

My Instruments

Release Notes

1 Introduction

This document describes the implemented changes, improvements, and corrected issues in the released versions of My Instruments.

2 Release 1.2.0

Improvements

- Color coding of the state of the instruments
- Update of chart for Chromatogram

Corrected defect

- When refreshing the main page, all the system states turn orange for one second.
- Connectivity symbol and state information mismatch.
- Utilization popup close button not working properly on smaller screen sizes.
- Unable to click **Close** when My Instruments is in tablet mode.
- Status of the systems never changes to **Unknown** when disconnects occur.

Known issues

- Curve data from previous run is still presented even though the run has ended.
- Connectivity symbol constantly red (not connected) in My Instruments . This can be caused by a faulty component in the instrument. Removing the faulty component might solve this issue.
- Browsing not secure icon appears on Android™ 10 and 11 even if the certificate is installed and the browsing is secured.

- My Instruments server is not always automatically started after the restart of the lab computers.
- The search field clears spaces only after clicking the **Clear** button.
- On the dashboard, the Utilization icon for a system in less detailed view seems clickable even though it is not.
- Some parts of the software may not always fit in the screen of smaller screen size devices, such as smart phones.
- When viewing the systems in thumbnail view, there can be big gaps between systems.
- UNICORN™ certificate installer creates certificates with expiry date not based on the root certificate expiry date.
- Method information pops up eventhough no method run is performed.
- Wrong color RED for Warning when it should be ORANGE.
- Run data is enabled even though no run is in progress.

3 Release 1.1.2

Improvements

- Added image of ÄKTA™ go.
- Installation report files are updated to reflect actual status of signing.
- The version number of My Instruments is visible on the logon page and on the **About** page.

Corrected defects

My Instruments extensions successfully start on the Instrument Server even if UNICORN online is installed in the same computer.

4 Release 1.1.1

Improvements

- All required server side certificates are automatically generated and installed during installation.
- The **Control** button is now called **UNICORN**.

Known issues

Upgrade from an older version is not supported.

5 Release 1.1.0

Improvements

- A new field named **Remote Control** is added in the **System settings** tab. This field is used to access a web-based system control software using an access URL.
- Due to security hardening, now only encrypted communication is possible.
- Removed all occurrences of **... Preview** in the **Extensions** description.
- Now, the **Curve** extension refreshes itself after every 30 seconds without any user interactions.
- A new function named **Control** is added. The user requires elevated privilege to access the **Control** function.

Corrected defects

- Corrected an issue in the **System settings** tab.
- Corrected the issue regarding the presence of the REST attributes ("/app#/dashboard") in the URL when a login session has become invalid.
- Corrected a connection issue between the Dashboard Extension and the My Instruments Gateway at the startup. Now, the extension retries to connect until it succeeds.
- Changed the internal boot sequence for the adapters to avoid time-outs. This resolved the intermittent startup problem with My Instruments Gateway.

Known issues

An upgrade of My Instruments will overwrite any existing user in the **DSADashboard.dll.config** and **CADashboard.dll.config** files. It is recommend to save a backup of these files and if necessary, restore the users after the upgrade. The files are located here:

- For CA,C:\Program Files (x86)\GE Healthcare\My Instruments\Bin\Adapters\CADashboard
- For DSA,C:\Program Files (x86)\GE Healthcare\My Instruments\Bin\Adapters\DSADashboard

6 Release 1.0.2

6.1 Corrected Defects

- Some unessential prerequisites (e.g., MSXML4) have been removed from the installer.

- An issue with the End-user license agreement displaying unsupported characters is fixed.

Page intentionally left blank



cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

ÅKTA and UNICORN are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Windows, Microsoft .NET, and Windows Server are registered trademarks of Microsoft Corporation.

All other third-party trademarks are the property of their respective owners.

© 2020-2021 Cytiva

UNICORN © 2020-2021 Cytiva

Any use of UNICORN is subject to Cytiva Standard Software End-User License Agreement for Life Sciences Software Products. A copy of this Standard Software End-User License Agreement is available on request.

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

29326793 AE V:4 02/2021