



GET THE MOST FROM YOUR PERKINELMER INSTRUMENTS

Training options for every stage of the instrument lifecycle

From new user courses in a traditional classroom setting to a customized training session with our experienced scientists in the comfort of your lab -- PerkinElmer has training options to help you gain the knowledge and insights you need to unlock everything your instruments offer.



For More Information on the Training Course

LEARN MORE



The Introduction You Need to Get Started. Provided with each instrument installed, Service Engineers deliver *Familiarization Training*.



Learn at Any Time, from just about Anywhere! Online training courses offer flexibility to learn at your own pace to help increase productivity and instrument performance.



Virtual Instructor Led Training (VILT) emulates the classroom experience right from your computer with small class sizes, group exercises, and discussions.



Get A Hands-on Experience in a Dedicated Laboratory Environment. Check out our *Classroom Training* schedule held at one of our PerkinElmer Centers of Excellence.



Basic Training is provided at your location for small groups and delivered by our certified Service Engineers. We focus on software operation, instrument maintenance, optimization guidance and more!



Advanced Training offers consultative training solutions for method optimization with PerkinElmer's Application Specialists.

The PerkinElmer Training Experience



Familiarization Training

- ✓ Training occurs during the instrument installation
- ✓ How to operate the main functions of the instrument
- ✓ Learn how to set up a simple method and perform quantitative analysis
- ✓ How to run the instrument software



Online Training

- ✓ Learn anytime from anywhere, at your own pace
- ✓ Save travel time and costs
- ✓ Individual modules and discounted learning bundles available
- ✓ Revisit courses and learn one topic at a time
- ✓ Online courses available for applications, tips and tricks, and more
- ✓ Gain valuable maintenance and operational training to limit downtime



Virtual Training

- ✓ 1-3 days of Virtual Instructor Led Training (VILT)
- ✓ Learn from just about anywhere
- ✓ Access to live instructors providing consultative training with results analysis and online Q&A
- ✓ Group exercises and discussions conducted using simulation software to emulate laboratory setting



Classroom Training

- ✓ 1-5 days of training conducted at PerkinElmer Centers of Excellence by subject matter experts
- ✓ Courses ranging from the basics to advanced applications and methods
- ✓ Hands-on experience in a laboratory environment
- ✓ Peer networking and collaboration
- ✓ Small class size to ensure individual attention
- ✓ Trainees can practice with instruments without affecting their lab setups



Basic Training

- ✓ Training sessions delivered at your location for up to 4 users, over 1-2 days
- ✓ System overview and basic principles
- ✓ Operational techniques
- ✓ Instrument optimization and maintenance
- ✓ Software operation and data acquisition



Advanced Training

- ✓ 1-5 days of training; up to four users per session
- ✓ Training conducted by specialized application scientists at your location
- ✓ Customized hands-on training
- ✓ In-depth instrument and software operations
- ✓ Method optimization
- ✓ Application-specific, consultative training with results analysis

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 ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.