SVILCIEVS

Validation report Active Dashboard 2.0.3

2020-12-03 16:21

	Role	Name	Date
Issued by:	Software quality	Lisa Gabrielsson	2020-11-25
Revised by:	Program manager	Therese Ringvall	2020-11-30
	Product manager	Jonas Elfving	2020-12-03
Approved by:	Head of Development	Annika Finck	2020-12-03
	Head of Quality	Andreas Norén	2020-12-03

Content

1	Introduction	.2
2	Validation report summary	.2
2	.1 Validation package content	2
3	Validation task results	.2
	.1 Graphical comparison	2
4	Verification of installed software	.2
5	Source code	.2
6	Routines	.2
7	Bug handling	.2
8	Validation conclusion	.2

SVILCIEVS

1 Introduction

The purpose of the **Validation report** is to summarize and document the found differences that require corrective actions from the validation activities performed.

The scope of the validation tasks performed are described in paragraph 6.4 in the Validation plan.

This patch validation complements the full validation of Active Dashboard 2 (version 2.0.0.3697).

Approving this document includes approval of all subdocuments and results referred to in this document.

2 Validation report summary

The purpose of the **Validation report** is to summarize the validation activities performed, listed in the Validation plan, and document the found differences that need corrective actions.

The graphical validation of Active Dashboard 2.0.3.22146 was done versus Active Dashboard 2.0.0.3697.

2.1 Validation package content

The validation package includes files and folders as follows:

- Active Dashboard 2.0.3.22146 validation documentation pdf, a compilation of validation documents including this document, the Validation report.
- Bugs folder Lists details for the bugs referenced in the validation package, if any.
- Graphical validation folder Documents containing the compared graphs, lists and tables.

3 Validation task results

3.1 Graphical comparison

The graphical validation was done using the PI Data Collector and the SIMCA-online 16.1 Data Collector, and the results compared to results from Active Dashboard 2.0.0.3697.

No differences that require corrective actions were found in the graphical comparison.

4 Verification of installed software

To verify that your license of the software has been correctly installed follow the instruction here:

- 1. In Active Dashboard, click "I" (About) and under the header about Umetrics Active Dashboard verify that the version is Active Dashboard 2.0.3.22146.
- 2. Follow the instructions in Active Dashboard Installation Guide and execute the section Test your installation.

5 Source code

All source code for the final version of a full release is transferred to electronic media and kept both at the Umeå office as well as in the safe of a local bank.

6 Routines

The relevant routines are stored in TFS in the QualityManual and QualityManagementSystem folders.

7 Bug handling

Work items describing bugs found during development are stored electronically in the database TFS. Bugs that require a corrective action are listed in the tables in paragraph 3.

8 Validation conclusion

All bugs found between the release of Active Dashboard 2.0.0.3697 and the release of Active Dashboard 2.0.3.22146 that remain not fixed were considered unimportant and therefore not fixed.



SVILCIEVS

All differences that require a corrective action are listed under paragraph 3, and the WIs referenced to are stored in TFS and available in the Bugs-folder.

None of the found differences are serious. The used routines together with the validation ensure that Active Dashboard 2.0.3.22146 gives correct results and is reliable.