

Introduction to application training courses

ThermoFisher
SCIENTIFIC

Increase your working knowledge of applications and instrument operation, including hands-on lab experience, with Application Training Courses, available at our Customer Training Centres located throughout Europe.

Our centres are state-of-the-art installations for training and collaborative research. We offer interactive instructor led training, a fully staffed working laboratory and the unique opportunity to collaborate with our application scientists.

Spend some time in our lab to get the most out of the time you spend in yours

We offer hands-on and practical training for a wide range of applications and all experience levels:

- **Basic introductions** to sequence analysis, real-time PCR, flow cytometry, and cell biology applications
- **Advanced applications** such as forensics, pathogen detection, and whole-genome analysis

Our classes are designed to help you get the most out of your experiments.

Courses are delivered across the region, with our main sites in the UK, Germany, France and Dubai offering the full range of classes. The most popular courses are also available in Spain, Italy, Sweden and South Africa.

Benefits of attending one of our courses:

- **Learn complete workflows**
- **Access the latest instrumentation and applications**
- **Interact with other researchers** working on similar projects

Courses are delivered by our award-winning* team of application scientists, who have extensive collective experience using, training on, and troubleshooting the entire Life Technologies™ portfolio.

Their expertise, plus our world-class facilities, make us the ideal choice for all of your training needs.

For more information on courses available, and to register, go to

learn.lifetechnologies.com/europe

Email the training team at

learn-europe@lifetech.com



- Customer Experience Centres
Glasgow, Darmstadt, Dubai and Paris
- Customer Training Labs
Madrid, Milan, Stockholm
- Customer Demo Labs
Johannesburg

*Best Customer Service & Technical Support by The 2011 Life Science Industry Awards™

Popular application training courses include:

- **Ion PGM™ & Ion Proton™ system workflow (3 days)**
 - Introduction to library prep
 - System workflow
 - Template preparation/setup
 - Enrichment
 - Initialisation of the instrument
 - Introduction to the Torrent Server
 - Sample loading and starting a run
 - Validation and review of the data
 - Introduction to Ion Reporter™ Software
 - Application-related data analysis
 - Instrument maintenance
- **Ion DNA Library preparation & essentials (1 day)**
 - Introduction to system workflow
 - Utilising Ion Xpress™ Plus and/or Library Builder™ systems
 - Library construction
 - Library qualification
- **RNA Library construction (3 days)**
 - Working with RNA
 - Small RNA with poly(A) enrichment system workflow
 - Whole-transcriptome system workflow
 - Evaluation of generated libraries
- **Real-time PCR (1 day / 2 days)**
 - Real-time applications:
 - Real-time (qPCR) instrument training
 - Gene expression workflow and data analysis
 - TaqMan™ genotyping workflow and data analysis
 - Introduction to pathogen detection
 - High resolution melting (HRM) training
 - Protein thermal shift training
 - Introductory CE-IVD real-time PCR training
 - Introductory QuantStudio™ with OpenArray™ system training
 - Gene silencing using RNAi
- **Digital PCR training (1 day / 2 days)**
 - Principle of digital PCR
 - Sample preparation
 - Data analysis using DigitalSuite™ software
- **CE – Capillary Electrophoresis (1 day / 2 days / 3 days)**
 - CE instrument training course
 - Sequencing applications and software training course
 - Fragment analysis applications and software training course
 - Introductory CE-IVD CE sequencing training
- **Molecular & cell biology (1 day / 2 days / 3 days)**
 - Basic molecular biology
 - Basic cell culture
 - Gene silencing using siRNA
 - Practical cloning course in Gateway™ technology
 - Attune™ Acoustic Focusing Cytometer and Autosampler training
 - Pluripotent stem cell workshop
- **Bioinformatics education for life scientists (1 day)**
 - Relational databases – MySQL
 - Introduction to Perl or Python
- **Bioinformatics for molecular biologists (4 days)**
 - Introduction to programming
 - Relational databases
 - Practical project (creating a knowledge base from NGS data)

Individual, customised training and consultation

While learning in a group is often the most productive and effective method, sometimes this simply isn't practical.

With that in mind we can arrange to deliver all of these courses, or parts of these courses, as well as consultancy services in your facility.

Find out more at lifetechnologies.com/europe

ThermoFisher
SCIENTIFIC