

My Instruments 1.2 Service Pack 3 Release Notes

1 Introduction

This document describes the main functionality of the released version of My Instruments 1.2 Service Pack 3 (SP3), which has been verified using UNICORN™ versions 7.6 to 7.10.1 and My Instruments 1.2.

SP3 is required when using My Instruments 1.2 in combination with UNICORN version 7.9 and later versions of UNICORN.

SP3 is not required when using My Instruments 1.2 in combination with UNICORN version 7.6-7.8.

Note: *SP3 consists of new versions of two separate extensions that need separate downloads from the Cytiva Marketplace. The extensions are referred to as **My Instruments 1.2 Service Pack 3 – D** (containing the **Dashboard** extension) and **My Instruments 1.2 Service Pack 3 – DC** (containing the **DashboardConfiguration** extension).*

For installation instructions, refer to the *My Instruments Installation Guide*.

Note: *My Instruments 1.2 Service Pack 3 replaces and contains the functionality of My Instruments 1.2 Service Pack 2 and Service Pack 1.*

Prerequisite

UNICORN (version 7.6 or a later version) and the My Instruments 1.2 package need to be installed before downloading and installing SP3.

For detailed instructions, refer to the *My Instruments Installation Guide*.

2 Corrected defects

- Fixes an issue with certificate validation of included extension files.

- Contains new versions of the **Dashboard** extension (version X) and the **Dashboard-Configuration** extension (version X), which replace the previous versions of the extensions.

Page intentionally left blank

**Give feedback on this document**

Visit cytiva.com/techdocfeedback or scan the QR code.



cytiva.com

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

UNICORN is a trademark 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.

© 2020–2023 Cytiva

UNICORN © 2020–2023 Cytiva

Any use of software may be subject to one or more end user license agreements, a copy of, or notice of which, are available on request.

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

29743728 AA V:1 10/2023