

Sartorius Lab Academy Customized Training for Laboratory Professionals



Increase the Quality, Efficiency and Safety of Your Workflows

Professional and customized training is the basis for successful, reliable and efficient work, and is instrumental in boosting your motivation and job satisfaction. Systematic training will enable you to successfully meet the challenges posed by rapid evolving technologies, complex regulatory requirements and the need to make constant improvements in productivity.

This is why Sartorius offers a comprehensive training program, which focuses on making the steps in your daily lab work more efficient and even more productive while increasing quality awareness. Our courses and seminars provide a high percentage of hands-on training to extend your techniques and to give you valuable advice and tips from the experts. In this brochure, we will introduce our program of concise training sessions, called Lab Academy.



Sartorius Lab Academy – Why It's Worth Attending

This is what you can expect:

- Intensive professional training that combines theory and practice
- Theoretical part for increasing background knowledge and gaining industry updates
- Hands-on session to put theory into practice and rehearse work situations
- Extensive course material provided for later reference

Your advantages:

- Individual knowledge customtailored to your requirements and routine procedures
- Industry updates about the most recent changes to the relevant standards and regulations and about trends
- Solutions for common problems based on actual work situations and challenges
- First-hand tips from knowledgeable and experienced experts to facilitate your daily work

Program

Sartorius | filtration academy



Sartorius | wateracademy



Sartorius | pipetting academy



Sartorius | weighing academy



Target Group

Laboratory professionals who would like to enhance the efficiency of their workflows and receive valuable tips on maximizing the benefits of using Sartorius technologies.

Trainers

You will be trained by Sartorius industry experts. They are professionals with extensive experience and are specialists in their fields.

Location

We will conduct training sessions either at your site, at another local facility of your choice or at one of the many Sartorius Training Centers located around the world.

Certificate

After the training, each participant receives a certificate from the Sartorius Lab Academy.

Register here:



www.sartorius.com/lab-academy

Sartorius | filtration academy

Filtration is an initial and essential task in a large number of applications in the lab. Getting this step right by choosing the adequate tool is essential for achieving reliable and reproducible results. The Sartorius filtration academy will give you the background knowledge and practical insight to choose the right filtration devices, use them as ergonomically as possible and speed up your work.

Modules

Cleanliness Analysis:

- Overview on Technical Cleanliness and Cleanliness Analysis
- Information on how to achieve the most accurate results in the filtration step by selecting the right filter
- Guidance on which balance attributes are essential for the gravimetric analysis step and selecting the right model
- Basic trouble shooting in filter weighing

Filtration of Small Volumes:

- Practical selection guide for choosing the correct membrane and device types for different applications
- Key factors to consider in terms of device quality, integrity and chemical compatibility
- Essential things to consider for the filtration of samples prior to analytics like HPLC and ICP/MS
- Integrity testing on device after sterile filtration
- Ergonomic factors for working with syringe and syringeless filters

Protein and Virus Concentration:

- Introduction to ultrafiltration terminology and definitions
- Achieving highest sample recovery after protein concentration with the suitable device and the correct MWCO
- Introducing crossflow as a versatile concentration method in labs
- Buffer exchange and diafiltration using ultrafiltration



Sartorius | pipetting academy

Technique is key for pipetting. With the right technique, you get the most out of your precious samples and make the best use of your valuable time while pipetting without any pain. The Sartorius pipetting academy will give you the background knowledge and practical insights to work efficiently and safely, choose the right material and ensure efficient and compliant calibration – for best results even in long pipetting series.

Modules

Pipetting Techniques:

- Understanding the influence of pipetting techniques on pipetting results
- Identification and effect of different error sources on pipetting results
- How to select optimum pipetting tools and techniques for different liquids and conditions
- How to achieve the best possible pipetting results

Ergonomics:

- Recognition and minimization of ergonomic risk factors in pipetting and lab work
- Selecting the right tools and working postures
- Application of best ergonomics to make your work more enjoyable

Pipetting in Cell Culture Applications:

- How to avoid contamination
- Best practices for cell culture work
- Recognition of the factors affecting your results

Pipetting Standards and Maintenance:

- How to work according to ISO standards
- Handling of pipettes as precision instruments
- Ensuring the working condition of pipettes



Sartorius | wateracademy

Impurities are found in all laboratory water and have a negative influence on your scientific analyses. With the Sartorius water academy, you learn where impurities come from and how to purify water to the standard you need for your specific application. In order to save resources, you'll gain a theoretical and practical understanding of how to work most efficiently with your lab water system.

Modules

Water Impurities:

- Types of impurities in drinking water
- Effect of impurities on water purification systems and measurements
- Testing and removing impurities (purification technologies)

Water Quality:

- Classification of water
- Definition of water qualities
- Regulatory affairs and standards

Lab Water Systems:

- Identifying the ideal lab water solution for a given application
- Best practices for efficiency, correct quality and quality consistency



Sartorius | weighing academy

Especially in pharmaceutical laboratories, compliance with international regulations and quality standards is mandatory. The Sartorius weighing academy provides you with comprehensive theory and practical tips to comply with regulatory requirements. Furthermore, we make sure that you know how to avoid typical mistakes and achieve the most accurate weighing results.

Modules

General Weighing Facts:

- Terminology
- Influence of operators and environmental factors on weighing results
- Handling samples and minimizing the influence of static electricity and hygroscopic samples

Balance Calibration and Qualification:

- Calibration and servicing of your laboratory balances
- Calibration weights
- Balance qualification
- Pass | fail parameters to monitor the performance of a balance

Quality Management and Regulations:

- Complying with regulations and quality requirements in the pharmaceutical laboratory (GLP, ISO17025, USP, 21CFR Part 11 etc.)
- Operating range according to USP regulations



Do You Already Know Our Intensive Sartorius Training?

You agree to our Lab Academy Program and perhaps you already participated? Then you will surely be interested in our detailed intensive training courses. Each unit lasts 1–3 days and is held in Goettingen, Germany; Shanghai, China; Bangalore, India; Bohemia, United States and Singapore. These training courses focus on sharing information with other colleagues in your field, as well as on interactive group work closely related to your applications.

For more information, visit: www.sartorius.com/training



Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen, Germany

Phone +49.551.308.0 Fax +49.551.308.3289

