruments.

N020-0522 | 3 Days | Tuition \$2,050

### LC/ICP-MS Instrumentation-Metal Speciation

This three-day advanced course will cover the concept and implementation of speciation analysis by coupling the analytical techniques of inductively coupled plasma emission mass spectrometry (ICP-MS) and liquid chromatography (LC). This course will include a one-day primer on LC, a review of ICP-MS optimization and the interfacing of the two techniques.

Also included is basic training on how to use the Chromera® software. Laboratory experiments will be conducted to show how to set up this technique and collect data. This data will be collected using two of the available modes of LC. All lab experiments will be conducted on the current models of the ICP-MS, LC pump and autosampler.

#### Prerequisite

Students should be very familiar with the ELAN series of ICP-MS instrumentation and should have a strong working knowledge of the ELAN software before attending this course. Ideally students should have attended the basic ICP-MS course (course number N020-0008).

#### Lectures and Labs

Introduction to Databases  
 Fundamentals of HPLC  
 Interfacing HPLC to ICP-MS  
 HPLC Hardware Familiarization  
 Selenium Speciation  
 Review of ICPMS Optimization  
 Software (Chromera)  
 HPLC Troubleshooting  
 Computer Based Software Exercises  
 Arsenic Speciation



## Liquid Chromatography

N020-0501 | 2 Days | Tuition \$1,450

### Fundamental Liquid Chromatography

For the beginner, this two-day introduction to liquid chromatography provides an overview of the principles and practice of high-performance liquid chromatography (HPLC). The topics covered are basic chromatographic terminology, modes of HPLC (normal and reversed-phase, size exclusion, and ion exchange chromatography), instrumentation, peak identification and method development, a brief applications review, and routine maintenance and troubleshooting.

#### Prerequisite

The student should have spent at least one month becoming familiar with the instrument.

#### Lectures and Labs

- History and theory of LC
- Modes of HPLC
- Hardware components
- Peak identification and method development
- Applications overview
- Troubleshooting
- Basic chromatographic parameters
- Hardware familiarization
- Peak identification
- Method development
- Routine maintenance

## Chromatography Data Systems (CDS)

N022-3025 | 3 Days | Tuition \$2,050

### Empower Quickstart

This three-day lecture and lab course Presents users of Empower 3 Workstation with a fundamental knowledge of the Quickstart interface. This course is for key operators who need formal product training or operators who need more comprehensive training than that supplied by on-site familiarization courses.

#### Prerequisite

A fundamental understanding and/or experience using Microsoft® Windows® 7 operating system is strongly recommended for this course. 20 to 60 days of product usage before the class is recommended.

#### Lectures and Labs

- Operation of the software, basic maintenance, and troubleshooting.
- Discussion and hands-on exercises emphasize acquiring data
- Developing quantitative methods
- Data processing,
- Basic report generation

N020-1043 | 3 Days | Tuition \$2,050

### TotalChrom Principles of Operation

This three-day lecture and lab course presents users of TotalChrom™ Workstation and Client/Server with a fundamental knowledge of the theory and operation of the software, basic maintenance, and troubleshooting. Discussion and hands-on exercises emphasize acquiring data, developing quantitative methods, and reprocessing data. This course is for key operators of laboratories who need formal product training or operators who need more comprehensive training than that supplied by on-site familiarization courses.

#### Prerequisite

A fundamental understanding and/or experience using the Microsoft® Windows® 95, 98 or 2000 operating systems are strongly recommended for this course. Thirty to sixty days of product usage before the class is recommended.

#### Lectures and Labs

- System overview
- PerkinElmer Intelligent Interface
- Software configuration
- Peak detection and identification
- Generation and calibration of method files
- Data acquisition
- Graphic reprocessing
- Comparing chromatograms
- Batch
- Generation of sequences



## Gas Chromatography

N020-0407 | 3 Days | Tuition \$2,050

### Fundamental Gas Chromatography

This three-day course provides an overview of the basic principles, hardware, and operational techniques used in gas chromatography, along with column overview.

#### Prerequisite

The student should have spent at least one month becoming familiar with the Clarus instrument.

*Note: Students in need of detailed instruction using TotalChrom software should attend the appropriate Chromatography Data Systems course.*

TotalChrom Principles will not be included in this course.

#### Lectures and Labs

Basic principles of GC  
Starting a chromatograph  
Hardware, gases, injectors,  
column installation, flow measurements  
detectors, etc.  
Setup of GC and data handling  
Qualitative and quantitative analysis  
Principles and optimization of  
data-handling parameters

## GC Mass Spectrometry

N020-0417 | 4 Days | Tuition \$2,700 (if not included w/instrument purchase)

### GC-MS Operator

This four-day course introduces new Clarus® GC/MS and users to gas chromatography/mass spectroscopy (GC/MS) and gives them an orientation to the hardware, software, and experimental requirements to accomplish successful system operation. Laboratory exercises provide an opportunity to carry out the main operations described in the lecture material.

#### Prerequisite

Students are required to have completed one of the PerkinElmer gas chromatography training courses, or to certify a history of extensive GC experience. The student will also be required to have at least one month of experience with an installed and operating Clarus MS system.

#### Lectures and Labs

System maintenance and software  
troubleshooting procedures  
Connecting the GC and MS  
Spectral data processing and cold starting  
the instrument  
Library searching  
Tuning the instrument  
Developing a method for construction and  
data analysis  
Instrument control and data acquisition  
Developing a Clarus MS method for  
quantization and a GC control method  
General maintenance

## GC Mass Spectrometry



N020-9010 | 3 Days | Tuition fee \$2,050

### Torion GC/MS Basic Operations

This three-day class will cover basic operations of the Torion T-9 Portable GC/MS system. Topics will include:

- Setting up the instrument for laboratory or field use
- Start up and preparation
- Conducting Performance Validation runs to ensure correct operation of the system
- Basic SPME sampling techniques for Gas and liquid samples
- Basic data viewing and analysis using Chromion Software
- Very basic maintenance for the T-9 GC/MS and SPME Custodion

N023-9011 | 5 Days | Tuition \$3,150

### Torion GC/MS Advanced Operations

This five-day class will cover advanced operations of the Torion T-9 Portable GC/MS system. Topics will include:

- Advanced data analysis using Chromion software
- Building custom target lists and unknown libraries in Chromion
- Air Sampling using the Custodion Needle Trap and Clarion air pump



## LC Mass Spectrometry

N023-3019 | 3 Days | Tuition \$2,050

### AxION 2 DSA

#### Prerequisite

ChromAcademy LC and LCMS online training.  
Understanding of basic MS theory and operation  
(1-2 years) recommended.

#### Lectures and Labs

Use the ToF Driver Software to Tune and Calibrate the ToF MS  
Use Autotune to refine Tune files and detector voltages  
Connect/configure the system with the AxION DSA to perform Direct Sample Analysis  
Build Acquisition methods for the DSA/ToF MS  
Run Samples as a DSA/ToF using the DSA controller  
View and Analyze DSA ToF data using the ToF Driver and AxION Solo

N023-3020 | 5 Days | Tuition \$3,150 (if not included w/instrument purchase)

### AxION 2 LC/DSA

#### Prerequisite

Fundamental LC training course, or a minimum of  
one year experience running an LC instrument.  
ChromAcademy LC and LCMS online training.  
Understanding of basic MS theory and operation  
(1-2 years) recommended.

#### Lectures and Labs

Use the ToF Driver Software to Tune and Calibrate the ToF MS  
Use Autotune to refine Tune files and detector voltages  
Build Acquisition methods for the LC/ToF MS and for the DSA/ToF  
Connect/configure the system with the Flexar LC system to run the system as an LC/ToF  
and with the AxION DSA to perform Direct Sample Analysis  
Run Samples as an LC/ToF using Chromera and as a DSA/ToF using the DSA controller  
View and Analyze LC/ToF and DSA ToF qualitative and quantitative data

N023-3022 | 3 Days | Tuition \$2,050

### AxION 2 LCMS

#### Prerequisite

Fundamental LC training course, or a minimum of  
one year experience running an LC instrument.  
ChromAcademy LC and LCMS online training.  
Understanding of basic MS theory and operation  
(1-2 years) recommended.

#### Lectures and Labs

Use the ToF Driver Software to Tune and Calibrate the ToF MS  
Use Autotune to refine Tune files and detector voltages  
Build Acquisition methods for the LC/ToF MS including the use of EIC data  
channels and LockMass Calibration  
Connect/configure the system with the Flexar LC system in order to run the  
system as an LC/ToF  
Run Samples as an LC/ToF using Chromera  
View and Analyze LC/ToF data using both the ToF Driver (qualitative) and  
Chromera (quantitative)

N023-3042 | 5 Days | Tuition \$3,150

### QSight® Triple Quad LC/MS/MS

#### Prerequisite

Students are required to have basic knowledge on  
mass spectrometry and chromatogram. It would be  
ideal to have LC-MS working experiences including  
LC and MS method development. The students are  
also required to have at least one month of experience  
on the installed and operating QSight LC-MS system.

#### Lectures and Labs

Overviews including the fundamentals of mass spectrometry, QSight technology,  
Simplicity software, basic applications, and basic maintenance and troubleshooting.  
Hands -on practices including instrument tuning and calibration, QSight Acquisition  
and Simplicity software familiarization, method development, sample preparation and  
data analysis, and basic maintenance and troubleshooting.



## Infrared Spectroscopy

N020-0126 | 2 Days | Tuition \$1,450

### Introduction to FT-IR and Spectrum Software

This two-day fundamental course provides instruction in theory, instrumentation, maintenance, data acquisition software, sample preparation and the use of various accessories, enabling students to acquire good quality spectra by either reflectance or transmission. The course will cover instrument maintenance, calibration, and the preparation of solids, liquids, and pastes for both reflectance (Universal Attenuated Total Reflectance UATR), and transmission studies. The use of the UATR accessory will be the main focus on day two along with additional techniques for sample preparation.

#### Lectures and Labs

IR theory and advantages of IR spectroscopy  
 Instrument parts, maintenance and validation  
 Software operation  
 Sample preparation using the UATR  
 Casting films on IR transparent windows  
 KBr pellet making  
 Gas cell preparation  
 Troubleshooting

N020-0152 | 2 Days | Tuition \$1,450

### IR Spectral Interpretation for Beginners

This two-day infrared interpretation course provides students with basic knowledge of the location of absorption bands and the chemical bonds present that are the cause of these absorptions. Various examples of infrared (IR) spectra representing a wide variety of chemical classifications both aliphatic and aromatic will be shown. Each sample spectra discussed will focus on the chemical bonds present and the corresponding infrared absorption. The student will learn the most direct way to interpret an IR spectrum and how to compare an unknown spectrum with known reference spectra. Exercises will focus on a displayed spectrum and the chemical bonds present that cause an infrared absorption in a particular region of the spectrum.

N020-0188 | 2 Days | Tuition \$1,450

### AssureID Software

This two-day workshop includes the use of AssureID software for the collection and evaluation of Mid Infrared and Near Infrared product spectra. Method Explorer, Method Developer, Analysis Setup and Results Browser modules will be demonstrated, and IQ/OQ/PQ setup and System Suitability testing will be covered. Extensive time will be devoted to hands on method development, including spectra processing, workflow choices, and final reports, as well as evaluating and manipulating the model results.

## Thermal Analysis



N020-0607 | 2 Days | Tuition \$1,450

### Introduction to Differential Scanning Calorimetry (DSC 4000, 6000 or 8000, Diamond DSC and Jade)

This course is for students who have recently started using the DSC 4000, 6000 or 8000, Diamond™ DSC or Jade instruments. The practical aspects of operation and maintenance as well as introductory theory are covered in the course. In addition, the common functions of the Pyris™ software and their application to the use of the various power compensated DSC instruments will be demonstrated.

#### Prerequisite

The student should have spent at least one month becoming familiar with the instrument.

#### Lectures and Labs

- System configuration/overview
- Setup
- Baseline optimization
- Calibration
- Maintenance and troubleshooting
- Practical laboratory experiments
- Generic applications
- Data handling
- Operating variables
- Safety

*\*The Pyris Player and the TGA Autosampler will not be covered in this course.*



N020-0605 | 2 Days | Tuition \$1,450

### Introduction to Thermogravimetric Analysis (TGA 4000, TGA7 and Pyris 1 TGA)

This course is designed for people who have recently started using TGA 4000, TGA7 and Pyris™ 1 TGA. The practical aspects of operation and maintenance as well as introductory theory are covered.

In addition, the course covers the common functions of the standard Pyris software and applications of the TGA7 and Pyris 1 TGA.

#### Prerequisite

The student should have spent at least one month becoming familiar with the instrument.

#### Lectures and Labs

- System configuration/overview
- Setup
- Calibration
- Sample handling
- Maintenance and troubleshooting
- Practical laboratory experiments
- Generic applications
- Data handling
- Operating variables
- Safety

*\*The Pyris Player and the TGA Autosampler will not be covered in this course.*

## Training Locations - Life and Analytical Sciences

### Atlanta, Georgia

PerkinElmer, Inc.  
11695 Johns Creek Parkway  
Suite 150  
Johns Creek, GA 30097

### Chicago, Illinois

PerkinElmer, Inc.  
2651 Warrenville Road  
Suite 100  
Downers Grove, IL 60515

### San Jose, California

PerkinElmer, Inc.  
75 Nicholson Lane  
San Jose, CA 95134

### Shelton, Connecticut

PerkinElmer, Inc.  
710 Bridgeport Avenue  
Shelton, CT 06484-4794  
Tel: 800-762-4000 x4  
Fax: 203-944-4902

### Toronto, Canada

PerkinElmer, Inc.  
501 Rowntree Dairy Road  
Unit 6  
Woodbridge, Ontario L4L8H1

## Customer Training/How to Enroll

Please mail or fax the completed enrollment form to:

### For Canada/US Enrollment

PerkinElmer, Inc.  
710 Bridgeport Avenue  
Shelton, CT 06484  
Tel: (800) 762-4000 option 4  
Fax: (203) 944-4902  
For more information please call (800) 762-4000 x4  
(outside US) or fax (203) 944-4902.

Please carefully review all payment terms and conditions listed below.

### Registration

Purchase order or credit card payments must accompany the registration. A confirmation email will be sent within 48 hours of receipt of enrollment. If the class you have requested has been filled, a training representative will contact you to advise you of alternate class availability and place you on a waiting list to fill any late cancellations.

### To register by phone:

Contact the Training Center (800) 762-4000, option 4 with course request and payment information.

## Course Agreement/ Refund Policy

### Cancellation

PerkinElmer may cancel or reschedule courses up to 14 days prior to the start date if the minimum enrollment is not met. A full refund of the tuition fee will be provided if this happens. If air travel is required, we discourage the purchase of non-refundable tickets.

PerkinElmer is not responsible for any travel expenses incurred due to a class cancellation.

Our refund policy regarding a student's cancellation from a class is as follows:

Up to 14 days prior to the course	Full Refund
Less than 14 days	No Refund ( <i>customer may schedule alternate date</i> )

All cancellations must be made with the training administration office directly. For all training inquiries or cancellation, please contact us at (800) 762-4000, option 4.

### Payment

Pricing is quoted in US Dollars. Payment is required upon enrollment. Purchase order and credit cards accepted. If paying by purchase order, a faxed copy of the purchase order must be provided during course enrollment.

PerkinElmer's OneSource® laboratory services leverage deep scientific knowledge and expertise in the development of the most comprehensive suite of scientific laboratory services, from instrument maintenance and enterprise solutions to scientific consulting services to optimize laboratory efficiency. Discover our integrated approach and take advantage of an ideal set of tools to help empower your science and drive your business.

For more information, please visit [www.PerkinElmer.com/OneSourceTraining](http://www.PerkinElmer.com/OneSourceTraining)

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2017-2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

013799E\_01

PKI

# CLASSROOM TRAINING COURSE ENROLLMENT FORM

Email application to: [ustraining@perkinelmer.com](mailto:ustraining@perkinelmer.com)

PerkinElmer will not guarantee seating without purchase order or credit card for payment. PH: (800) 762-4000, option 4 (US/Canada only) / FAX: (203) 944-4902

NAME \_\_\_\_\_

PHONE \_\_\_\_\_ FAX \_\_\_\_\_ EMAIL \_\_\_\_\_

COMPANY \_\_\_\_\_

ADDRESS \_\_\_\_\_

CITY/TOWN \_\_\_\_\_ STATE/COUNTY \_\_\_\_\_ ZIP/POSTAL CODE \_\_\_\_\_

EXPERIENCE WITH PERKINELMER PRODUCTS \_\_\_\_\_ # OF YEARS \_\_\_\_\_

DIETARY REQUIREMENTS \_\_\_\_\_

\*PerkinElmer reserves the right to cancel a course if the minimum requirement is not met. Two (2) weeks minimum notice will be given.


**METHOD OF PAYMENT: \*\*Payment must be submitted UPON registration.**  
Hard copy purchase orders must be submitted to 203-944-4902 (001 203-944-4902 Europe). PerkinElmer cannot guarantee seating without payment.

\_\_\_\_\_ BILL MY COMPANY USING PURCHASE ORDER NUMBER \_\_\_\_\_  
(Copy of purchase order *MUST* be included with registration in order to guarantee a seat in the class)

\_\_\_\_\_ CREDIT CARD (circle one)    **MC**    **VISA**    **AMEX**

Exp.Date: \_\_\_\_\_ Number: \_\_\_\_\_

(Credit card information **MUST** be supplied upon registration).

Name as it appears on card: \_\_\_\_\_

\_\_\_\_\_ TRAINING WAS INCLUDED WITH SALE OF INSTRUMENT

(Please provide PerkinElmer sales order number) \_\_\_\_\_

**PerkinElmer, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2017-2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.