



Sartorius Lab Academy

Customized Training for
Laboratory Professionals

Simplifying Progress

SARTORIUS

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 filtration academy

- Basic Filtration
- Filtration Devices
- Ultrafiltration Devices

Sartorius pipetting academy

- Pipetting Techniques
- Ergonomics
- Pipette Calibration

Sartorius water academy

- Water Impurities
- Water Quality
- Lab Water Systems

Sartorius weighing academy

- General Weighing Facts
- Balance Calibration and Qualification
- Quality Management and Regulations

Register here:

<https://www.sartorius.com/en/knowledge/trainings>



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



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.

Costs and Registration

The price for attending the Sartorius Lab Academy is just as individual as your training. Explain to your local Sartorius partner which content you would like to include in your individual training plan and where the training should take place. We will then send you a non-binding offer.

Our Method

- Three hours of intensive professional theoretical and practical training
- 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

Individual Training

We want you to gain maximum benefit from your training. The Sartorius Lab Academy has a modular structure so you can customize your training courses in a way that best suits your needs:

- Select content from one or more Sartorius Lab Academy programs.
- Combine individual modules or even individual topics
- Decide whether you prefer theoretical or practical training or a combination of both

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.

Your Advantages:

- You will learn to choose the right filtration medium for your application from the comprehensive range available on the market
- You will learn about ergonomic handling and find out how to accelerate your processes and achieve maximum sample recovery
- You will learn how to choose the right filtration system as well as the best ultrafiltration unit and method for concentrating your sample

Modules

Basic Filtration:

- Overview of basic filtration media (paper and membrane filters, glass and quartz microfilters)
- Basic knowledge of membrane structures, materials and pore sizes
- Choosing the right material for various applications
- FAQs (for example: How can I compare paper types from different suppliers with one another?)

Filtration Devices:

- Practical guidelines for choosing the right membrane and unit for sample preparation
- Key factors relating to the quality and integrity of the unit and chemical compatibility
- Integrity tests on a unit after filtration
- Ergonomic factors when working with syringe filters

Ultrafiltration Devices:

- Introduction to ultrafiltration methods
- Efficient concentration of initial sample volumes between 0.5 l and 5 l with crossflow units
- Maximum sample recovery after protein concentration using the right unit and correct MWCO
- Typical applications involving ultrafiltration units





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.

Your Advantages:

- You will discover the capabilities of your pipette as well as when to use each individual pipetting technique
- You will learn how to calibrate your pipette correctly and comply with international quality standards at work
- You can teach your colleagues the most ergonomic posture to adopt when pipetting

Modules

Pipetting Techniques:

- Influence of pipetting technique and environmental factors on your results
- Best practice pipetting methods that reduce costs and increase efficiency
- The right pipetting techniques for different liquids
- Accurate, reproducible pipetting results – irrespective of the user

Ergonomics:

- Ergonomic risks of handling liquid and how to avoid them
- Ideal laboratory environment and posture for efficient, safe, and ergonomic pipetting

Pipette Calibration:

- Correct calibration for precision results
- International quality standards (ISO 8655)



Sartorius Water Academy

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.

Your Advantages:

- You will learn about different types of water contamination, how to test them and which filtration type is most suitable for subsequent treatment
- You will learn everything there is to know about current laboratory standards and the best water quality for your scientific applications
- You will receive tips on how to reduce operating costs without compromising on water quality

Modules

Water Impurities:

- Types of impurities in drinking water
- Effect of impurities on water purification systems and measurements
- Testing and removing impurities (purification technologies)
- Minimizing the risk of recontamination and bacterial growth

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.

Your Advantages:

- You will learn how to avoid typical weighing errors and achieve precision results
- You will discover how to safely adhere to international regulations
- You will learn how to adjust and qualify your laboratory balances



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



 Find out more
For more information, please visit
www.sartorius.com

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further contacts, visit
www.sartorius.com