



Pipetting Academy

ISO Certified Trainings to
Enhance the Quality
of Your Work

Simplifying Progress

SARTORIUS

Pipetting Academy

ISO Certified Trainings to Enhance the Quality of Your Work

Reach mastery over pipetting and liquid handling through high quality training from certified Pipetting Academy trainers. The ISO 9001 certified Pipetting Academy Program guarantees that the laboratory personnel will be trained with the pipetting knowledge and skills necessary to meet quality requirements. Each participant will receive a certificate after successfully completing a training. Thousands of laboratory professionals around the globe have been trained in the Pipetting Academy since 1994.



Pipetting Academy Modules

Basics of Pipetting

It is an excellent package for anyone who is new to pipetting or feels a need to polish pipetting practices. In the Basics of Pipetting module, you will learn:

- to choose the right equipment
- the basics of pipetting techniques
- the best practices in lab ergonomics
- daily maintenance and checking of pipettes

Pipetting in Cell Culture Applications

Working in a cell culture lab has special requirements in terms of techniques and equipment.

In the Cell Culture module, you will learn:

- how to avoid contamination
- the best pipetting practices for cell culture work
- to recognize the factors affecting your results

Best Practices for Cell-Based Sample Preparation

Have you ever wondered if sample preparation could be the source of variance in your cell-based assays?

In this module, you will learn:

- The aspects that affect your cell-based assays with Incucyte or IQue screener
- How to reduce variability between experiments
- How to speed up the workflow and improve your cell health with smart pipetting practices

Best Practices for Working With Proteins

To achieve the best results, protein samples should have a high concentration and maintain a high biological activity.

Working with proteins teaches you:

- Best pipetting practices for protein work
- How to ensure your sample quality and yield

Contact Sartorius to Set Up Your Pipetting Academy

- Basic and advanced modules covering various aspects of pipetting
- Emphasis on hands-on training
- Learn mastery of pipetting practices

www.sartorius.com/pipetting-academy



Best Pipetting Practices for Nucleic Acid Work

DNA and RNA extraction and qPCR assays are the cornerstones of a large part of research in life science. In Best Pipetting Practices for Nucleic Acid Work, you will learn how to:

- Maintain sample quality and purity
- Avoid contamination
- Recognize the factors affecting your results

Ergonomics

The ergonomics module is intended for laboratory personnel and occupational and health professionals who are interested in improving the work ergonomics, minimizing ergonomic risk factors and preventing musculoskeletal disorders.

In the Ergonomics Module you will learn:

- to recognize the ergonomic risk factors in pipetting and lab work
- to select the right tools and working postures
- to apply ergonomics and make your work more enjoyable

Pipetting Standards and Maintenance

Regular pipette maintenance and calibration ensures reliable results and conformance to regulation. In the Pipetting Standards and Maintenance module, you will learn:

- to define your pipette maintenance program according to ISO 8655 standards
- to handle pipettes as precision instruments
- to ensure the reliable performance of your pipettes

Pipetting Techniques

Pipetting is a precision task influenced by multiple factors. In the Pipetting Techniques module, you will learn:

- to understand the influence of pipetting techniques on pipetting results
- to identify the influence of different error sources on pipetting results
- how to select pipetting tools and techniques for different liquids and conditions

Sales and Service Contacts

For further contacts, visit
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