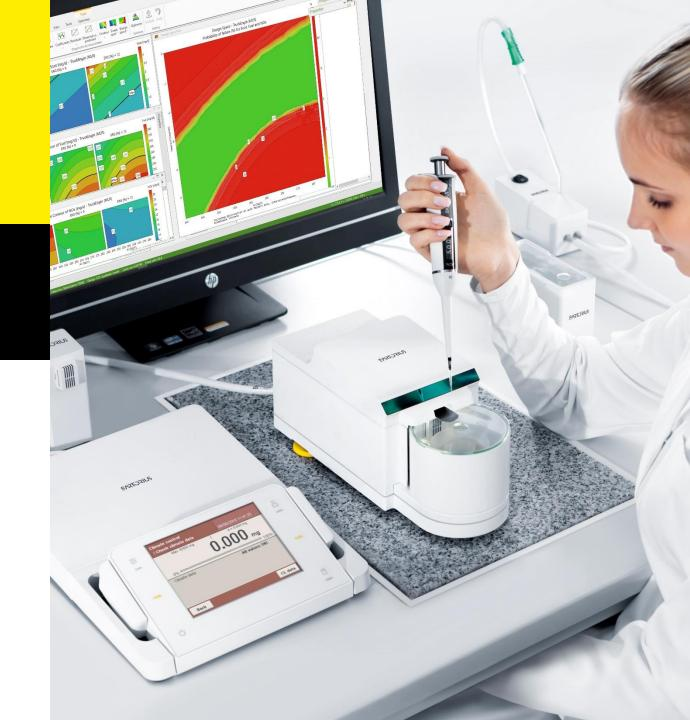
# SARTURIUS

Simplifying Progress

MODDE® 13 – What's New?

Release March 2, 2021



#### MODDE® 13 Overview

- Umetrics® suite MODDE® is focused on delivering a full design of experiment solution, from creating an investigation plan to analyze results and support decision making base on scientific principles.
- MODDE® 13 is focused on improving Design Selection and making Optimization easier and more powerful.
- With MODDE® 13 you get the new Optimization Wizard that complements the reworked Design Wizard and updated Analysis Wizard.
- The most frequently used function and customizations can now be accessed in the Properties pane making it easier than ever to succeed with design of experiments.



## MODDE® 13 Highlights

- Design Selection
  - New response Objectives and Conditions
  - Factor setting for Normal Operating Range
  - Detailed design Power per factor and response
  - Optimal selection of replicated design points
  - Interactive design selection view
  - New investigation objectives
- Model and Analysis
  - Model verification in Analysis Wizard
  - Visualization of desirability

- Optimization
  - Optimization Wizard
  - Optimization within design space
  - Setpoint comparison
- Other Improvements
  - Access to most used functions in new Properties panes
  - Define favorite setpoint
  - Create new plot from with existing settings for contour like plots
  - Performance



## Learning What's New in MODDE® 13

■ In the following slides you will get an overview of the changes and additions made in MODDE® 13

■ Please also check out more videos on MODDE® and other Umetrics® suite products by looking up Sartorius Data Analytics on YouTube



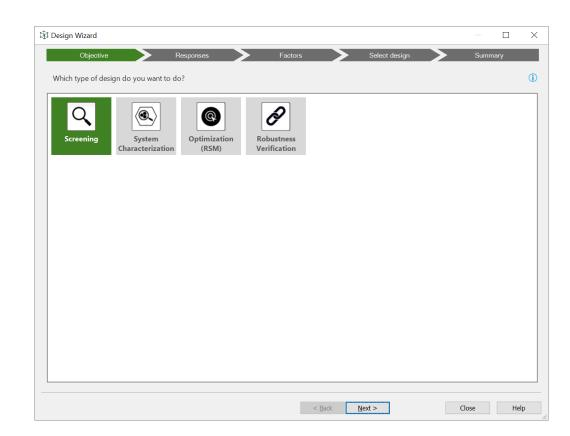
Simplifying Progress

MODDE® 13 What's New Design Selection



#### Design Wizard - Investigation Objective

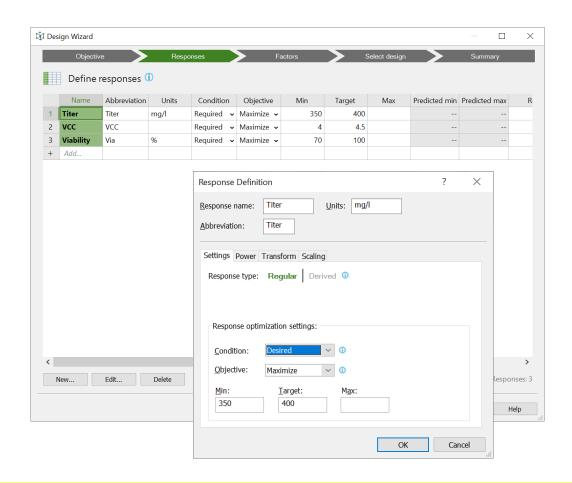
- Screening, first stage of an investigation when little is known
- System characterization, investigate the influence of the most influential factors, including their two-factor interactions and quadratic effects
- Optimization (RSM), optimization using the most influential factors and focus on low prediction error
- Robustness verification, investigate the system's sensitivity to changes in certain critical factors





#### Design Wizard - Responses definition dialog

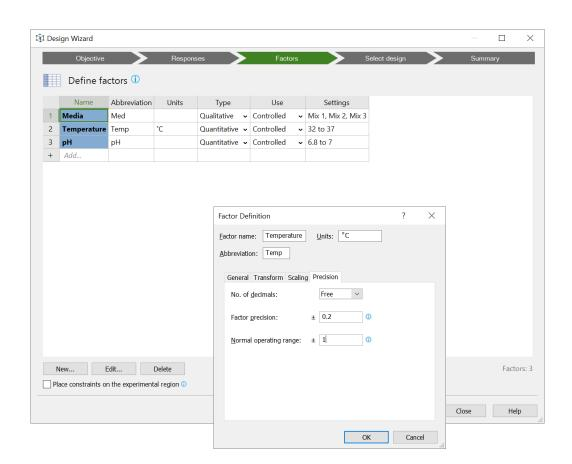
- Condition importance of response
  - Required responses are quality attributes with hard limits that must be fulfilled. They define sweet spot, design space and are used in robust setpoint calculations
  - Desired responses are influencing optimal setpoint
  - Observed responses are predicted but do not influence design space or optimization
- Objective what you want to do with the response.
  - Minimize the response, often with max limit
  - Maximize the response, often with min limit
  - Target set the response close to a value
  - Inside, not influencing optimal setpoint
- Power to detect a coefficient as significant
  - Signal to noise ratio





#### Design Wizard - Factor definition dialog

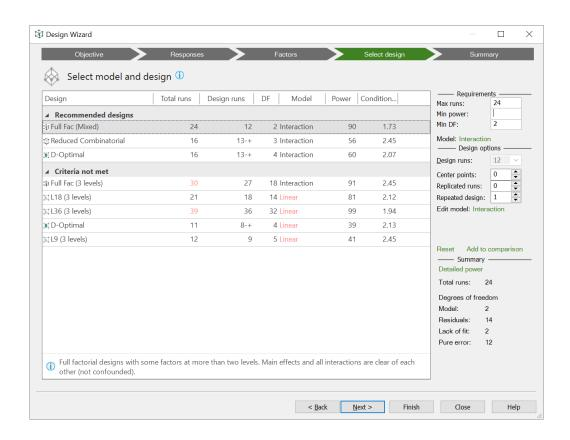
- Precision is the uncertainty in measurement of the factor setting
  - If set, it complements the prediction uncertainty range in Design Space estimation and Setpoint Exploration estimation
- Normal Operating Range, NOR, is defined as the common range during daily use of the application
  - Use the model to simulate response distributions
  - NOR can be used in used in Setpoint Comparison and Setpoint Exploration





#### Design Wizard - Select Design

- Requirements
  - Used to filter the design alternatives
  - Max number of runs, Design Power and Degrees of freedom
  - Model complexity
- Design options for selected design
  - Replicated runs, add best replicated design points
  - Edit model of a design
  - Add to comparison to add the current design to Compare designs section
- Other improvements
  - Create D-Optimal design directly from select design page
  - Detailed Power for each factor and response combination
  - Sort design on Design power or I-optimality







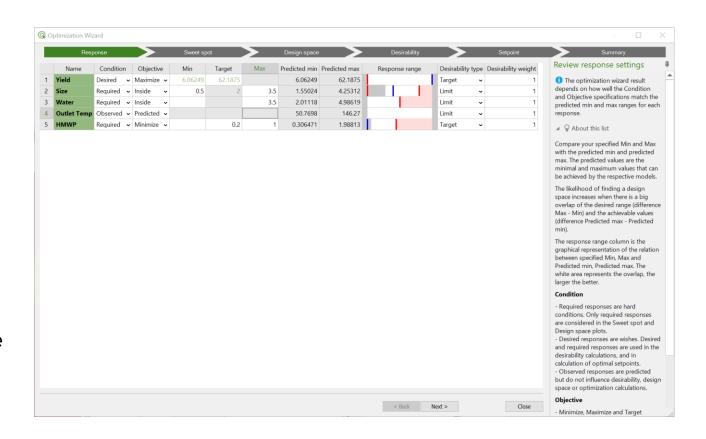
Simplifying Progress

MODDE® 13 What's New Optimization



#### Optimization Wizard - Inspect Response Settings

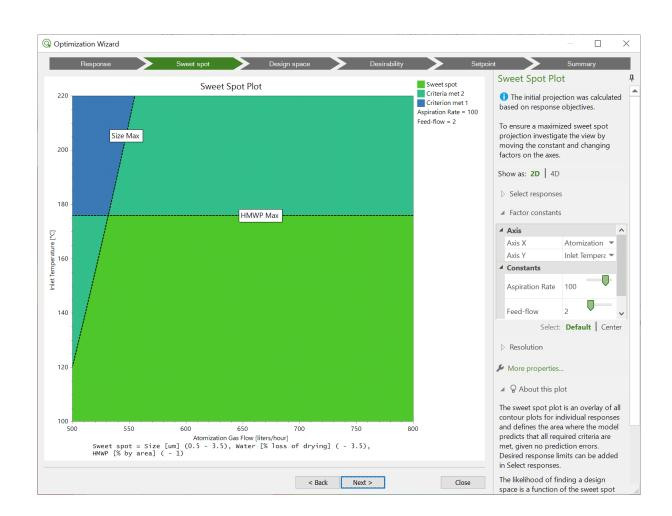
- Verify that response Condition, Objective and limits are set correctly
- Response condition, objective and limits are automatically fetched from the response definition. Green values are tooltips that indicate the range of values valid when considering the model
- Set desirability type
  - Target, reach a solution close to target
  - Limit, reach a solution within limits
- Weight can be used to limit the influence of some responses in order to find a compromise between many.





#### Optimization Wizard - Sweet spot

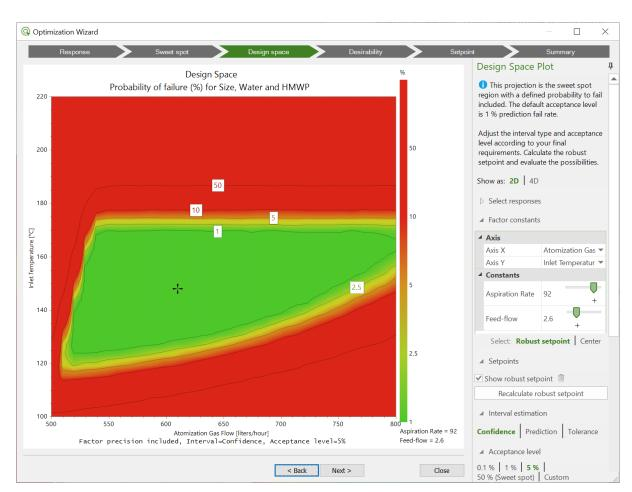
- Automatically sets factor constants to display the largest sweet spot.
- When no sweet spot can be found there is guidance in pane to the right
- Sweet spot plot is shown when there is at least one response with Condition: Required





#### Optimization Wizard - Design Space

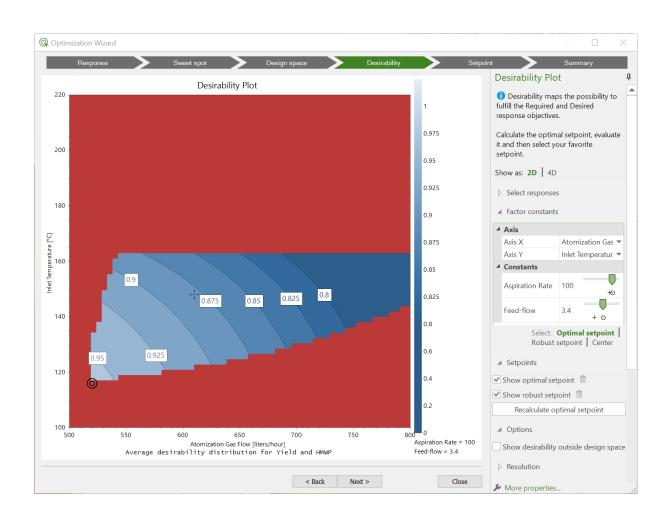
- Calculate the robust setpoint by clicking "Find robust setpoint"
  - Factor constants are adjusted to match the robust setpoint
  - Robust setpoint is marked by a crosshair
- Customization of Design Space calculation in the properties pane
- When no Design Space can be found, there is guidance in pane to the right





#### Optimization Wizard - Desirability

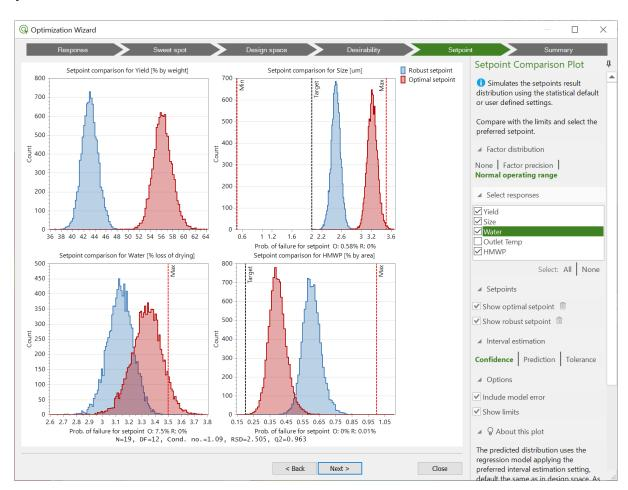
- Interpreting the desirability plot:
  - Red: outside the design space
  - Blue: inside the design space. The lighter the blue, the higher the desirability
- Calculate optimal setpoint by clicking "find optimal setpoint"
  - Factor constants are adjusted to match the optimal setpoint
  - Optimal setpoint is marked by a double circle
- When Design Space can be found there is guidance in pane to the right





#### Optimization Wizard - Setpoint Comparison

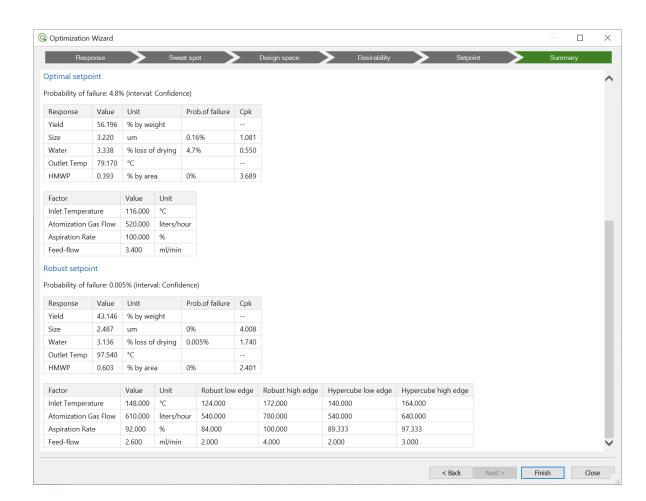
- Compare predicted distribution of selected responses. Can be used to simulate future process output
- Probability of failure for each setpoint for all responses with Objective: Required
- Adjustment of factor distribution and Interval estimation settings in pane



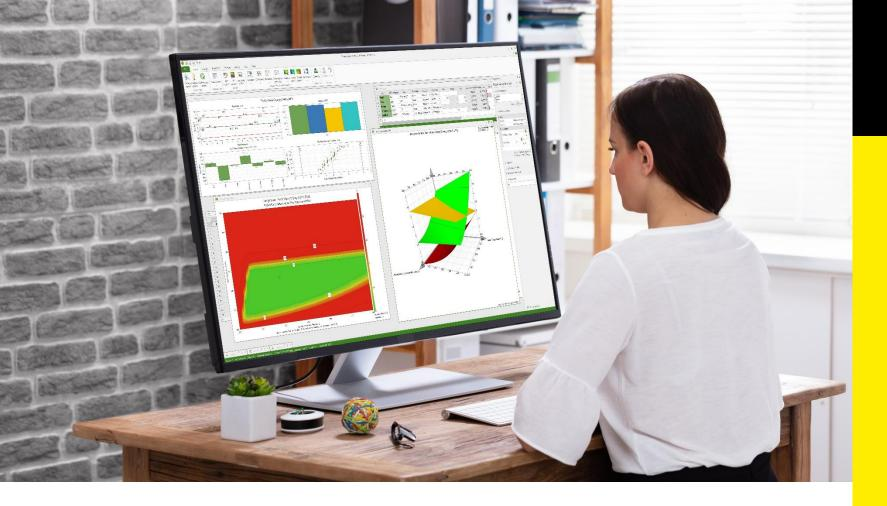


#### Optimization Wizard – Summary

- Summary of setpoint characteristics
  - Response settings
  - Factor settings
  - Predicted value
  - Probability of failure
  - Cpk
- Proven acceptable range calculated for robust setpoint
  - Robust low/high edge
  - Hypercube low/high edge







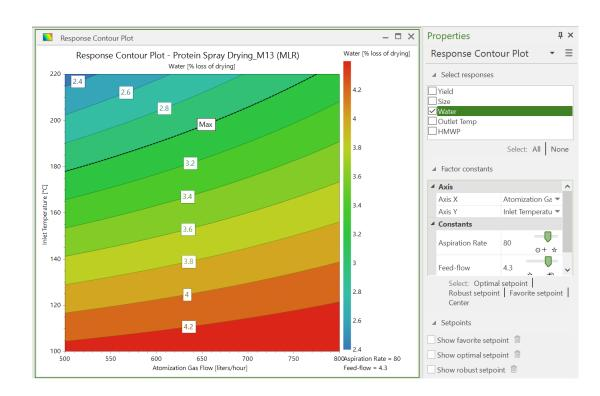
Simplifying Progress

MODDE® 13 What's New Features and improvements



#### User interface

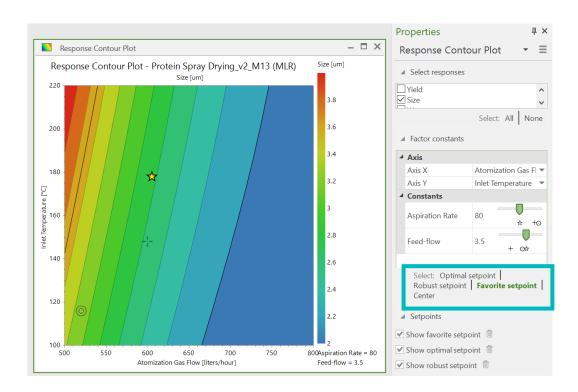
- Properties pane with most used settings
  - Available for all plots
  - Select or deselect responses
  - Select axis and adjust constants with slide bars
  - In the pane to the right specific settings for current contour plot are available and alternatives can be activated by point and click.
- Right-click any contour plot and select "... from this" creates the selected sweet spot, contour plot desirability or designs space plot with the current axis and constant settings





#### Robust, Optimal, and Favorite setpoint

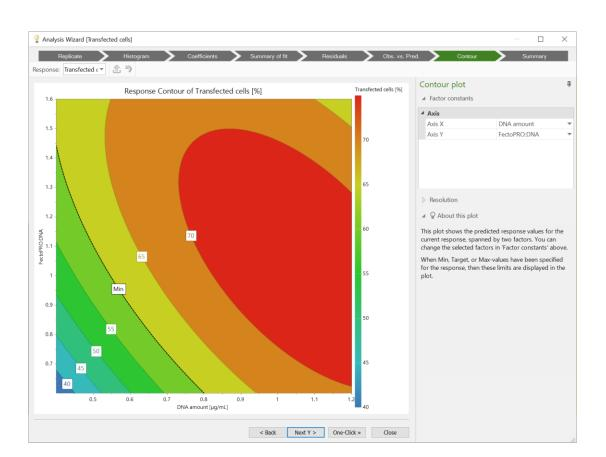
- MODDE 13 have three different setpoints and can be used in all types of 2D and 4D contour like plots
- Robust setpoint crosshair symbol
  - Calculate in design space plot and optimizer
- Optimal setpoint double circle symbol
  - Calculate in Desirability plot or optimizer
- Favorite setpoint identified by star symbol
  - Position your own favorite setpoint (right-click, Set as favorite setpoint)
- Select factor constants by clicking a setpoint in Properties pane
- When factor constants don't match exactly for a setpoint, that setpoint symbol is displayed in grey





#### Analysis Wizard

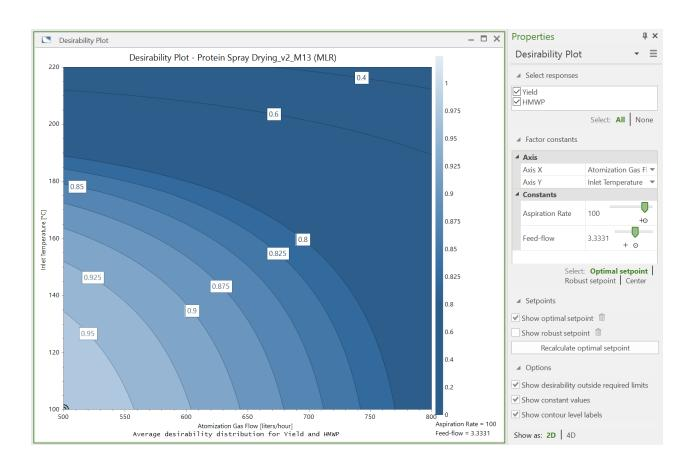
- Contour plot added in Analysis wizard for model verification by user
  - Is the model in line with expectations and prior knowledge
- 4D contour plot available, customize axis in properties pane
- Improved Interaction test-dialog for reduced factorial designs





#### Desirability plot

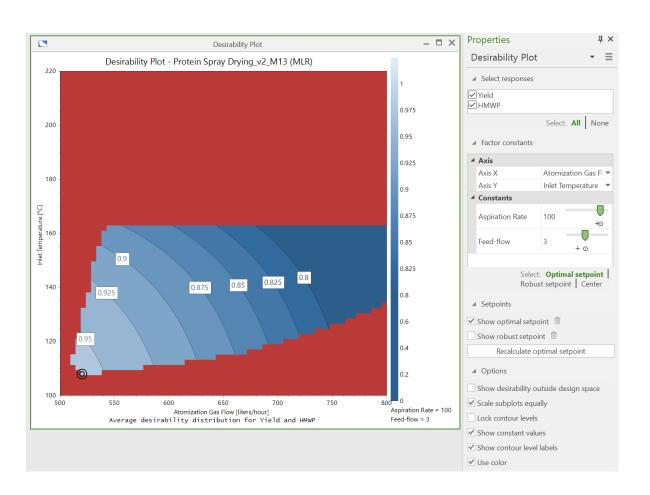
- The desirability plot shows how well the response objectives are fulfilled
  - Visualization of desirability function
  - Used to identify optimal setpoint
  - Find optimal setpoint function in properties pane to identify the optimal setpoint
  - Combine with sweet spot or design space if there are required responses





#### Optimization within design space

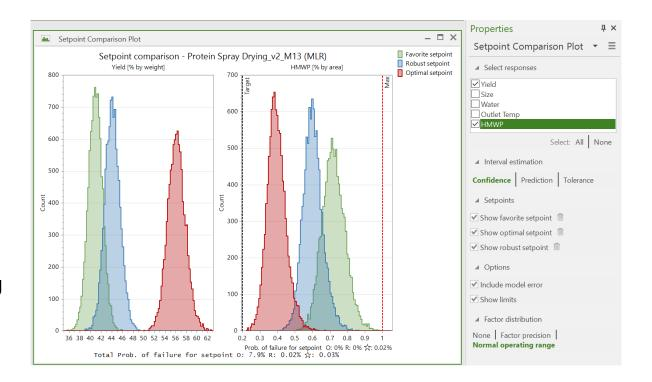
- When a design space has been calculated, MODDE can identify the factor combination with best desirability within the design space.
  - Optimal setpoint marked with double circle symbol
  - For full view of desirability, select "Show desirability outside of design space" in Properties pane
  - Create desirability plot from a sweet spot plot to calculate optimal point inside the Sweet spot.





#### Setpoint comparison

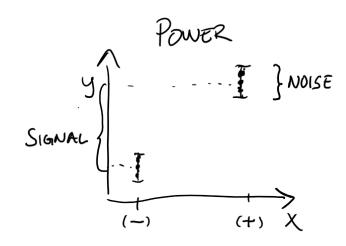
- Setpoint comparison plot shows histograms of simulated process output profiles based on factor distribution
- All defined setpoints can be visualized in the same histogram
- Total Probability of failure and individually for each response are shown for selected setpoints
- Normal Operating Range, NOR, is the default setting

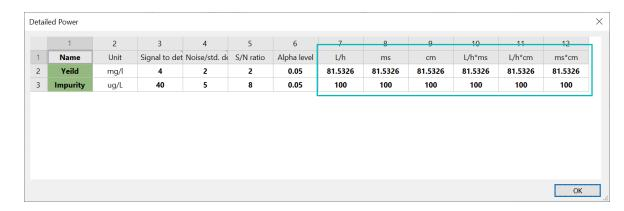




## Design Wizard - Detailed power in Responses definition dialog

- A measure of the DOE's ability to detect an effect as significant. Power is expressed in percent.
- Ratio between expected noise and the size of the effect to be detected
- Intended mainly for screening investigations
- MODDE provides detailed power for each factor per response







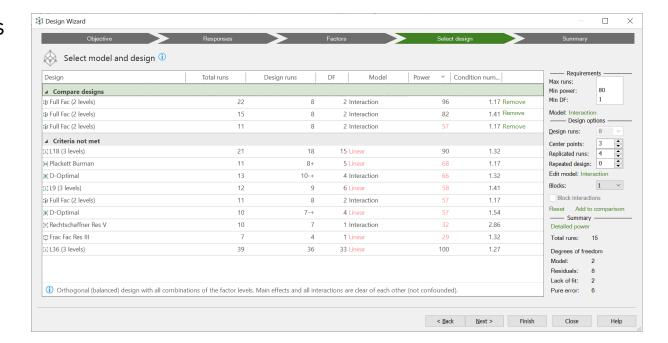
#### Design Wizard - Optimal selection of replicated design points

- Repeated design creates a DOE with 22 experiments and Power of 96
- Add replicated design runs to create a DOE with sufficient power

Original design (full fac) : 11 runs Power 57

■ Repeated design : 22 runs Power 92

Optimal replicated runs :15 runs Power 82





# Thank You for Your Interest in MODDE® 13

Don't forget to check out the instructional videos in Sartorius Data Analytics YouTube channel

