



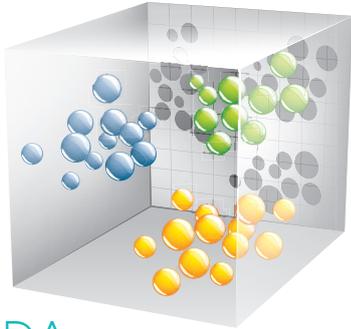
MVDA EduPack for Multivariate Data Analysis

We help you see what
others don't using SIMCA®

Simplifying Progress

SARTORIUS

Reduce Complexity and Learn from Your Data



MVDA
Multivariate
Data Analysis

Multivariate Data Analysis (MVDA) are methods to reduce complexity in your data and reveal information hidden in the numbers. The information is transformed into graphs that are easily interpreted. Multivariate methods are useful for presenting an overview of the data, identify reasons for deviations and predict the properties of the data that you want to investigate. The value of using MVDA is thorough understanding of historical data enabling you to learn from the past and improve the future.

Visualize Information in Your Data

MVDA is valuable when you don't know where to start analyzing all of your historical data. Applying MVDA methods will reduce noise in the data and provide information rich graphs where important patterns are easily detected. These patterns can be investigated further by the use of interactive graphics to understand the source of pattern. With this insight, you can visualize and explain your historical data.

DOE EduPack & MVDA EduPack

We provide EduPack with two complementary methodologies for gaining knowledge from data, Design of Experiments (DOE) and Multivariate data analysis (MVDA). DOE provides experimental series that reveals cause and effect relations between input variables and process output. MVDA can be used to extract information from big data sets with historical data.



“Build knowledge and understanding the smart way - to learn the most from your data”

Education to Master your Data

The education package starts with an introduction to MVDA and how the methods can be applied to solve problems and gain information. With EduPack students get the best education in MVDA for generating good results.

MVDA EduPack is designed to give your students hands-on skills to solve problems and learn:

- How to explore data and look deeply into how changing parameters can affect outcomes
- How to visualize trends and clusters using an easy-to-understand graphical interface
- How to analyze, define and identify problems
- How to use sophisticated modeling techniques to create accurate predictions about product quality
- How to interpret and gain in-depth understanding of your data and learn from the data

Get Started with MVDA

The MVDA EduPack will help you get started with all the educational material you need for lectures and hands-on training. Slides for classroom teaching are supported with 12 recorded lectures that will cover theory and examples. The slides are structured, so theory and exercises go together.

MVDA EduPack includes:

- Complete slide deck for classroom teaching
- 12 recorded lectures
- Exercises with relevant examples from industry
- Software: SIMCA® Multivariate Data Analysis Solution including one license with 20 activations valid for three years



Sartorius Data Analytics – Change a Little. Grow a Lot.

Sartorius Data Analytics are leading data analytics experts with multiple applications in high-tech industries worldwide. We enable customers at both the individual and enterprise levels to maximize the value of their data, enabling better decision-making and process excellence. Through our complete solutions in the Umetrics® Suite of Data Analytics Solutions, which include state of the art software, academy training and education, consulting, and OEM services, we help our customers to grow their businesses by maximizing profitability and minimizing time to market.

To learn more about how Sartorius Data Analytics helps you grow your business please visit www.sartorius.com/umetrics and follow us on LinkedIn.



Sartorius Data Analytics
Phone: +46.40.664.258.0
E-mail: umetrics@sartorius.com
www.sartorius.com/umetrics