

# MODDE

Design of Experiments Solution

## Readme MODDE Pro 13.0.1

18 May 2021

---

Thank you for your interest in MODDE® 13.

MODDE Pro 13.0.1 is a minor release that introduces a new licensing system. This release also 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 13.0.1.

## 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. Due to the change of licensing system, a previous MODDE 13 license will not work.

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](mailto:umetrics_support@sartorius.com).