



2024 Europe CUSTOMER TRAINING

January - December



PerkinElmer brings you the best laboratory and applications service in the industry, the quality consumables you need to optimize performance, and **stellar customer training** – when and where you need it.



CENTER OF
EXCELLENCE

Course	Course #	Virtual Training	Rodgau, Germany
TotalChrom Software – ONLINE	N0233204	April 22-23 October 28-29	Available Upon Request
SimplicityChrom CDS Software			Available Upon Request
LC Fundamentals and Operator Training	N0200535		Available Upon Request
LC-MS/MS with Simplicity Software	N0233060		Available Upon Request
GC Fundamentals and Operator Training	N0200430		March 12-14 August 27-29
GC Fundamentals Operator Training – ONLINE	N0233205	May 13-14 December 2-3	Available Upon Request
GC 2400 Platform User Training			Available Upon Request
GC Headspace Operator Training– ONLINE	N0233206	September 30	
GC Headspace Fundamentals Operator Training	N0200431		April 9-10 November 5-6
Headspace - Trap Workshop			April 11 November 7
GC-MS Fundamentals Training	N0232038		January 22 October 7
GC-MS Software Training	N0232039		January 23-24 October 8-9
GC-MS Operator Training	N0232040		January 25 October 10
GC-MS Fundamentals and Operator Training	N0200432		January 22-25 October 7-10
GC-MS Operator Training - ONLINE	N0233207	June 10-12 November 25-27	

**REGISTER FOR TRAINING
IN 3 EASY STEPS AND
SECURE YOUR SPOT!**

- 1 [Click to go to the PerkinElmer Education and Training Portal](#)
- 2 Register for required course and date
- 3 Happy training



2024 Europe

CUSTOMER TRAINING

January - December



PerkinElmer brings you the best laboratory and applications service in the industry, the quality consumables you need to optimize performance, and **stellar customer training** – when and where you need it.



CENTER OF EXCELLENCE

Course	Course #	Virtual Training	Rodgau, Germany
IR Fundamentals and Operator Training	N0200128		April 16-17 October 1-2
IR Fundamentals – ONLINE	N0233212	February 19 September 16	
IR Spectrum Operator Software Training	N0200131		June 18-19
IR - Spectrum Software Training – ONLINE	N0233213	March 11 October 21	
IR - Interpretation - Module 1-4 – ONLINE	N0233214	May 13-14	
IR - Interpretation - Module 5 – ONLINE	N0233215	May 15	
IR Spectral Interpretation	N0200129		November 4-6
FTIR - Microscopy Spotlight Operator Training	N0237031		Available Upon Request
DSC Fundamentals and Operator Training	N0200632		June 24-26
DSC Fundamentals and Software – ONLINE	N0233218	November 11-12	
UV WinLab Fundamentals Software Training	N0200211		July 1-3
UV WinLab Advanced Software Training	N0235003		July 4
UV WinLab Software – ONLINE	N0233219	March 4-5 December 2-4	
Quality Management UV WinLab - ONLINE	N0233220	March 6 December 5	

**REGISTER FOR TRAINING
IN 3 EASY STEPS AND
SECURE YOUR SPOT!**

- 1 [Click to go to the PerkinElmer Education and Training Portal](#)
- 2 Register for required course and date
- 3 Happy training



2024 Europe

CUSTOMER TRAINING

January - December



PerkinElmer brings you the best laboratory and applications service in the industry, the quality consumables you need to optimize performance, and **stellar customer training** – when and where you need it.



**REGISTER FOR TRAINING
IN 3 EASY STEPS AND
SECURE YOUR SPOT!**

- 1 [Click to go to the PerkinElmer Education and Training Portal](#)
- 2 Register for required course and date
- 3 Happy training

			CENTER OF EXCELLENCE
Course	Course #	Virtual Training	Rodgau, Germany
AA Flame Familiarization Training	N0200920		April 9-11 September 24-26
AA Operator Training Flame	N0231008		February 6-7
AA Operator Training Graphite Furnace	N0231009		October 22-23
THGA Method Development	N0200924		May 14-15
AA Fundamentals and Software - Flame – ONLINE	N0233200	March 25 September 30	
AA Fundamentals and Software – Graphite Furnace – ONLINE	N0233201	March 26 October 1	
WinLab / Syngistix Software for AA – ONLINE	N0200925	September 9	
ICP-OES Fundamentals and Operator Training	N0200078		February 19-22 September 16-19
ICP-OES – Fundamentals and Software – ONLINE	N0233208	March 4-5 November 18-19	
ICP-OES Advanced Operator Training	N0200079		April 3-4 November 26-27
WinLab / Syngistix Software for ICP – ONLINE	N0200035	April 15	
ICP-MS Fundamentals and Operator Training	N0200194		Jan 29-Feb 2 March 18-22 June 3-7 November 11-15
ICP-MS Advanced Operator Training	N0236005		February 27-28 July 2-3 October 15-16
ICP-MS Fundamentals and Software – ONLINE	N0233210	March 11-13 October 7-9	
Training – General Lab Practices - ONLINE	N0209777	April 29	



2024

CUSTOMER TRAINING

Course Descriptions and Objectives



PerkinElmer brings you the best laboratory and applications service in the industry, the quality consumables you need to optimize performance, and **stellar customer training** – when and where you need it.



Atomic Absorption

Atomic Spectroscopy courses provide the analyst with the knowledge and skills needed for operating, optimizing and troubleshooting including:

- Technical introduction/features
- Setting up element parameter file
- Background correction
- Optimization and calibration
- Flame emissions analysis
- Interferences and Matrix Modifiers
- Maintenance and Quality Control

Inductively Coupled Plasma Optical Emission Spectrometry

ICP-OES courses provide analysts with emission theory and the knowledge to develop a method for analysis of complex samples. Courses include the following topics:

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

ICP-MS Inductively Coupled Plasma Mass Spectrometry

Courses will cover the NexION series of ICP-MS along with the following topics:

- Theory and method development
- Hardware overview
- NexION software training
- Optimization and instrument setup
- Universal Cell Technology (UCT)
- Data acquisition, visualization, reporting
- Troubleshooting and maintenance
- Reporting and data exporting
- Interference and correction tools

Gas Chromatography

GC courses provide an overview of the basic principles, hardware, and operational techniques including:

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

GC/MS

GC/MS courses provide users with orientation to the hardware, software, and experimental requirements including:

- System maintenance and software troubleshooting procedures
- Connecting the GC and MS
- Spectral data processing and cold starting the instrument
- Library searching, Instrument Tuning
- Method for construction and data analysis
- Instrument control, data acquisition
- MS method for quantization and GC control method

Liquid Chromatography

LC courses provide an overview of the principles and practice of high-performance liquid chromatography including:

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

LC/MS

LC/MS courses cover an overview of the fundamentals of mass spectrometry along with the following:

- QSight technology
- SimplicityChrom™ software
- Basic applications, maintenance, technology
- Creating LC/MS/MS methods
- Mass calibration
- Setting up batches
- Acquiring data, data analysis
- Demonstration data using SimplicityChrom™ software

CDS Chromatography Data Systems

CDS courses provide users with a fundamental knowledge of the theory and operation including:

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

Molecular Spectroscopy

Molecular Spectroscopy courses provide instruction in theory, instrumentation, and maintenance for both reflectance and transmission studies along with the following:

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



2024

CUSTOMER TRAINING Course Descriptions and Objectives



PerkinElmer brings you the best laboratory and applications service in the industry, the quality consumables you need to optimize performance, and **stellar customer training** – when and where you need it.



PerkinElmer[®]

Thermal/Elemental Analysis

Thermal/Elemental Analysis courses are designed to help users quickly become proficient on their own with DSC systems including:

- Basics of thermal analytics
- DSC principles and troubleshooting
- Influence on the measurement data
- Calibration, method editor, data acquisition, multi-tasking, and data optimization and evaluation
- Pyris software and Pyris Player
- Measurement and interpretation of data, reporting

UV-VIS Spectroscopy

UV-VIS Spectroscopy courses provide analysts with the ability to configure and use the software, create methods, create the equations, (process), and realize the report formats including:

- Understanding the structure of the software
- Different methods (wavelength scanning, wavelength programming, quantitative analysis with calibration curves and kinetic method)
- Method creation and report design
- Database maintenance and queries