

Change log

Date	Prepared by	Description
2019-05-29	Andreas Norén	List of fixed bugs with reference to knowledge base article

Contents

Change log	1
1 Bugs fixed in SIMCA 16.0.1.....	2



1 Bugs fixed in SIMCA 16.0.1

The table below list the fixed bugs in SIMCA 16.0.1.

ID	Title	State	System Info
25119	Incorrect predictionset for batch model with excluded observations in BEM	Closed	https://umetrics.com/kb/batch-level-predictionset-including-workset-batch-excluded-observation-may-be-incorrect
25069	Predictionset not stable	Closed	https://umetrics.com/kb/batch-level-predictionset-including-workset-batch-excluded-observation-may-be-incorrect
25047	Newsfeed makes SIMCA crash	Closed	https://umetrics.com/kb/simca-may-crash-due-firewall-settings

