

Risk assessment for software updates Cubis® MCA

Table of content

General Information	2
Firmware package 09-03-01.17.04.....	3
Firmware package 09-03-01.18.12.....	4
QAPP-Center 09-06-03.00.00.....	5
QAPP-Center 09-06-03.00.03.....	6
Notes on the software update procedure.....	7

General Information

Overview

In case that a new software version is released a risk assessment for re-qualification is executed. The objective is to evaluate all software changes regarding the question, whether they require a repetition of the qualification of this instrument which ran with a previous software version.

This document provides the information and recommendation for all software updates for Cubis MCA.

The software for MCA is delivered in two packages: the "Firmware package" and the "QAPP-Center". This document covers information for both.

Scope

This document is valid for all Cubis® Models MCA.

Risk assessment

All software changes (bug fixes, improved functionality and new functionality) implemented in the new version will be classified as following:

Classification	Algorithm changed	Effect on result	Example	Re-qualification
major change	yes	yes	The measurement algorithm was changed, so that the numerical results could be affected.	recommended
minor change	yes	no	Improvement of measurement parameters and reliability	Whether a re-qualification is recommended depends on the current users application; usually not necessary
announcement	no	no	Adding a new language.	not recommended

Firmware package 09-03-01.17.04

Release date: 2019-03-08

Summary:

This is the first public firmware release.

Firmware package 09-03-01.18.12

Release date: 2019-04-05

Summary of changes from previous version:

This update contains fixes, stability enhancements and no functional extensions.

List of changes from previous version:

Sartorius ID	Brief Change Description	Comment	Classification
Cubis-14003	Fix: sporadic display issues	improved initialization routines	minor
Cubis-15226	Fix: module names for audit trail	improve readability of audit trail	announcement
Cubis-15171	Fix: no filter set in audit trail	shows all entries in audit trail with no filter	announcement
Cubis-15176	Fix: electronic signature	added missing signature with PC-Direct	minor
Cubis-14737	Fix: return from sleep under LOW	prevents lock of balance	minor
Cubis-15194	Fix: sort order of audit trail	corrects sort order on website	announcement
Cubis-15184	Fix: translation of search field	missing translation added	announcement
Cubis-15106	Fix: backup MD5 file	MD5 was unreadable in some cases	minor
Cubis-15383	Fix: missing translation	added translations to several screens	announcement
Cubis-15557	Fix: sporadic ERR235 after restart	internal communication improved	minor
Cubis-15838	Fix: isoCAL menu change needs reboot	changes now effective immediately	minor
Cubis-15840	Fix: sensitivity of motion sensors	new default setting: LOW	announcement
Cubis-15845	Fix: missing translation in weighing	missing translation added for identifiers	announcement

Recommendations:

The update is recommended.

A re-qualification is not recommended for the update from the previous version.

QAPP-Center 09-06-03.00.00

Release date: 2019-03-07

Summary of changes from previous version:

This is the initial public QAPP-Center release.

QAPP-Center 09-06-03.00.03

Release date: 2019-03-07

Summary of changes from previous version:

This update contains fixes, stability enhancements and no functional extensions.

List of changes from previous version:

Sartorius ID	Brief Change Description	Comment	Classification
Cubis-15235	Improve: QAPP-Library and initialization of QAPPs	Faster startup of QApps	minor
Cubis-15145	Fix: Backweigher	Rounding for loss in % improved	minor
Cubis-15278	Fix: Flexible Formulation	App workflow with manual takeover	minor

Recommendations:

The update is optional.

A re-qualification is not recommended for the update from the previous version.

Notes on the software update procedure

It is recommended to perform the software update in the following order:

1. Make a backup of you device data on a USB-stick before starting the update procedure.
2. Make sure that the power supply is not interrupted during the update.
3. First, install the firmware package update.
4. Second, install the QAPP-Center update.
5. Finally, create new tasks with the new QAPPs as needed.

Note: Tasks remain unchanged by the update and always use the QAPP version when the task originally created.