

EDUCATION SERVICES TO HELP ACCELERATE YOUR SCIENCE

OneSource
Laboratory Services

Invest in Instrument Training

You are under tremendous pressure to meet tight budget constraints while performing more experiments than ever. Due to these dynamics, instrument training can seem like a luxury versus a necessity, and important benefits can be overlooked.

PerkinElmer Education Services enables you to:

- Keep your instrument running at peak performance with less downtime.
- Discover new instrument capabilities.
- Advance your knowledge base and skill set.
- Overcome the unique challenges of your application.

Choose the Option That's Right for You:

Classroom Training

Conducted by faculty at PerkinElmer Centers of Excellence

- Courses ranging from the basics to advanced techniques
- Hands-on experience in a laboratory environment
- Peer networking and collaboration
- Small class size to ensure individual attention
- 1-5 day options

Basic Training: Operation Essentials

Conducted by a certified Service Engineer in your lab

- For the beginner to intermediate user
- Courses adapted to meet your lab's needs
- System overview and basic principles
- Hands-on training using your instruments
 - Operational techniques
 - Instrument optimization/maintenance
 - Software operations/data acquisition
- 1-3 day options; up to four users per session

Advanced: Optimization Training and Applications

Conducted by a specialized Field Application Scientist in your lab

- For the experienced user
- Courses designed to address your unique challenges
- Customized hands-on training
 - In-depth instrument and software operations
 - Method optimization
 - Application-specific, consultative training
- 1-5 day options; up to four users per session

For more information including custom training options:

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


PerkinElmer
For the Better

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.

013531A_01

PKI