

RUN YOUR INSTRUMENT AT PEAK EFFICIENCY

OneSource
Laboratory Services

Basic and Advanced Training for Inductively Coupled Plasma Optical Emission Spectroscopy (ICP-OES)

Basic Onsite Training for Operational Excellence

Learn tips and techniques to keep your ICP instrument up and running at peak performance. Receive expert, hands-on instruction from a certified training service engineer right in your lab. Courses range from one to three days, and you can choose from the following modules:

Theory and Hardware/Software Overview

Covers ICP-OES theory and chemistry, including how to identify optimal plasma emission zones (ionic/atomic emission) for analysis. Review ICP-OES hardware and sample introduction options, as well as software options and functionality.

Setup and Sample Analysis Review

Learn how to perform basic method setup with manual and automated analysis options. Covers instrument hardware checks plus optimization and performance checks. Review the importance of background equivalent concentration (BEC) and how to calculate it. Gain an understanding of detection limits. Also covers data viewing and reporting features.



Avio® 200 ICP-OES



Avio® 500 ICP-OES

Method Editor

Detailed review of how to create and edit methods. Includes how to set up analytes, how to set up sample acquisitions, and how to optimize spectrometer acquisition. Learn how to set up QC checks (periodic, spike recovery, etc.), plus review options for result output. Covers data process functions and options.

Interferences

Learn about types of interferences and typical correction techniques. Covers background correction for basic spectral interference, plus Multicomponent Spectral Fitting (MSF) and Inter-element Correction (IEC) for advanced spectral interference correction.

Maintenance and Troubleshooting

Reviews the maintenance required to ensure consistent instrument performance plus troubleshooting and problem-solving techniques.

Data Maintenance

Includes Syngistix™ file structures, how to manage/maintain data and method databases, and time-dependent QC charting. Learn how to create a custom report and options for data export.

Advanced Optimization Training and Applications

Looking to solve your lab's specific challenges or focus on accelerated topics such as method optimization? Contact us to discuss our customized advanced training packages.

For more information:

Email: OnSiteTraining@PerkinElmer.com | Visit: 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



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

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