

MODDE

Design of Experiments Solution

Technical features for MODDE 13

27 March 2024

Design of Experiments (DOE) is the most effective method to achieve product and process efficiency and optimization. MODDE® is a state-of-the-art design of experiments software package that is used by scientists, engineers, and statisticians alike to help understand complex processes and products. **Highlighted rows are new in MODDE 13 or 13.1.**

Feature

Design generation

Design Wizard guides the design generation

Four design objectives: Screening, **System Characterization**, Optimization (RSM), **Robust Verification**

Split objective

Up to 48 factors

Factor ranges (scaling) can be updated retroactively

Factor types: Quantitative, quantitative multilevel (24 levels) factors, qualitative factors (24 levels),

Factor types: Formulation (mixture), Filler

Constant and uncontrolled factors

Linear constraints on factors

Combination of process and formulation factors

128 responses possible

New response objective and condition

Linear, Log, Neglog, Logit, Exp and Power transformations of factors and responses

Detailed design power estimation

Optimal selection of replicated design points

A wide variety of classical designs: Fractional factorial, Full factorial (2 levels, 3 levels and mixed), L9, L18, L27, L36, CCF, CCC, CCO, Reduced CCF and CCC, Box Behnken, Rechtschaffner designs in 2 and 3 levels, Doehlert designs, regular and Super saturated Plackett Burman designs. Definitive screening designs.

Axial (reduced, normal, and extended), Cubic centroid (Mod, Mod w/face, Special and Full) designs

Reduced combinatorial designs (J2)

Generalized subset designs - optimal and balanced multilevel designs

Stability testing designs

Rectangular Experimental Designs for Multi-Unit Platforms, RED-MUP. Supports designs for up to 4 plates with sizes 8x12 and 16x24, with 32x48 size plate. Includes RED-MUP specific designs

D-Optimal designs using state of the art algorithm

Blocking of classical and D-Optimal designs

Inclusions can be imported and edited

Candidate sets can be read from file

Feature

Import design data from external files

Paste design to import it

Complementing designs, using classical and D-Optimal approaches

Onion designs from scores generated in SIMCA

Onion design in ordinary factors, both with imported candidate set and candidate set generated by MODDE

Analysis of worksheet including Scatter Plots, Histogram, Descriptive Statistics, Correlation Matrix, Replicate Plots and condition number

Qualitative factors with missing levels supported

Export and open in SIMCA

Analysis and modelling

Fit with MLR or PLS

Cox and Scheffé Mixture models

Handles process and mixture models and their combinations

Cross validation of models

Indication of confounded model terms for fractional factorial designs

Analysis guidance

Analysis wizard guides the user through the analysis step by step allowing model customization from the graphs

One-Click analysis feature, including automatic outlier detection, transformation and model tuning

Automatic Square and Interaction tests in the Analysis wizard

Advisor pane which explains analysis plots and results and advises you on what to do next

Reviewing the model

Multiplots and lists displaying selected responses

Summary of the model fit plot and list with Q², R², Model validity (LOF) and Reproducibility

Customizable model overview multiplot

ANOVA plots and lists

Residual vs Run Order, Predicted, Variable plots and lists

Normal Probability of residuals, Observed vs Predicted and Distance to Model plots

Coefficient plots and lists

Effects and Interaction plots

Variable importance (VIP) plots and lists

Score and Loading plots

Box Cox plot

Refining the model

Interactive pruning of model terms with automatic model fitting and updating of all open plots and lists

Automated model tuning feature

Separate model for each response with all fit methods

Predictions

Feature

Contour, Sweet Spot and Prediction plots wizards for simple generation of plots

Design Space plot wizard to find design space and robust setpoint

Visualization of desirability

2D, 3D (mixture) and 4D plots make it possible to display up to 5 factors simultaneously.

4D plots with qualitative factors on the outer axes

Contour surface with multiple responses

Option to lock contour levels in Contour plot

Prediction plot interval estimates include confidence, prediction and tolerance options

Prediction plots display raw data

Overlay prediction plots for multiple responses

Factor effects plot including confidence intervals

Prediction Scatter plot updated with changes in the Predictions spreadsheet

Transformed factors by default displayed in original units in prediction plots

Optimization guidance

Optimization wizard guides the user through the optimization step by step

Visualization of simulated process output

Favorite setpoint defined by user

Summary of alternative setpoints and statistics

Optimizer

Uses a multidimensional Simplex method

Customizable desirability functions

Possible to set target values and optimization criteria

Optimizer predicts possible ranges for all responses

Weighting according to the importance of the responses

Optimization of multiple responses, regular or derived

Risk analysis of the optimal setting

Option to set response limits as absolute in Optimizer

Robust optimization feature presenting the most robust setpoint

Response correlation effect optionally included in Design Space calculations

Optimization within design space

Design Space Explorer

Design space explorer plot expansion with hypercube to facilitate communication of the Proven Acceptable Range (PAR)

Export of complete Design Space as a data matrix

Setpoint validation

Statistical robustness validation of the investigated system

Interactive GUI and automatic functions for robust range establishment

Setpoint comparison histogram

Feature

Plots and lists

Contextual properties pane for easier access to plot properties

Predefined plot sizes when copying to various presentation types

Create list from plots

Color coding in lists to highlight suspicious values

Plots can be customized, and templates saved

Reports

Customizable report generator for fast and standardized documentation

Report integrated in the MODDE *.mip file

Experimental Design sharing

Share DOE investigations in json file format
