

Data Analytics eLearning: Level 2 Intermediate Training on Design of Experiment (DOE) and Multivariate Data Analysis (MVDA)

These Level 2 intermediate courses focus on either Design of Experiments (DOE) or Multivariate Data Analysis (MVDA). Design of Experiments (DOE) is a rational and cost-effective approach to practical experimentation that allows the effects of critical process parameters to be assessed using only the minimum of resources. MVDA is the key to unlocking the information residing in complex process datasets and presenting the results in a simple graphical format. Both DOE and MVDA empower scientists to better understand, develop

and optimize processes more efficiently, leading to better product quality and reduced development costs.

These Level 2 training courses are a direct continuation of the Level 1 courses for DOE and MVDA. For interested participants new to DOE or MVDA or feel that a knowledge refresher is needed, it is recommended that the respective Level 1 training is completed first before embarking on the Level 2 course.

Cost and Conditions

The Level 2 Intermediate course fee is 500 € per person per course. Please contact your local sales representative to receive a quote.



Course Objective

The objective of the Level 2 DOE course is to provide a deeper understanding of DOE, by focusing on more elaborate design protocols as well as specific application designs supported by MODDE® to handle constrained situations and non-standard demands such as the D-optimal design family, which includes Mixture design common in formulation application, REDMUP® designs for DOE in microtiter plates, generalized sub-set design for accomplishing time distributed, sequential DOE protocols, and stability testing DOE for ensuring or establishing product shelf-life.

The objective of the Level 2 MVDA course is to provide a deeper understanding of MVDA and cover additional functionality that comes with SIMCA®, such as model exploration tools (What-If, Score Space Explorer, and OPLS® Solver), specific model types and applications, and further theory. In addition, the batch evolution model (BEM), the batch level model (BLM), the multi-block orthogonal components (MOCA) model, and the hierarchical model framework will be covered in this course.

Who Should Participate?

The Level 2 course is an intermediate course, ideally suited for existing MODDE® or SIMCA® users from all sectors of industry and academia, and intended for those who have already completed the respective Level 1 DOE and | or MVDA courses.

Language Information

The eLearning courses are given in English.

Schedule

A self-paced, on-demand online training for both courses is available for a period of 3 months from the time of activation.

Prerequisite Software, Course Material and Certification

The course material will be provided within the training platform. Each participant will need SIMCA® (for MVDA), and | or MODDE® (for DOE) software installed on their own computer with an active license as a prerequisite before taking these training courses. The software and licenses will not be included with the Level 2 courses. Each course ends with a Final Test. This test can be taken multiple times. A course certificate for each course is awarded after passing the test.

Order Information

Article Number	Description	Price
UT-OC-2496	Data Analytics eLearning MVDA (Level 2)	500 €
UT-OC-2497	Data Analytics eLearning DOE (Level 2)	500 €

Please contact your local sales representative to receive a quote.

Also, in the Data Analytics eLearning Series:

SIMCA®-online Operator Training Course

Are you confident that your process is on track to meet quality targets? Learn how SIMCA®-online uses multivariate data analysis (MVDA) modeling to eliminate the complexity of traditional process monitoring so you can get clear, actionable insights – exactly when you need them, no matter where you are. Find out how you can better understand typical use cases for real-time monitoring, increase process visibility from end-to-end, accurately implement end point prediction, forecasting, early fault detection and warning of process deviations. This self-paced eLearning is intended for end users (operators) of SIMCA®-online. There is no course fee required to access this training. An insight into the basics of MVDA is recommended, but not required. If you would like to find out more about SIMCA®-online, please contact your local Sartorius sales representative.

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178



For more information, visit
[sartorius.com](https://www.sartorius.com)