

Tecan Training 2018.

THE KEY TO SUCCESS



Tecan's comprehensive training program enables customers to build up the necessary knowledge to optimize and maximize the productivity of Tecan devices. These training courses offer the trainees a thorough learning experience. The participants will benefit from Tecan training resources and obtain at first hand, the skills needed in their daily work.

BENEFITS

- maximized profitability justifies your training investment
- highly effective learning environment without the distractions of daily business
- certified and highly skilled trainers
- opportunities to ask questions and challenge what you are learning
- learn from experiences of other participants
- time for practicing
- ongoing access to trainers and product experts
- hands-on training using the instruments
- small training groups



TRAINING TERMS AND CONDITIONS

- **Training dates and course fees are published on the corresponding training webpages under www.tecan.com/training_schedule.**
- Course fees include full documentation, refreshments and lunch. A social event is included if course duration is more than three days.
- The courses require a minimum number of three participants.
- Registration is required four weeks prior to the start date of any course.
- Cancellation with less than 10 working days notice incurs a 50 per cent fee. Hotel fees arising due to cancellations with less than three days notice will be charged.
- Each training course includes a final written test. Successful participants will be awarded a training certificate. Participants who fail the test will be awarded a training confirmation. To take part in the test, full attendance of the course is mandatory.
- Customers should register via their local Tecan subsidiary or distributor.
- The charge for courses held at the customer's site will be communicated on request. Accommodation and travel expenses for the trainer will be charged separately.
- It may be necessary for reasons beyond our control to alter the venue, timetable or content of a training course. Fees will be refunded should any event be cancelled by Tecan. We accept no liability for any other costs incurred.
- Special terms and conditions apply for the Tecan Academy.



CONTENTS

Training structure overview4

Applications and special courses

● Freedom EVOware® Te-Chrom™ Wizard
– on request5

Software training for liquid handling products

● Freedom EVOware Standard6
 ● Freedom EVOware Plus7
 ● EVO Logic™ application software, basic8
 ● Freedom EVOLution™ application software for
Freedom EVOlyzer® and
Freedom EVO® Clinical, basic9
 ● Freedom EVOware
Sample Tracking – on request10
 ● Freedom EVOware
Sample Oriented – on request11
 ● Freedom EVOware
Normalization Wizard – on request12
 ● Freedom EVOware
Hit Picking Wizard – on request13
 ● Freedom EVOLution application software for
Freedom EVO Clinical, – on request14
 ● FluentControl™15

Dedicated liquid handling training courses

● Basic Liquid handling for Freedom EVO® 16
 ● Air LiHa Liquid handling for Freedom EVO 17
 ● Liquid handling with low volume option
for Freedom EVO – on request 18
 ● MCA96 Liquid handling for Freedom EVO 19
 ● MCA384 Liquid handling for Freedom EVO20
 ● Pressure Monitored Pipetting (PMP™)
for Freedom EVO – on request21
 ● QC Kit for Freedom EVO – on request 22
 ● FCA Liquid Handling for Fluent® 23
 ● MCA384 Liquid Handling for Fluent24
 ● Infinite® multimode reader product training –
on request25
 ● Basic Magellan™ for
Infinite F50 and Sunrise™26
 ● Advanced Magellan™
for high end Infinite Readers27

Separation training courses

● ProfiBlot™ – on request28
 ● HydroSpeed™ washer – on request29
 ● HydroFlex™ washer – on request30

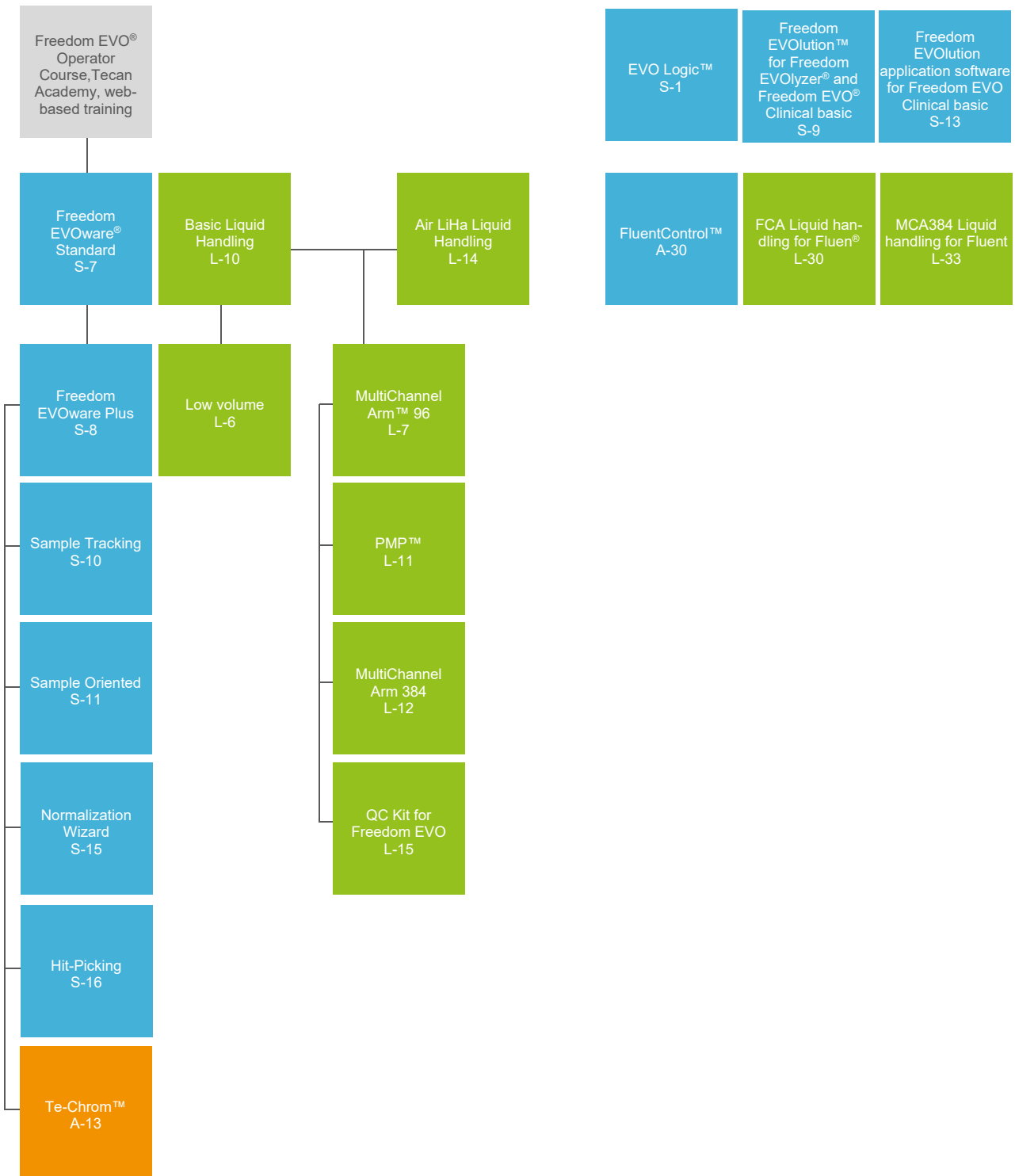
Microarray training courses

● Array suite (HS, LS, PS, ArrayPro) –
on request31

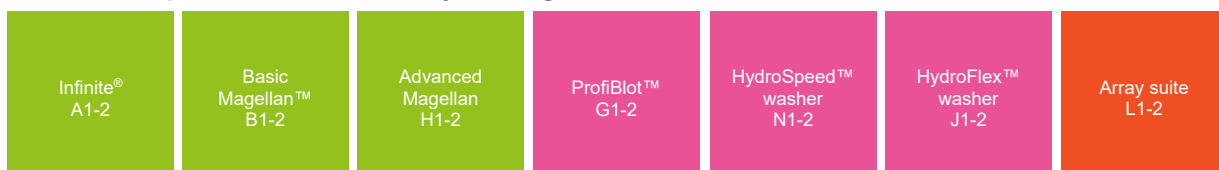


CONTENTS

Training courses for liquid handling products



Detection, separation and microarray training courses*



* these courses are only available for Tecan Distributors and Tecan Sales force



COURSE
NO. A-13

FREEDOM EVOWARE® TE-CHROM™ WIZARD

On request

Description

This course includes theory with a strong focus on hands-on training and will provide participants with convenient step-by-step instructions for implementing and optimizing processes using the Freedom EVOware Te-Chrom Wizard with Te-Chrom hardware. The course aims to develop knowledge, helping users to gain experience and become confident in using the Freedom EVOware Te-Chrom Wizard to control the instrument.

Target audience

This course is ideal for those who will run and support Te-Chrom applications on Freedom EVO instruments. Successful completion certification of this training course qualifies the application specialists/sales oriented staff/users to setup and optimize the Freedom EVOware Te-Chrom Wizard, including hardware installation.

Prerequisites

S-7 Freedom EVOware Standard

Duration

2 days

Language

English

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- software concept/structure
- terminology
- integration of Te-Chrom hardware and Te-Chrom Wizard
- importance of SnS setup and Freedom EVOware Te-Chrom-specific RoMa vectors
- installing and configuring the specific Te-Chrom drivers
- how to use the Wizard
- how to implement different example configurations of Te-Chrom
- overview of the different Te-Chrom wizard liquid classes
- error handling
- exercises
- collaboration and interface between Tecan and Atoll
- integration of the Te-Chrom Wizard and Freedom EVOware Sample Tracking software
- final test and certification

Upon course completion

At the end of the course, trainees will be able to implement and optimize processes on Freedom EVO instruments using Freedom EVOware Te-Chrom Wizard



COURSE
NO. S-7

FREEDOM EVOWARE® STANDARD

Description

This course includes some theory and has a strong focus on hands-on training. This training will provide participants with convenient step-by-step instructions for implementing or programming scripts, as well as optimizing existing applications with Freedom EVOware software. The course aims to develop knowledge, helping users to gain experience and become confident in using Freedom EVOware Standard software. This course is a prerequisite for all instructor based classroom training.

Target audience

This course is ideal for those who need to implement and optimize applications on Freedom EVO® Standard software.

Prerequisites

Freedom EVO Operator Certificate, basic web-based training course provided at the Tecan Academy. Visit www.tecan.com/academy for more information. Microsoft Windows® and Microsoft Office experience

Duration

4 days

Language

English

Training dates

Please refer to www.tecan.com/training or email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf, Switzerland
Tecan Germany GmbH, Crailsheim, Germany
Tecan UK, Reading, UK
Tecan US, Durham, NC, USA

Course outline

The following topics are covered in this module:

- introduction to Freedom EVO hardware
- installation of Freedom EVOware software
- system configuration
- the software components (editors) and their interaction
- settings and data structure
- configuration of worktable, carriers and labware with LiHa and RoMa arms
- defining applications using a wide variety of commands
- introduction to liquid handling
- how to use worklists
- data export
- defining and modifying RoMa vectors
- application optimization / troubleshooting
- maintenance procedures
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- implement and optimize applications on Freedom EVO instruments with Freedom
- EVOware Standard software understand the importance of maintenance and operating safety for perfect performance of the instrument



COURSE
NO. S-8

FREEDOM EVOWARE® PLUS

Description

This course includes some theory and has a strong focus on hands-on training. This training will provide participants with convenient step-by-step instructions for implementing and optimizing processes with the Freedom EVOware Plus software. The course aims to develop knowledge, helping users to gain experience and become confident in using Freedom EVOware Plus software to allow full control of the instrument.

Target audience

This course is ideal for those who need to implement and optimize processes on Freedom EVO instruments.

Prerequisites

S-7 Freedom EVOware Standard

Duration

4 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland
Tecan Germany GmbH, Crailsheim,
Germany
Tecan UK, Reading, UK
Tecan US, Durham, NC, USA

Course outline

The following topics are covered in this module:

- overview of the Freedom EVOware software concept
- creating, prescheduling and running processes
- plate handling including lid handling
- implementation of devices
- optimizing processes with Freedom EVOware Plus scheduling capabilities
- setting up variable steered processes
- worklist handling
- application optimization/troubleshooting
- barcode handling
- final test and certification

Upon course completion

At the end of the course, trainees will be able to implement and optimize processes on Freedom EVO instruments using Freedom EVOware Plus software.



COURSE
NO. S-1

EVO LOGIC™ APPLICATION SOFTWARE, BASIC

Description

This course enables the trainee to implement and program methods and assays in EVO Logic on Freedom EVO instruments. It will provide participants with convenient step-by-step instructions to help them in their day-to-day work. The course includes some theory and has a strong focus on hands-on training. The aim is to develop skills and techniques, helping users to gain experience and become confident in using Logic software to control instruments.

Target audience

This course is ideal for those who wish to develop the skills to implement and set up tests and methods with EVO Logic software, and want to optimize and maximize the productivity of their applications.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

4 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- instrument hardware introduction
- overview of the EVO Logic concept
- set-up configuration
- how to create assays
- organizing tests in a method
- liquid handling concept with liquid LiHa and Air LiHa
- working with the database
- EVO Logic applications: sample distribution and pooling
- error handling within EVO Logic
- working in batch, joblist or ASTM mode
- maintenance procedures
- operation and safety
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- set-up a configuration in EVO Logic
- implement tests and methods in EVO Logic
- set up a range of different applications in EVO Logic like pooling and sample preparation for ELISA assays.
- understand the importance of maintenance /operation safety in EVO Logic



COURSE
NO. S-9

FREEDOM EVOLUTION™ APPLICATION SOFTWARE FOR FREEDOM EVOLYZER® AND FREEDOM EVO® CLINICAL, BASIC

Description

This course includes some theory and has a strong focus on hands-on training. This training will provide participants with convenient step-by-step instructions for installing, setting up and running assays in Freedom EVOLution software for the Freedom EVOLyzer and the EVOClinical. The course aims to develop knowledge, helping users to gain experience and become confident in using Freedom EVOLution software to allow full control of the Freedom EVOLyzer and the EVOClinical instrument.

Target audience

This course is ideal for those who need to program and optimize assays on Freedom EVOLyzer and the Freedom EVOClinical instruments. Upon request, we also provide separate "Freedom EVOLution application software for Freedom EVOClinical" (S-13, 5 days) or separate "Freedom EVOLution application software for Freedom EVOLyzer" trainings (9 days).

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

10 days

Language

English

Training dates

Please refer to www.tecan.com/training or email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf, Switzerland

Course outline

The following topics are covered in this module:

- introduction to the Freedom EVOLyzer and Freedom EVOClinical hardware and Freedom EVOLution application software
- installation of the application software
- link between instrument set-up and service software, and the application software
- the software components (editors) and their interaction
 - configuration builder
 - consumable data editor
 - process definition editor
 - HydroControl™ (washer software, as part of the EVOLution for EVOLyzer software package)
 - Magellan™ Tracker (reader software, as part of the EVOLution for EVOLyzer software package)
 - Freedom EVOLution studio
 - additional tools
 - programming ELISAs
- working modes, file structure and output files
- result files and evaluation with Magellan Tracker
- working with schedulers

- basics of liquid handling
- basics of troubleshooting and error handling
- advanced application features
- maintenance procedures
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- implement and optimize applications on Freedom EVOLyzer and EVOClinical instruments with the Freedom EVOLution software
- understand the importance of maintenance/operating safety for perfect performance of the instrument



COURSE
NO. S-10

FREEDOM EVOWARE® SAMPLE TRACKING

On request

Description

This training will provide participants with convenient step-by-step instructions for implementing and optimizing EVOware scripts and processes with Freedom EVOware Sample Tracking. The course aims to develop knowledge, helping users to gain experience and become confident in using Freedom EVOware Sample Tracking.

Target audience

This course is ideal for those who wish to develop the skills to implement and optimize Sample Tracking applications on Freedom EVO® instruments.

Prerequisites

S-7 Freedom EVOware Standard

Duration

2 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- software concept /structure
- terminology
- installation
- barcode handling
- Freedom EVOware integration (labware registration, labware rotation and print report)
- data interpretation
- set up customized report queries with the Report Designer
- Sample Tracking Web Reporting Tool
- Sample Tracking Quick Viewer
- error handling
- final test and certification

Upon course completion

At the end of the course, trainees will be able to implement and optimize processes on Freedom EVO instruments using Freedom EVOware Sample Tracking.



COURSE
NO. S-11

FREEDOM EVOWARE® SAMPLE ORIENTED

On request

Description

This course includes both theory and hands-on training. This training will provide participants with convenient step-by-step instructions for implementing and optimizing processes with Freedom EVOware Sample Oriented. The course aims to develop knowledge, helping users to gain experience and become confident in using Freedom EVOware Sample Oriented.

Target audience

This course is ideal for those who wish to develop the skills to implement and optimize sample oriented applications on Freedom EVO instruments.

Prerequisites

S-7 Freedom EVOware Standard

Duration

2 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

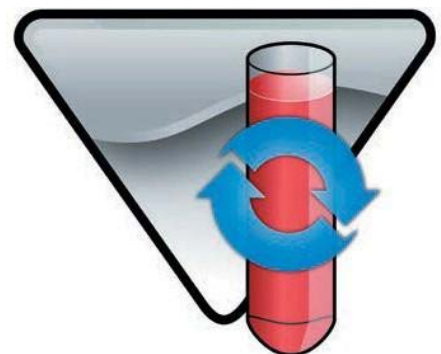
Course outline

The following topics are covered in this module:

- overview of Freedom EVO platform
- overview of Freedom EVOware Sample Oriented
- software structure and terminology
- wizards
- dilution schemes
- input /output files
- joblists
- assay implementation
- final test and certification

Upon course completion

At the end of the course, trainees will be able to implement and run scripts and processes on Freedom EVO instruments using Freedom EVOware Sample Oriented.



COURSE
NO. S-15

FREEDOM EVOWARE® NORMALIZATION WIZARD

On request

Description

This course will provide participants with convenient step-by-step instructions for implementing and optimizing processes with the Freedom EVOware Normalization Wizard. It includes some theory, and has a strong focus on hands-on training, and aims to develop knowledge, helping users to gain experience and become confident in using the Freedom EVOware Normalization Wizard.

Target audience

This course is ideal for those who wish to develop the skills to implement and optimize applications on Freedom EVO® instruments using the Freedom EVOware Normalization Wizard.

Prerequisites

S-7 Freedom EVOware Standard

Duration

1.5 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

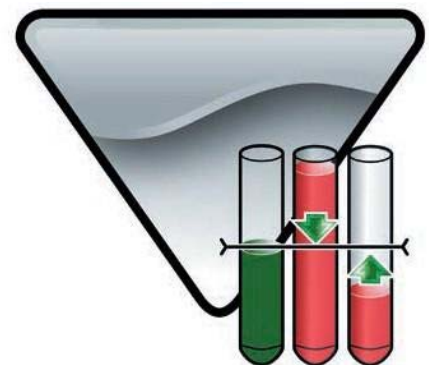
Course outline

The following topics are covered in this module:

- software concept /structure
- terminology
- installation
- integration of Quantitation Wizard (sample dilution and distribution, control and standard addition, quantification)
- interface with Magellan™ and output files
- integration of the Normalization Wizard
- calculation report (including conversion of .xml to .csv format)
- error handling
- sample exercise
- final test and certification

Upon course completion

At the end of the course, trainees will be able to implement and optimize processes on Freedom EVO instruments using the Freedom EVOware Normalization Wizard.



COURSE
NO. S-16

FREEDOM EVOWARE® HIT-PICKING WIZARD

On request

Description

This course will provide participants with convenient step-by-step instructions and hands-on-training for implementing and optimizing processes with the Freedom EVOware Hit-Picking Wizard. The course aims to develop knowledge, helping users to gain experience and become confident in using the Freedom EVOware Hit-Picking Wizard to manage full control of the instrument.

Target audience

This course is ideal for those who wish to develop the skills to implement and optimize applications on Freedom EVO instruments using the Freedom EVOware Hit-Picking Wizard.

Prerequisites

S-7 Freedom EVOware Standard

Duration

1.5 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- software concept /structure
- terminology
- installation
- wizard configuration
- choosing the source and destination labware
- integrating a sample list file
- additional options (dilutions, reagent addition, use of controls)
- plate layout
- error handling
- sample exercise
- final test and certification

Upon course completion

At the end of the course, trainees will be able to implement and optimize processes on Freedom EVO instruments using the Freedom EVOware Hit-Picking Wizard.



COURSE
NO. S-13

FREEDOM EVOLUTION™ APPLICATION SOFTWARE FOR FREEDOM EVO® CLINICAL

On request

Description

This training will provide the participants with convenient step-by-step instructions for installing, setting up and running assays in the Freedom EVOLution software for the Freedom EVO Clinical. It includes both theory and a strong focus on hands-on training. We help to establish confidence, develop knowledge, enable powerful usage of the tools and gain appropriate experience with the Freedom EVOLution application software for the Freedom EVO Clinical.

Target audience

This course is designed as a basic training to get to know Freedom EVOLution software for the Freedom EVO Clinical and learn to program assays on Freedom EVO Clinical instruments.

To learn how to work with and setup tests on the Freedom EVOLution software for Freedom EVOlyzer instruments, please refer to course number S-9.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

5 days

Language

English

Training dates

Please refer to www.tecan.com/training or email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf, Switzerland

Course outline

The following topics are covered in this module:

- introduction to the Freedom EVO Clinical hardware and the Freedom EVOLution application software
- installation of the application software
- link between instrument set-up and service, and the application software
- the software components (editors) and their interaction
 - configuration builder
 - consumable data editor
 - process definition editor
 - Freedom EVOLution studio
 - additional tools
- work modes, file structure and output files
- work with scheduler
- basics troubleshooting and error handling
- advanced application features
- maintenance procedures
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- implement applications on Tecan Freedom EVO Clinical instruments
- understand the importance of maintenance/operating safety for optimal performance of the instrument



COURSE
NO. A-30

FLUENTCONTROL™

Description

This course includes both theory and hands-on training on the Fluent® Instrument using the FluentControl Application Software. It will provide participants with convenient step-by-step instructions to help them in their day-to-day work. The aim is to develop skills and techniques, helping operators to gain experience and become confident in using FluentControl.

Target audience

This course is ideal for those who need an introduction to FluentControl and wish to develop skills to understand applications on Fluent Instrument.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

4 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland
Tecan Germany GmbH, Crailsheim,
Germany
Tecan UK, Reading, UK
Tecan US, Durham, NC, USA

Course outline

The following topics are covered in this module

- Introduction to instrument hardware
- Introduction to FluentControl Application Software
- Set-up and configuration
- Workspace concept, how to configure worktable layout
- Overview of commands for FCA, RGA and MCA384
- Application Optimization and Troubleshooting
- TouchTools
- Introduction to Liquid Handling
- Test and Certification.

Upon course completion

At the end of the course, trainees:

- understand and optimize applications on Fluent instruments with FluentControl
- troubleshoot errors at runtime
- understand the importance of maintenance and operating safety for perfect performance of the instrument



COURSE
NO. L-10

BASIC LIQUID HANDLING FOR FREEDOM EVO®

Description

This course contains both theory and hands-on training with step-by-step instructions for Tecan instruments with liquid displacement system (liquid-filled tubing). The trainee will establish a broad knowledge of liquid handling skills for effective use of pipetting instruments. The course will focus on improving the quality of pipetting and maintenance. This course is the basis for advanced user courses L-6, L-7, L-11, L-12, L-14 and L-15.

Target audience

This course is ideal for those who operate a Tecan instrument with a liquid displacement system and would like to optimize and maximize the performance of their system.

Prerequisites

Familiarity with Tecan Freedom EVO instrument and S-7 Freedom EVOware® Standard

Duration

2.5 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland
Tecan Germany GmbH, Crailsheim,
Germany
Tecan UK, Reading, UK
Tecan US, Durham, NC, USA

Course outline

The following topics are covered in this module:

- overview of hardware relevant to pipetting
- the liquid system of Tecan instruments using standard fixed and disposable tips
- the liquid handling language of Tecan instruments
- wash and flush procedures
- importance of maintenance
- the liquid handling concept of the Tecan application software packages
- the influence of the various liquid handling parameters on pipetting quality
- how to implement a new liquid class
- single and multipipetting modes
- measuring precision with photometric tests
- troubleshooting
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand the default liquid classes
- create customized liquid classes
- implement the concept of liquid classes for their own applications
- improve the performance of their instruments with appropriate maintenance



COURSE
NO. L-14

AIR LIHA LIQUID HANDLING FOR FREEDOM EVO®

Description

This course combines theory and hands-on training, with step-by-step instructions for the use of Tecan instruments equipped with an **air displacement pipetting arm (Air LiHa)**. The trainee will acquire a broad knowledge of liquid handling skills to enable the effective use of the pipetting arm, as well as learn how to improve the quality of pipetting.

Target audience

This course is ideal for those who operate and maintain a Tecan Air LiHa on a Freedom EVO® instrument, and would like to optimize and maximize the performance of their system.

Prerequisites

S-7 Freedom EVOware® Standard and L-10 Basic Liquid Handling for Freedom EVO

Duration

2.5 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland
Tecan US, Durham, NC, USA

Course outline

The following topics are covered in this module:

- introduction to the Air LiHa hardware
- Tecan instruments' air displacement system, using disposable tips
- the liquid handling concept of the Tecan application software packages
- the influence of various liquid handling parameters on pipetting quality
 - free dispense and contact pipetting
 - aspiration/dispense speeds and delays
 - acceleration and deceleration
- single and multi-pipetting modes
- single-channel calibration
- overview of the Air LiHa MultiSense™ option (see L-11 Pressure Monitored Pipetting (PMP™))
- use of conductive (cLLD) and pressurebased (pLLD) liquid level detection
- measuring precision with photometric tests
- daily, weekly and preventive maintenance
- troubleshooting
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand the default liquid classes for the Air LiHa on the Freedom EVO
- create customized liquid classes
- implement the concept of liquid classes for their own applications
- improve the performance of their instrument by optimizing the parameter settings



COURSE
NO. L-6

LIQUID HANDLING WITH LOW VOLUME OPTION FOR FREEDOM EVO®

On request

Description

This course is based on basic liquid handling course L-10 and is dedicated to low volume pipetting (<3 µl). It contains both theory and hands-on training, with step-by-step instructions.

Target audience

This course is ideal for those who operate Freedom EVO instruments and would like to optimize and maximize pipetting of small volumes.

Prerequisites

L-10 Basic Liquid Handling for Freedom EVO

Duration

2 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- important hardware considerations for the low volume option
- the liquid handling system of Tecan instruments with low volume options
- influence of various liquid handling parameters on pipetting quality of small volumes
- use of low volume options with low volume standard fixed tips and 10 or 50 µl disposable tips
- liquid handling with and without pinch
- single and multi-pipetting modes
- measuring precision with photometric tests
- wash and flush procedures
- daily, weekly and preventive maintenance
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand the default liquid classes for low volume options
- improve the performance of their instruments with appropriate maintenance
- implement the concept of liquid classes for their own applications with the low volume option
- know how to improve the liquid handling of small volumes (<3 µl) with and without the pinch valve



**COURSE
NO. L-7**

MCA96 LIQUID HANDLING FOR FREEDOM EVO®

On request

Description

This course is designed to develop the ability to precisely, accurately and safely pipette with the MCA96 option on a Freedom EVO instrument. This training should develop knowledge and establish confidence for programming liquid handling with fixed and disposable tips using Freedom EVOware®. This course will provide participants with convenient step-by-step instructions to help them in their day-to-day work. It includes some theory with a strong focus on hands-on training.

Target audience

This course is ideal for those who operate a Freedom EVO instrument equipped with an MCA96 arm and would like to optimize and maximize the performance of their system.

Prerequisites

Familiarity with Tecan Freedom EVO instrument and L-10 Basic Liquid Handling for Freedom EVO

Duration

3 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- MCA96 overview
- set-up and configuration
- liquid handling with MCA96 options
- wash steps /disposable tips / fixed tips
- wash procedures
- liquid classes, with and without contact pipetting
- single and multi-pipetting modes
- aspiration and dispense speeds and delays
- precision and accuracy measurements (color performance test)
- daily, weekly and preventive maintenance
- service information
- results and conclusions
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand the liquid classes for standard and disposable tips with a MCA96 device
- create customized liquid classes with and without contact pipetting
- improve the performance of their instruments with appropriate maintenance



COURSE
NO. L-12

MCA384 LIQUID HANDLING FOR FREEDOM EVO®

Description

This course is designed to help users develop the skills to precisely, accurately and safely pipette with a MCA384 option. It aims to develop knowledge and establish confidence in programming liquid handling with fixed and disposable tips using Freedom EVOware®. The course will provide participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This course is ideal for those who operate a Freedom EVO instrument equipped with an MCA384 arm and would like to optimize and maximize the performance of their system.

Prerequisites

Familiarity with Tecan Freedom EVO instrument and L-10 Basic Liquid Handling for Freedom EVO

Duration

3 days

Language

English

Training dates

Please refer to www.tecan.com/training or email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf, Switzerland

Course outline

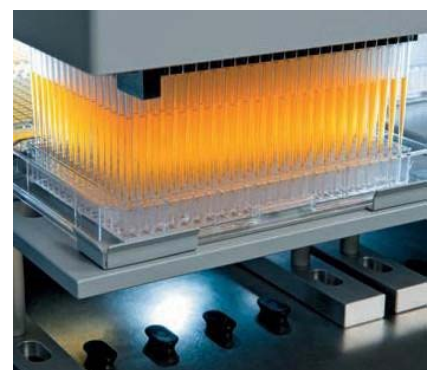
The following topics are covered in this module:

- MCA384 overview
- set-up and configuration
- liquid handling with the MCA384
- the influence of the various liquid handling parameters on pipetting quality
 - free dispense and contact pipetting
 - aspiration respectively dispense, speeds and delays
 - acceleration and deceleration
- liquid classes, with and without contact pipetting
- single and multi-pipetting modes
- working with flexible numbers of tips (8, 16, 24, 96 or 384)
- exchanging different adapter plates 'on the fly'
- wash procedures with fixed tips and disposable tips
- precision and accuracy measurements with photometric tests
 - colour performance test and QC Kit
- daily, weekly and preventive maintenance
- Troubleshooting
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand liquid classes for an MCA384 using standard and disposable tips
- create customized liquid classes with and without contact pipetting
- improve the performance of their instruments with appropriate maintenance
- optimize wash procedures
- perform precision and accuracy measurements with colour and the QC Kit performance tests



COURSE
NO. L-11

PRESSURE MONITORED PIPETTING (PMP™) FOR FREEDOM EVO®

On request

Description

This course contains both theory and hands-on training with step-by-step instructions. The trainee will establish knowledge about the function of the MultiSense™LiHa on a Freedom EVO instrument. The course will focus on pressure-based liquid level detection (pLLD) and pressure control of pipetting quality (PMP).

Target audience

This course is ideal for those who use the PMP option on their Freedom EVO instrument equipped with MultiSense™LiHa arm and would like to optimize and maximize the performance of their system.

Prerequisites

Familiarity with Tecan Freedom EVO instrument and L-10 Basic Liquid Handling for Freedom EVO

Duration

3 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- overview of the MultiSense LiHa hardware
- PMP concept for 200 µl and 1000 µl disposable tips
- configuring and testing the PMP option in service and application software
- the use of pLLD
- pLLD parameters in liquid class editor (pLLD or combined pLLD /cLLD)
- detection of substantial errors during pipetting (based on physical model algorithms)
- the influence of the various PMP parameters in the liquid class
- short introduction into the evaluation of pressure curves
- overview of the evaluation configuration manager, and the algorithm to detect pipetting errors
- daily, weekly and preventive maintenance
- troubleshooting
- final test and certification

Upon course completion

At the end of the course, trainees will be able to understand the use of pLLD, the function of PMP and create customized liquid classes for PMP.



COURSE
NO. L-15

QC KIT FOR FREEDOM EVO®

On request

Description

This course includes both theory and a strong focus on hands-on training, and is available as a standalone training course, bookable on request only. The training course will provide participants with convenient step-by-step instructions for implementing and operating the **QC Kit components** and software on Freedom EVO platforms.

Target audience

This course is ideal for those who operate and service the QC Kit on Freedom EVO instruments, and would like to optimize and maximize the performance of their system.

Prerequisites

S-7 Freedom EVOware® Standard

Duration

1 day

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

The following topics are covered in this module:

- introduction to QC Kit on Freedom EVO
- QC Kit Wizard
- QC Kit consumables, such as dye solutions and verification plates
- QC Kit hardware components; calibrator plate, barcode scanner, shaker and reader
- installation of the QC Kit devices
- link between QC Kit and the Freedom EVOware® application software, and their integrated operation
- using QC Kit in online or offline mode
- using different Tecan pipetting options with the QC Kit
- result evaluation with QC Kit
- recommended use of the Tecan QC Kit
- troubleshooting
- maintenance
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- implement and operate the QC Kit
- understand the importance of maintenance/operating safety for the precise performance of the Tecan instruments



COURSE
NO. L-30

FCA LIQUID HANDLING FOR FLUENT®

Description

This course contains both theory and hands-on training with step-by-step instructions for Tecan Fluent instrument with **liquid displacement system (FCA)** and/or with an **air displacement pipetting arm (Air FCA)**. The trainee will establish a broad knowledge of liquid handling skills for effective use of pipetting instruments. The course will focus on improving the quality of pipetting and maintenance.

Target audience

This course is ideal for those who operate a Tecan Fluent instrument with a liquid displacement system and would like to optimize and maximize the performance of their system.

Prerequisites

Familiarity with Tecan Fluent instrument and A-30 FluentControl™

Duration

2.5 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

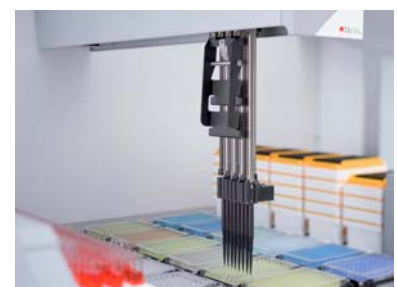
The following topics are covered in this module:

- overview of hardware relevant to pipetting
- the liquid system of Tecan Fluent instrument using standard fixed and disposable tips
- the liquid handling language of Tecan instruments
- wash and flush procedures
- importance of maintenance
- the liquid handling concept of the Tecan application software packages
- the influence of the various liquid handling parameters on pipetting quality
- how to implement a new liquid class
- single and multi-pipetting modes
- measuring precision with photometric tests
- troubleshooting
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand the default liquid classes
- create customized liquid classes
- implement the concept of liquid classes for their own applications
- improve the performance of their instruments with appropriate maintenance



COURSE
NO. L-33

MCA384 LIQUID HANDLING FOR FLUENT®

Description

This course is designed to help users develop the skills to precisely, accurately and safely pipette with a MCA384 option. It aims to develop knowledge and establish confidence in programming liquid handling with fixed and disposable tips using FluentControl™. The course will provide participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This course is ideal for those who operate a MCA384 instrument and would like to optimize and maximize the performance of their system.

Prerequisites

Familiarity with Tecan Fluent instrument and A-30 FluentControl™

Duration

2.5 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-ch@tecan.com

Venue

Tecan Schweiz AG, Männedorf,
Switzerland

Course outline

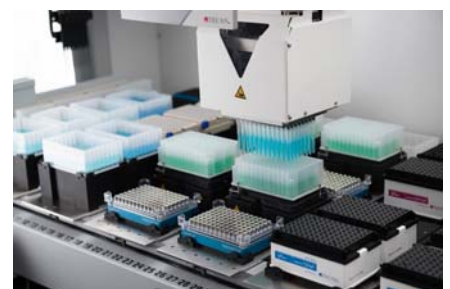
The following topics are covered in this module:

- MCA384 overview
- set-up and configuration
- liquid handling with the MCA384
- the influence of the various liquid handling parameters on pipetting quality
 - free dispense and contact pipetting
 - aspiration respectively dispense, speeds and delays
 - acceleration and deceleration
- liquid classes, with and without contact pipetting
- single and multi-pipetting modes
- working with flexible numbers of tips (8, 16, 24, 96 or 384)
- exchanging different adapter plates 'on the fly'
- wash procedures with fixed tips and disposable tips
- precision and accuracy measurements with photometric tests
 - colour performance test and QC Kit
- daily, weekly and preventive maintenance
- Troubleshooting
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- understand liquid classes for an MCA384 using standard and disposable tips
- create customized liquid classes with and without contact pipetting
- improve the performance of their instruments with appropriate maintenance
- optimize wash procedures
- perform precision and accuracy measurements with colour and the QC Kit performance tests



**COURSE
NO. A 1-2**

INFINITE® MULTIMODE READER PRODUCT TRAINING

On request

Description

The Infinite multimode reader product training course includes theoretical and practical knowledge and has a strong focus on hands-on training. The training goal is to develop skills and techniques to manage and fully control the instruments of the Infinite reader series and efficiently set up applications on every reader, alone or in combination with the Connect™ or built-in stacker. This training provides participants with convenient step-by-step instructions to help them in their daily work.

Target audience

This fundamental course is designed to develop skills for the efficient and powerful use of Infinite multimode readers, i.e. the Infinite 200 PRO, Infinite F500 and Infinite M1000. It is ideal for people who want to operate the Infinite instruments, to establish basic as well as advanced applications, and to maximize the productivity of their system.

Prerequisites

Understanding of basic measurement principles (e.g. absorbance, fluorescence, luminescence, fluorescence polarization), experience with Windows® and Microsoft® Office, esp. Excel for standard calculations (average, st. dev. CV, etc.)

Duration

5 days

Language

English

Training dates

Please refer to www.tecan.com/training or send an email to training-at@tecan.com.

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

The following products will be trained in the course:

- Infinite 200 PRO
- Infinite 200 PRO NanoQuant
- Infinite F500
- Infinite M1000
- NanoQuant Plate
- Connect, built-in Stacker
- injectors
- i-control™ software

The following topics are covered in this module:

- technical concepts of filter- and monochromator-based Infinite instruments (including Quad4 Monochromators™ technology)
- measurement techniques:
 - free absorbance
 - fluorescence
 - fluorescence polarization
 - luminescence
- advanced applications:
 - free RET (e.g. GeneBLAzer®)
 - TRF (e.g. Delfia®)
 - TR-FRET (HTRF®)
 - FP (e.g. PolarScreen™) hands-on experiments
- example assays and script set-up (e.g. DLR, HTRF, LanthaScreen™)
- optimization of measurement parameters (e.g. attenuation, gain, mirror and filter selection)
- final test and certification

Upon course completion

At the end of the course, trainees will:

- be able to operate the instruments of the Infinite multimode reader series with i-control software
- be able to operate the instruments of the Infinite multimode reader series in combination with the Connect stacker
- have a good general knowledge about the most common measurement techniques and advanced applications
- create, modify and optimize measurement processes
- receive a training certificate after performing the practical exercises and successfully passing a final test



**COURSE
NO. B 1-2**

BASIC MAGELLAN™ FOR INFINITE® F50 AND SUNRISE™

Description

This course includes some theory and has a strong focus on hands-on training. It aims to develop knowledge and skills, helping users to gain experience and become confident in managing and controlling the instruments. It teaches the skills required to efficiently set up applications on Magellan using an Infinite F50 or Sunrise instrument. This training will provide participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This fundamental course is ideal for those who operate Sunrise or Infinite F50 instruments via Magellan. It is recommendable for sales specialists and field service engineers, who want to gain basic knowledge of Magellan software

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

2 days

Language

English

Training dates

Please refer to
www.tecan.com/training or
email training-at@tecan.com

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

The following topics are covered in this module:

Basic measurement techniques:

- absorbance

Infinite F50:

- technical overview of instrument
- instrument options
- quality control plate and software: MultiCheck™ F50

Sunrise:

- technical overview of instrument
- instrument options
- quality control plate and software: QC Pac™
- hands-on: basic absorbance applications

Magellan software:

- Software overview
- Magellan wizards
- hands-on: qualitative and quantitative ELISAs
- Magellan Clinical/Tracker (for implementation of regulatory requirements of IVDD and FDA 21 CFR part 11)

Upon course completion

At the end of the course, trainees will be able to:

- operate Infinite F50 and Sunrise instruments with Magellan
- define Magellan methods for qualitative and quantitative ELISA's
- use Magellan Clinical/Tracker



COURSE
NO. H 1-2

ADVANCED MAGELLAN™ FOR HIGH END INFINITE® READERS

Description

This course is dedicated to setting up advanced Magellan applications using high end instruments of the Infinite series. This course includes some theory and has a strong focus on hands-on training, providing participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This course is designed for those who want to develop the skills required to become power users of Magellan software. It is recommendable for application specialists.

Prerequisites

Basic Magellan knowledge

(definition of plate layout, transformations, cut-off, QC validation)

Knowledge about measurement techniques (absorbance, fluorescence intensity, fluorescence polarization)

Microsoft Windows® and Microsoft Office experience

Duration

3 days

Language

English

Training dates

Please refer to

www.tecan.com/training or
email training-at@tecan.com

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

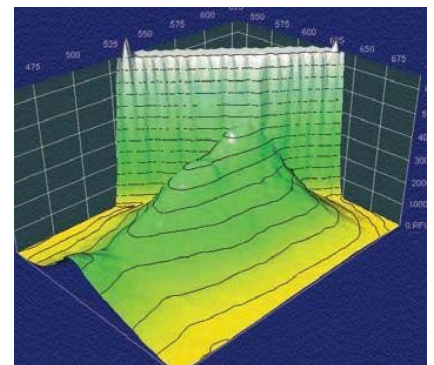
The following topics are covered in this module:

- Magellan applications:
 - kinetic evaluation
 - multi-label and advanced transformations
 - concentration and dilution series (ICx calculation)
 - standard curves
 - calculation for TR-FRET assays and HTRF®-calculation
 - calculation for FP assays (considering G-factor calibration)
 - spectra data reduction
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- modify and optimize existing processes
- set up high-end applications with the Magellan software
- receive a training certificate after successfully passing a final test and performing practical exercises



COURSE
NO. G 1-2

PROFIBLOT™

On request

Description

This course teaches the skills required to efficiently set up applications on ProfiBlot instruments. It includes some theory and a strong focus on hands-on training. The course aims to develop skills and techniques, helping users to gain experience and become confident in managing and controlling ProfiBlot instruments. This training provides participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This fundamental course is ideal for those who wish to optimize and maximize the productivity of their ProfiBlot system.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

2.5 days

Language

English or German (according to preference of the participants)

Training dates

Please refer to
www.tecan.com/training or
email training-at@tecan.com

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

The following topics are covered in this module:

- ProfiBlot overview
- use of instrument
- on-board programming
- BlotWare™ software with examples
- hands-on training
- programming procedures
- applications
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- set up ProfiBlot instruments and processes
- modify and optimize existing processes
- receive a training certificate after successfully passing a final test and performing practical exercises



COURSE
NO. N 1-2

HYDROSPEED™ WASHER

On request

Description

This course includes some theory and has a strong focus on hands-on training. The course aims to develop skills, helping users to gain experience and become confident in fully managing and controlling these instruments. This training provides participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This course is ideal for those who wish to optimize and maximize the efficiency of their HydroSpeed washer.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

2 days

Language

English or German

Training dates

Please refer to
www.tecan.com/training or
email training-at@tecan.com

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

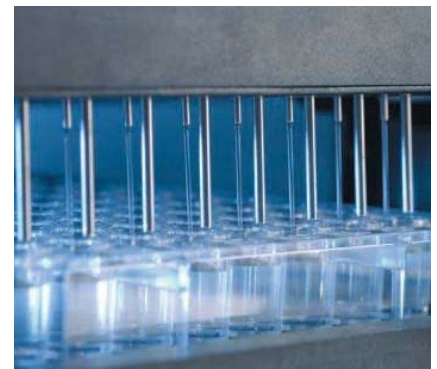
The following topics are covered in this module:

- overview of the HydroSpeed (plate washer)
- technical background (e.g. ELISA plate washing, ultrafiltration, magnetic bead separation, drip-mode cell washing)
- on board programming
- HydroControl™ software with examples
- hands-on training
- programming procedures
- applications
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- set up washer instruments
- modify and optimize existing programs and settings
- receive a training certificate after successfully passing a final test and performing practical exercises



COURSE
NO. J 1-2

HYDROFLEX™ WASHER

On request

Description

This course includes some theory and has a strong focus on hands-on training. The course aims to develop skills, helping users to gain experience and become confident in fully managing and controlling the instrument. This training provides participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This course is ideal for those who wish to optimize and maximize the efficiency of their HydroFlex washer.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

2 days

Language

English or German

Training dates

Please refer to
www.tecan.com/training or
email training-at@tecan.com

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

The following topics are covered in this module:

- HydroFlex (strip washer) overview
- technical background (e.g. ELISA plate washing, ultrafiltration, magnetic bead separation, drip-mode cell washing)
- on board programming
- HydroControl™ software with examples
- hands-on training
- programming procedures
- applications
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- set up HydroFlex washer instruments
- modify and optimize existing processes
- receive a training certificate after successfully passing a final test and performing practical exercises



**COURSE
NO. L 1-2**

ARRAY SUITE (HS, LS, PS, ARRAYPRO)

On request

Description

This course will provide participants with detailed technical information and step-by-step instructions to help them with their product installations. This course includes both theory and hands-on training. The course aims to develop skills and techniques, helping users to gain experience and become confident in managing and controlling these instruments. It teaches the skills required to efficiently set up protocols on the HS Pro series of hybridization stations, as well as the LS and PS series of laser scanner instruments.

The course provides a short overview showing how to use the sophisticated array analysis software ArrayPro and how to set up customized protocols. This training provides participants with convenient step-by-step instructions to help them in their day-to-day work.

Target audience

This fundamental course is designed to develop skills for basic usage of the HS 4800™ Pro, the HS 400™ Pro, the Power-Scanner™, LS Reloaded™ and ArrayPro products. It is ideal for those who operate such instruments and would like to optimize the productivity of their system.

Prerequisites

Microsoft Windows® and Microsoft Office experience

Duration

5 days: 2 days HS Pro, 2 days laser scanners (LS, PS), 1 day ArrayPro analysis demonstration – parts of the trainings can be booked separately.

Language

English

Training dates

Please refer to www.tecan.com/training or email training-at@tecan.com

Venue

Tecan Austria GmbH, Grödig, Austria

Course outline

The following topics are covered in this module:

- HS 4800 Pro, HS 400 Pro:
 - free background array methodology
 - key features of instruments
 - instrument overview (technical construction)
 - installation
 - operating software
- LS Reloaded:
 - instrument overview (technical construction)
 - image capture and analysis software
 - key features
 - hands-on training for set-up and QC of instrument
 - hands-on training slides and plate protocol design /scan
- PowerScanner:
 - instrument overview (technical construction)
 - image capture and analysis software

- key features
- hands-on training for set-up and QC of instrument
- batch run HD arrays with different protocols
- ArrayPro :
 - key features of image capture and analysis software
 - run template /experiment wizard
- final test and certification

Upon course completion

At the end of the course, trainees will be able to:

- set up HS 4800 Pro / HS 400 Pro and laser scanner instruments (LS and PS), as well as ArrayPro analysis software
- troubleshoot
- modify and optimize existing processes
- receive a training certificate after successfully passing a final test and performing practical exercises
- use the Freedom EVOware® Normalization Wizard



THE FOLLOWING TECAN SUBSIDIARIES OFFER CERTAIN USER TRAINING COURSES

Australia (on request)

Tecan Australia Pty Ltd.
21 / 3 Westside Avenue
Port Melbourne Vic 3207
Australia
T 1300 808 403 (free in Australia)
T +61 3 9647 4100
helpdesk-aus@tecan.com

Austria (on request)

Tecan Austria GmbH
Untersbergstrasse 1a
A-5082 Grödig, Austria
T +43 6246 8933 256
helpdesk-at@tecan.com

Benelux (on request)

Tecan Benelux bvba
Businesspark E19
Bettelsesteenweg 455
B 2800 Mechelen, Belgium
T +32 15 709 055 (french) or
T +32 15 709 054 (english) or
T +31 20 708 4773
helpdesk.benelux@tecan.com

China (on request)

Tecan (Shanghai) Trading Co.,
Ltd. Room 1802-1804, No.388,
Fushan Road, Pudong New Area
200122 Shanghai, P.R. China
T +86 21 2206 32 06 or
T +86 4008 213 888
helpdesk-cn@tecan.com

France (on request)

Tecan France S.A.S.
6, Avenue du Château de Gerland
F-69007 Lyon, France
T +33 820 88 77 36
helpdesk-fr@tecan.com

Germany

Tecan Deutschland GmbH
Werner-von-Siemens-Strasse 23
D-74564 Crailsheim, Germany
T +49 1805 8322 633
helpdesk-de@tecan.com

Great Britain

Tecan UK
Theale Court, 11-13 High Street,
Theale, UK-Reading RG7 5AH,
United Kingdom
T +44 118 930 0300
helpdesk-uk@tecan.com

Spain (on request)

Tecan Ibérica (Spain) Edificio
Mapfre
C/ de la Marina 16-18,
Planta 11a C-1
08005 Barcelona, Spain
T +34 93 40 91 237
helpdesk-sp@tecan.com

Italy (on request)

Tecan Italia S.r.l.
Via Brescia 39
I-20063 – Cemusco Sul Naviglio
(MI), Italy
T +39 800 11 22 91
helpdesk-it@tecan.com

Japan (on request)

Tecan Japan Co. Ltd Kawasaki
Tech Center 580-16,
Horikawa-cho, Saiwai-ku
Kawasaki, Kanagawa 212-0013,
Japan
T +81 44 556 7312 (Kawasaki)
T +81(0) 6305 8511 (Osaka)
helpdesk-jp@tecan.com

Scandinavia (on request)

Tecan Nordic AB
Taljegårdsgatan 11B
SE-43153 Mölndal, Sweden
T +46 317 54 40 00
helpdesk@tecan.se

Switzerland

Tecan Schweiz AG
Seestrasse 103,
CH-8708 Männedorf, Switzerland
T +41 44 922 82 82
helpdesk-ch@tecan.com

USA

Tecan US
9401 Globe Center Drive,
Suite 140
Morrisville NC 27560, USA
T +1 919 361 5200
Toll Free in the US:
T +1 800 TECAN US or
T +1 800 352 5128
helpdesk-us@tecan.com

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at <http://www.tecan.com/trademarks>. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

Tecan, Freedom EVO, Freedom EVOlyzer, Freedom EVOware and Infinite are registered and Freedom EVOLution, Gemini, FACTS, Logic, MultiChannel Arm, HID EVOLution, QC Pac, MultiCheck, Connect, MultiSense, Te-Chrom, HydroControl, i-control, Quad4 Monochromators, Sunrise, Magellan, ProfiBlot, HydroSpeed, HydroFlex, BlotWare, PMP, HS 4800 PRO, HS 400 PRO, PowerScanner and LS Reloaded are trademarks of Tecan Group Ltd., Männedorf Switzerland. Excel, Microsoft and Windows are registered trademarks of Microsoft Corporation. DLR is a trademark of Promega Corporation. HTRF is a registered trademark of Cisbio International. GeneBLAzer is a registered and PolarScreen and LanthaScreen are trademarks of Invitrogen Corporation, Carlsbad, USA. Delfia is a registered trademark of PerkinElmer, Inc., Wellesley, USA. © 2016 Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com



 **TECAN.**