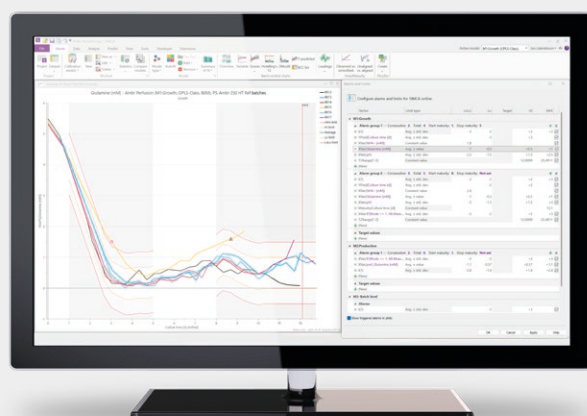


Umetrics® Digital Twin Build SIMCA® 18.2

Data Exploration Workflows,
Project Registry, and
Analytics Extensions



Product Information

SIMCA® 18.2 enhances data analytics with new features like data exploration workflows, allowing users to visualize and select data more effectively, and model lifecycle management, a model registry for tracking and managing models throughout their lifecycle.

The integration with SIMCA®-online ensures seamless predictive process control, optimizing yield and minimizing costs while fostering a user-friendly experience based on customer feedback.

Features and Benefits

- Seamlessly connect to and explore data in the cloud for data discovery
- Generate new variables to enhance data analysis and discovery
- Contextualize data for deeper understanding
- Track project versions and compare results directly in SIMCA® for greater control and accuracy

Introduction

Relevant Applications

- **Data exploration**— Understand, contextualize, select, and share data
- **Model lifecycle management registry**— Store, compare, version, and control created models

Relevant Process Steps

- Any upstream or downstream process in discovery, development, or manufacturing
- Supporting tools before model building
- Supporting tools for models from model-building

Overview

The Umetrics® Digital Twin AI Ecosystem equips you with end-to-end tools and support to accelerate results across your entire process. Users gain access to supporting documents, templates, training, and consultation to address their specific business challenges. Each year, our courses and webinars help more than a thousand professionals build expertise and confidence in extracting real value from their data.

A complete AI Ecosystem for Business Growth

The Umetrics® Digital Twin AI ecosystem is a family of proven data analytics, advanced process control, and application innovation solutions that work seamlessly together.

Umetrics® Digital Twin Build

Tools for empirical machine learning models in SIMCA®, for mechanistic models in cell Insights, and for generating and exploring system with design of experiments in MODDE®.

Umetrics® Digital Twin Deploy

Tools for visualizing the process, ensuring process consistency, and both predicting and forecasting the product quality through machine learning models.

Umetrics® Digital Twin Innovate

Tools for self-service rapid prototyping, creation, and deployment of streamlined workflows in Scibox® and built-in machine learning in SIMCA®-Q, and design of experiments in MODDE®-Q.

Umetrics® Digital Twin Elevate

Training and consulting services to support knowledge building, usage, and maximization of value from the full AI ecosystem.

These solutions give you control and confidence in your processes at every stage—from development to manufacturing.

Change a Little—Grow a Lot

The Umetrics® Digital Twin AI ecosystem supports organizations in generating value from their data. It provides tools for data analytics that can help identify key factors influencing research, development, and manufacturing outcomes. By enhancing process understanding and product consistency, the solution enables reduced risk and accelerated time to market. The offering includes software, training, support, and project management. As part of Sartorius, a global company with more than 10,000 employees, the ecosystem is backed by an international presence and infrastructure.

Technical Specifications

System Requirements

- SIMCA® runs on all modern PCs
- Any x64 computer hardware supported by Windows
- Supported versions of Windows 11 or Windows 10, Microsoft Windows Server 2022, 2019, or 2016
- Microsoft .NET Framework 4.6.2 or later. The latest 4.8 version is recommended
- 700 MB available hard disk space for a full installation

Ordering Information

Peripherals and Accessories

SIMCA® Extensions

- DTA

Consultancy Service Packages

Description	Order Number
GxP Validation Framework for SIMCA®-online (document templates)	UT-110060005
Direct Transition Analysis (DTA) for DSP (TBC)	UT-110060007
Spectro Model Building L1 Service Package	UT-110060017
Spectro Model Building L2 Service Package	UT-110060018
Spectro Model Building L3 Service Package	UT-110060019
SIMCA®-online Model Updates Package	UT-110060020
SIMCA®-online Model Updates Package	UT-110060021

Training Courses

Description	Order Number
DOE MVDA Level 1 Basic eLearning Course	UT-OC-2495
MVDA Level 2 Intermediate eLearning Course	UT-OC-2496
Multivariate Data Analysis (MVDA) Open Web Course	UT-SC-2132
Multivariate Data Analysis (MVDA) Client Onsite Training	UT-310015
Multivariate Data Analysis (MVDA) Client Web Training	UT-SC-2125
SIMCA®-online Admin Course	UT-OC-ADM
SIMCA®-online Operator eLearning Course	n/a

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



Experience the benefits for your business today

Find out how our solutions can help your business to grow, whatever industry you are in.

Visit sartorius.com/umetrics for details or to download a free 30-day trial.