

# Data Analytics User Meeting Workshop Agenda

Monday, September 19

Start	End	Topic	Instructors
1:00 p.m.	4:00 p.m.	<b>Scale-up   Scale-down in MODDE® and SIMCA®</b> You will learn how to use MODDE® and SIMCA® to support scale-up, scale-down, and tech-transfer activities. Most of the time will be spent hands-on in MODDE® and SIMCA®, allowing you to become familiar with functionality and applications to industry problems. Topics will include scale-up and tech transfer strategies for new and legacy processes, scaling parameter sensitivity analysis and optimization, and scale-down model qualification.	Bryan Looze and Ben Reed

Tuesday, September 20

Start	End	Topic	Instructors
9:00 a.m.	12:00 p.m.	<b>Python in SIMCA®</b> You will learn how to use the Python module in SIMCA®. Most of the time will be spent hands-on in SIMCA®, allowing you to become familiar with the process of adding and deploying a Python script from start to finish. Topics will include what functions are supported via scripting, how scripting can be used to enhanced model performance, and use cases.	Amin Ganjian and Greg Casee
1:00 p.m.	4:00 p.m.	<b>Spectroscopy Modeling in SIMCA®</b> You will learn how to use SIMCA® to model spectroscopy data. Most of the time will be spent hands-on in SIMCA®, allowing you to become familiar with the process of spectroscopy modeling from start to finish. Topics will include spectral analysis, transformations, and preprocessing, and PCA   PLS modeling.	Art Hamfeldt and Michael Roberto