

UNICORN™ software trending tool

Release Notes and Installation Instructions

1 Introduction

This document describes the main features and known issues in the released version of UNICORN™ software trending tool.

The document also describes how to install UNICORN software trending tool.

2 Release 2.1

Features

- UNICORN software trending tool 2.1 is compatible with UNICORN 7.10 and later versions. It requires a software license (the same license as for UNICORN software trending tool 2.0).
- UNICORN software trending tool 2.1 is equipped with new features, such as:
 - displaying deviations (alarms and manual interactions) directly in the trend
 - aligning UV peak area to certain fractions
 - adding certain peak parameters (peak height and asymmetry at 10%)

Some minor improvements and corrections of defects have also been implemented.

- Trends saved in earlier trending tool versions can be opened in the new trending tool version. If trends are saved in the new trending tool version, the trends cannot be opened in an earlier trending tool version.
- Results created with older UNICORN versions can be opened with UNICORN software trending tool 2.1 if the results are created with a method that is divided into different phases.

Known limitations

- Exiting UNICORN while a saved trend is open in edit mode causes the trend to become locked. To unlock a locked result:

1. Go to the **Administration** module.
 2. Select **Database Management**.
 3. Select the **Release Objects** tab.
- After repeatedly moving the legends in **Quick View**, UNICORN **Evaluation** sometimes crashes. Close and restart UNICORN.
 - After saving a snapshot in **Quick View** as an image and then opening **Quick View** in a new trend, UNICORN **Evaluation** sometimes crashes. Close and restart UNICORN.

3 Release 2.0

Features

- UNICORN software trending tool 2.0 is compatible with UNICORN 7.7 and requires a software license.
- UNICORN software trending tool 2.0 is equipped with new features, such as:
 - saving trends
 - printing trends
 - adding legends
 - result **Quick View** functionality
- Results created with older UNICORN versions can be trended as long as they are created with a method divided into different phases.

Known limitations

- Exiting UNICORN while a saved trend is open in edit mode will cause the trend to become locked. To unlock a locked result:
 1. Go to the **Administration** module.
 2. Select **Database Management**.
 3. Select the **Release Objects** tab.

4 Release 1.0

Features

- UNICORN software trending tool is a new extension to the UNICORN **Evaluation** module that is used to trend the UV peak area and the maximum pressure of multiple results, for instance, from protein purifications performed using cycling on Fibro devices.
- UNICORN software trending tool 1.0 is only compatible with UNICORN 7.6. It is available as an additional tab in the **Evaluation** module, not in evaluation classic.

- Results created with older UNICORN versions can be trended as long as they were created with a method divided into different phases. Results created with predefined methods on ÄKTA™ avant, ÄKTA go™, or ÄKTA pure™ with UNICORN 6.3 or later contain phases.

Known limitations

- The **Trending Tool** tab disappears if the **Evaluation** module is closed using the  in the upper right corner while a single result is open. Open a single result and close only that result to get the **Trending Tool** tab back.
- Right-clicking a result in the **Results** browser and then selecting many results induces very slow response times. Log out and log in again into UNICORN to return the response time to normal.

5 Installation Instructions

Install UNICORN software trending tool

1. Open UNICORN.
2. In the **Administration** module, click **Extension Management**.
3. Click **Install** and locate the UNICORN software trending tool extension software file.

Activate UNICORN software trending tool

On the UNICORN computer, where the **Trending Tool** is to be used:

1. Open UNICORN.
2. In the **Administration** module, click **Extension Management**.
3. Tick the box beside the **Trending Tool** extension, and click the **OK** button.
4. Log out and log in again into UNICORN.
5. Go to the **Evaluation** module in UNICORN. If you see evaluation classic, switch to the **Evaluation** module. After successful installation, you will see a tab called **Trending Tool** next to the **Results** tab.

6 Privacy and Security Capabilities

UNICORN includes a broad assortment of capabilities to enable privacy and security. For more details about this release, please refer to UNICORN 7.10 Privacy and Security Manual.



cytiva.com/unicorn

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

ÄKTA, ÄKTA go, ÄKTA pure, and UNICORN are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

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

© 2021–2023 Cytiva

UNICORN © 2021–2023 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.

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

29697532 AC V:3 07/2023