



sartorius stedim
biotech

Training Courses 2019 Southeast Asia



turning science into solutions

Our application-centered training courses are designed for Sartorius customers who work in research and development, production and quality assurance.

The theoretical part will help you better understand the background on your application and Sartorius products. In the practical part, you will work intensively in small groups so you can apply directly what you have just learned and quickly assimilate procedural methods.

Our approach is solution- and process-oriented to accommodate your routine work. Your procedural steps and processes will be simplified; you will gain confidence in performing procedures while increasing job satisfaction. One nice side effect: you'll get new contacts and can share experiences with other colleagues in your field.

"The only person who is educated is the one who has learned how to learn ... and change."

Carl Rogers



Welcome to our training program 2019.

You can choose from a broad range of training courses and seminars in the following areas:

Fermentation | Cell Culture Technologies page 6

Downstream Processing page 7

Single-Use Process Systems page 8

Filtration page 10

Software page 11

Our training courses focus on methodology. They are open to all interested parties and give you objective and unbiased perspectives.

It is our hope that the time you spend with us will be an excellent experience for you. We have put together these training courses in order to better enable you to meet your training objectives.



Sartorius

Overview of Our Training Courses

Fermentation | Cell Culture Technologies

Basics of Fermentation 6

Downstream Processing

B1 Use of Crossflow Filtration in DSP (Module 1) 7

B2 Use of Crossflow Filtration in DSP (Module 2) 8

Single-Use Process System

C1 Single-Use Technologies 9

Filtration

D1 Basics of Filters and Integrity Test 10



Software

E1 Multivariate Data Analysis (MVDA) 11

E2 Design of Experiments (DOE) 12

Lab Academy 13

Trainers

Inhouse Training

Your Training Facilities

About Us

Registration

Information on How to Register



Fermentation | Cell Culture Technologies

Basics of Fermentation

Who should attend: Technicians and scientists without any or basic practical knowledge or cultivation of microorganisms or cell culture

Course Date | Venue:
21 - 22 March 2019
Singapore Polytechnic

This course gives you an overview about the fundamentals of fermentation including typical applications, organisms, fermentation techniques and process control. The theoretical and practical knowledge will help you better understand your own fermentation processes and to apply fermentation knowledge for the operation of bioreactor equipment

Course Duration:
2 Days

Start First Day: 10:00am
End Last Day: 5:00pm

Trainer:
Application | Training Specialist

This course covers:

- Chemical and biological principles of fermentation
- Organisms and expression systems used for production
- Basics principles of bioreactor technologies:
 - what is a bioreactor, design of bioreactors and operations
- Different fermentation modes (batch, fed-batch, continuous fermentation)
- Basic principles of process control: sensors and automation
- Performing experiments (model solution) on bioreactor

Course Fee:
Please contact your local representative



Downstream Processing

Use of Crossflow Filtration in DSP (Module 1)

Who should attend: Technicians and scientists with or without previous knowledge in the field of crossflow filtration.

This course is designed to provide you with knowledge about the use of crossflow filtration in the biopharmaceutical industry. It consists of two different modules, which can be booked separately.

Module 1 focuses on the fundamentals of crossflow filtration and to apply crossflow filtration knowledge for concentration and diafiltration.

Theoretical aspects include:

- Crossflow filtration theory
- Membrane characterization | membrane selection
- Factors influencing performance
- Operating conditions
- Cleaning-in-place (CIP)
- Applications in biotechnology

Practical exercises:

- Operational set-up of the systems
- Determination of clean water flux
- Cell retention by microfiltration (model solution)
- Concentration of a protein solution by ultrafiltration
- Removal of low-molecular weight contaminants by diafiltration
- Cleaning

Course Date | Venue:

19 – 20 June 2019

Singapore Polytechnic

Course Duration:

2 Days

Start First Day: 10:00am

End Last Day: 5:00pm

Trainer:

Application | Training Specialist

Course Fee:

Please contact your local representative

B1

Downstream Processing



Downstream Processing

Use of Crossflow Filtration in DSP (Module 2)

Who should attend: Technicians and scientists with basic knowledge in the field of crossflow filtration.

Module 2 focuses on aspects of integrity testing of crossflow cassettes and efficient product recovery.

Theoretical aspects include:

- Scale-up
- Integrity testing
- Optimal recovery process: Harvesting of product and draining of the system after filtration
- Multi-use vs. single-use
- Use of cassettes and holders of different suppliers

Practical exercises:

- Integrity testing
- Efficient harvesting and draining of system

Course Date | Venue:

21 June 2019

Singapore Polytechnic

Course Duration:

1 Day

Start First Day: 10:00am

End Last Day: 5:00pm

Trainer:

Application | Training Specialist

Course Fee:

Please contact your local representative



Single-Use Process Systems

Single-Use Technologies

Who should attend: Technical staff with limited experience with single-use systems and consumables..

Establishing a new production process requires high initial cost and can take years to set up. At each step, from biosynthesis to formulation, materials must be validated to eliminate contamination risks.

Single-use technologies eliminate cleaning validation steps, reduce significant time and cost, provide higher flexibility and reliability.

Theoretical aspects and practical exercises:

- Bag assemblies: bag types and material, design options and handling (bag unpacking & insertion into tanks)
- Mixing technology and efficiency: magnetically coupled vs. levitated impeller
- Sensor technology (types, location, calibration, accuracy)
- Integrity testing of filters and bag leakage testing in single use systems
- Aseptic connections and disconnections of tubing
- Aseptic transfer of product across cleanroom barriers

Course Date | Venue:

19 - 20 September 2019

Singapore Polytechnic

Course Duration:

2 Days

Start First Day: 10:00am

End Last Day: 5:00pm

Trainer:

Application | Training Specialist

Course Fee:

Please contact your local representative

C1

Single-Use Process Systems



Filtration

Basics of Filters and Integrity Test

Who should attend: Technicians and scientists with little or no previous knowledge in the field of filtration.

Course Date | Venue:
04 - 05 December 2019
Singapore Polytechnic

In this training course, you will acquire the necessary knowledge and practical skills in handling the filters employed for sterilizing filtration.

Course Duration:
2 Days

Theoretical aspects include:

- Basic principles of filtration
 - Depth filters | membrane filters
 - Hydrophobic | hydrophilic filters
 - Retention mechanisms
- Integrity testing of membrane filters
 - Regulatory requirements
 - Integrity test methods
- Physical principles of the steam sterilization of filter lines

Start First Day: 10:00am
End Last Day: 5:00pm

Trainer:

Application | Training Specialist

Course Fee:

Please contact your local representative

Practical exercises:

- Manual determination of bubble point | diffusion
- Automated integrity testing
 - Bubble point test | diffusion test
 - Integrity testing of hydrophobic filters using the water intrusion test (WIT)
- Trouble shooting



Software

Multivariate Data Analysis

Who should attend: Researchers, scientists and engineers working with collection and interpretation of data.

Using the latest multivariate techniques, participants will learn how to interpret complex data quickly and confidently. Discover the secrets of overviewing data tables and also learn how to build robust predictive models that turn data into decisions. The course is composed of lectures, software demonstrations and computer exercises in software SIMCA, based on real-life datasets.

Course Objectives: To show how to convert data tables to robust and reliable models and figures that can be easily interpreted and overviewed.

Topics include:

- Organization, visualization and treatment of different types of data
- Multivariate data modeling and diagnostics
- Understanding PCA, PLS and OPLS
- Model interpretation
- Possibilities for on-line process monitoring

Course Date | Venue:

14 -15 March 2019

04 - 05 June 2019

05-06 September 2019

Also available on request
Sartorius Office, Singapore

Course Duration:

2 Days

Start First Day: 9:00am

End Last Day: 5:00pm

Trainer:

David Wang

Course Fee:

Please contact your local representative

Condition:

You are required to bring your own computer for the exercises

E1

Software



Software

Design of Experiments (DOE)

Who should attend: Researchers, scientists and engineers from all sectors of industry and academia.

Typical applications include product development, process improvement, optimization, validation and quality control. No prior knowledge of statistics is assumed.

Design of Experiments (DOE) is the most effective method to achieve product and process efficiency and optimization. MODDE is a state-of-the-art design of experiments software package that is used by scientists, engineers, and statisticians alike to help understand complex processes and products.

Course Objectives:

On completion, participants will know how to:

- Create efficient experimental designs to match the objectives
- Analyze experimental data using sound statistical principles
- Use graphical and statistical tools for understanding
- Communicate the results in graphical format
- Use numerical and graphical optimization tools
- Understand limitations and possibilities in conflict situations
- Create follow up experiments when applicable

Topics include

- Understanding the DOE concept
- Data modeling and diagnostics
- Create a solid base for decisions
- Finding the optimal setting

Course Date | Venue:

11-12 March 2019

06 - 07 June 2019

03 - 04 September 2019

Also available on request

Sartorius Office, Singapore

Course Duration:

2 Days

Start First Day: 9:00am

End Last Day: 5:00pm

Trainer:

David Wang

Course Fee:

Please contact your local representative

Condition:

You are required to bring your own computer for the exercises

Sartorius Lab Academy

Customized Training for Laboratory Professionals

In addition to our detailed intensive training program, Sartorius also offers concise and targeted training sessions at your company, called Lab Academy. These sessions are ideally suited to get an overview about specific applications and to enhance the efficiency of your workflows. Our experts provide a high percentage of hands-on training to extend your techniques and to give you valuable advice and tips for your daily work and challenges.

Choose from several course topics about lab water purification systems, pipettes, filtration and balances.

What you can expect:

- 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 - Custom-tailored 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

The Program

- Sartorius | Pipetting Academy
- Sartorius | Weighing Academy

For more information
or to register for a training session,

contact
caryn.chai@sartorius.com

Trainers

Sherwin Ting is responsible for the fermentation business of Southeast Asia, focusing on product proposals and application support. Prior to joining Sartorius, Sherwin was in the R&D industry for 8 years, focusing on upstream process development for stem cells. Sherwin has completed his Bachelor Degree in Biomedical Engineering from Boston University, United States.

Eileen Sim graduated from University of Queensland, Australia with Bachelor of Biotechnology (IIA). Prior to joining Sartorius, Eileen was in the R&D industry for 7 years, focusing on upstream process development for stem cells and also vaccine development.

In her current role, her primary responsibility will be to develop FRT business in South East Asia focusing on product proposals and application support.

Vicky Neo is responsible for the purification business of Southeast Asia. Prior to joining Sartorius, Vicky worked at Lonza (UK and Singapore) for 5 years in both Development Services and Manufacturing divisions. Her specialties include downstream process optimisation for scale up, membrane chromatography, technology transfer across sites and process improvement projects. She has an Honor's Degree in Chemistry and Biological Chemistry.

Lucas Loy graduated from Nanyang Technological University, Singapore, with a Bachelor Degree in Biological Sciences. Prior to joining Sartorius, he worked in Amgen for 4 years and took part in the tech transfers of 2 molecules from Amgen US to Singapore and was the Process Lead for Bulk Drug Filling and Harvest process. In his current position with Sartorius, Lucas aims to help emerging and existing biopharma companies to integrate single-use technologies into their processes.



Trainers

Tan Ming Wei is responsible for the filtration business within SEA, providing filter related training & technical support, identifying opportunities for filtration optimization and streamlining of processes. He started off his career at Baxter Bioscience as a Biotechnologist in Upstream Cell Culture | Media Prep | Inoculation.

Over the years, Ming Wei progresses to a Process Engineer, involved largely in tech transfer and critical investigation relating to product quality and product yield.

Caryn Chai completed her Bachelor Degree in Biomedical Science from University of Western Australia. Prior to joining Sartorius, Caryn was in the R&D industry for 10 years, focusing on cancer science and cardiovascular disease, with the goal of identifying novel targets for therapeutic intervention.

Caryn is a ACTA certified trainer.

David Wang is a Principal Data Scientist at Sartorius-Stedim Data Analytics based in Singapore. He mainly works on improving efficiency, optimality, and profitability for product development, process development, and manufacturing processes by applying advanced data analytics technology, where the work is focused on data-driven modelling, process monitoring, advanced process control, design of experiments, PAT and QbD. He has delivered a large number of data analytics training courses and industry projects across Asia-Pacific region.



Inhouse Training

We'll be happy to hold any of our on-going training courses and seminars inhouse at your own facilities. Of course, we can also put together complete new training courses for you! Let us know how we can help you to develop and implement a training concept that is right for you.

The benefits of our inhouse training include:

1. You get personalized training courses – designed according to your wishes and tailored to meet your needs.
2. You choose the dates and times for your training courses and seminars that suit.
3. A group of several colleagues can take part in the courses – all at the same time, without any added travel expenses.
4. The courses will be held using your own equipment and materials (e.g. cell lines). The knowledge gained can be put right into practice – on site!

Training Laboratories



Registration



Course Participant:

Name
Institute Company
Address
City
State
Phone
Email

Sartorius Office
10 Science Park Road, The Alpha
#02-13/14
Singapore Science Park II
Singapore 117684

Singapore Polytechnic
T11A504C
500 Dover Road
Singapore 139651

Invoice Address (if different from the one above):

Name
Institute Company
Address
City
State

Phone +65.6431.0530
+65.9782.2258
Email caryn.chai@sartorius.com

Invoice Address (if different from the one above):

Name
Institute Company
Address
City
State

Areas of Interest | Questions | Remarks

Date Signature

Information on How to Register

Registration is subject to our General Terms and Conditions of Business.

Registration

Registrations must be submitted in writing. A written confirmation of the booking will follow. The fees will be invoiced after completion of the seminar.

Cancellations

Cancellations will be accepted if made in writing up to 30 workdays prior to the beginning of the course. After this time, the course fee is payable in full. However, registered participants always have the option to appoint a substitute to attend.

Changing the Scope of Services

We reserve the right to change the content and schedule of the seminar and the trainer while preserving the overall nature of the seminar.

We reserve the right to cancel the seminar or to offer an alternative date. If we cancel the seminar due to illness of the trainer, acts of God, or other adventitious events, the registrants have no claim to having the seminar conducted.

Copyrights

Seminar materials and folders are governed by copyright laws and under no circumstances and at no time may they be reproduced by photographic or electronic means. These materials are intended for the sole personal use of the course participants.

2019 Training Calendar

January

February

March

11 - 12 March
Design of Experiments (DOE)

14 - 15 March
Multivariate Data Analysis

14 - 15 March
Basics of Fermentation

April

May

June

04 - 05 June
Multivariate Data Analysis

06 - 07 June
Design of Experiments (DOE)

19 - 20 June
Crossflow Filtration (Module 1)

21 June
Crossflow Filtration (Module 2)

July

August

September

03 - 04 September
Design of Experiments (DOE)

05 - 06 September
Multivariate Data Analysis

19 - 20 September
Single-Use Technologies

October

November

December

04 - 05 December
Basics of Filters and Integrity Testing

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Southeast Asia Main Offices

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Road,
The Alpha, #02-13/14
Singapore Science Park 2
Singapore 117684
Phone +65.6872.3966
Fax +65.6778.2494
Email info.singapore@sartorius-stedim.com

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4 Technology Park
Malaysia Bukit Jalil
57000 Kuala Lumpur
Phone +60.3.8996.0622
Fax +60.3.8996.0755
Email info.sea@sartorius-stedim.com

Thailand

Sartorius Thailand Co. Ltd.
129 Rama 9 Road, Huaykwang
Bangkok 10310
Phone +66.2643.8361-6
Fax +66.2643.8367
Email info.sea@sartorius-stedim.com

Vietnam

Sartorius Vietnam Co., Ltd
Bitexco Financial Tower
Unit 4631, Level 46,
2 Hai Trieu Street, Ben Nghe ward, District 1,
Ho Chi Minh City, Vietnam
Phone +84.8.6287.5951
Email info.sea@sartorius-stedim.com

Southeast Asia Distributors

Indonesia

PT Sartonet Filtrasi Indonesia
Jl. Kemanggisan Ilir Raya No.22, RT.19/RW.8,
Kemanggisan, Palmerah, Kota Jakarta Barat,
Daerah Khusus Ibukota Jakarta 11480, Indonesia
Phone +62.21.5366.9338/41/42/43
Fax +(62-21)53669339
Email sartonet@cbn.net.id

Philippines

Membrane Separation Tech. (MST) Trading, Inc.
1208 Richville Corporate Tower
Madrigal Business Park
1107 Alabang-Zapote Rd. Ayala Alabang
Muntinlupa City 1780
Phone +632.8098594 / 7763607
Fax +632.8093653
Email julibell_dupo@msttradinginc.com

Thailand

Renown Technical Co., Ltd.
1194 Soi Wachirathamsathit 57 Sukhumvit 101/1
Road Bangchak Prakanong
Bangkok 10260, Thailand
Phone +66 (0) 2185 4333 ext 2
Fax +66 (0) 2332 9390
Email rt_marketing@spcgroup.co.th

Vietnam

Indochina International JSC,
HanoiOffice
No.27-29 Gamuda Gardens 2-2 Str,
Tran Phu Ward, Hoang Mai Dist,
Ha Noi City, Vietnam.
Phone 04.6291.7941/04.6291.7942
Email trang.dinh@ibiotech.com.vn



◀▶ www.sartorius.com