

MODDE®

Design of Experiments Solution

Readme MODDE® Pro 13.0.2

November 16, 2021

Thank you for your interest in MODDE® 13.

MODDE® Pro 13.0.2 is a minor release that addresses some stability issues found in MODDE® 13.

MODDE® includes state-of-the-art functionality for creating designs and analyzing the results with confidence, supporting the Quality by Design initiative through a set of useful tools.

New features in MODDE® 13:

- Optimization wizard (Pro only)
- Contextual properties pane. Most frequently functions easily accessible.
- Design recommendation based on Design Objective and specific requirements
- Favorite setpoint defined by user
- New response objective and condition
- Optimal selection of replicated design points
- Detailed design power
- Optimization within design space (Pro only)
- Visualization of desirability
- Setpoint comparison (Pro only)
- Definitive Screening Design improvement
- Sweet spot histogram

Detailed information about functionality in MODDE® is available in “Technical Features for MODDE 13” and at our [website](#).

Installation

Administrator permissions are necessary to install MODDE® Pro 13.0.2.

Licensing and Activation

To run MODDE® you need to have a valid license that has been activated. The first time MODDE® is started, if the activation key is entered in the installation wizard, MODDE® automatically attempts to activate your license and download it from a web service. The licensing system was updated in MODDE® 13.0.1. This means that if you have a license for 13.0.1 it works for 13.0.2 as well. A license from an earlier version does not work for 13.0.2.

If the automatic activation fails, please read “Manual product activation.pdf” located in the Program folder. With a license server, you need information from your local IT department on where the license server is located.

Virtual Machine

To run MODDE® Pro on a virtual machine, your license must specifically allow this. Contact your sales representative for more information.

Trial Limitations

MODDE® Pro can be run as a trial for 30 days.



Minimum System Requirements

- CPU: Intel Core i5-4460, 4:th generation, (3.20GHz) or AMD FX-6300
- RAM: 4 GB
- Disk space: 400 MB available disk space
- OS: Microsoft Windows 8.1 or 10, 64-bit only
- Video card: Nvidia GeForce GTX 460, ATI Radeon HD 4850, or Intel HD Graphics 4400

Support

The latest information and support questions can be found in [Sartorius Data Analytics Knowledge base](#).

For technical support, contact umetrics_support@sartorius.com.