



sartorius

Sartorius | pipetting academy

ISO Certified Trainings to Enhance
the Quality of Your Work



- Basic and Advanced Modules covering various aspects of Pipetting
- Emphasis on hands-on training
- Useful Tips & Tricks

Contact Sartorius to set up your
Pipetting Academy:

www.sartorius.com/pipetting-academy

turning science into solutions

Sartorius | pipettingacademy offers insight and training for pipetting related topics. It's a comprehensive combination of theory and practice that provides you with useful knowledge and tips for your daily work. Thousands of laboratory professionals around the globe have been trained in the Pipetting Academy over the past decades.

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 basic 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

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

Contact Sartorius to set up your Pipetting Academy:

www.sartorius.com/pipetting-academy

ISO Certified Training Program

Sartorius | pipettingacademy is an ISO 9001 certified training program. You will get high quality training from the certified Pipetting Academy trainers. The ISO certified Pipetting Academy guarantees that the laboratory personnel will be trained with the pipetting knowledge and skills necessary to master their tasks and meet the quality requirements. Each participant will receive a certificate after successfully completing the training.

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen, Germany
Phone +49.551.308.0
www.sartorius.com

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG.
Printed in the EU on paper bleached without chlorine.
Publication No.: LLH1512-e180501
Order No.: 98649-024-07
Ver. 05 | 2018