

# Cytiva data bridge

## Release Notes

### Introduction

This document describes the major implemented changes, improvements, and corrected issues in released versions of Cytiva data bridge.

### Release 1.0

#### Functionality

- eCoA and eCoA + lot genealogy data for Cytiva cell culture media products shared through a secure Snowflake database.
- Dedicated Data-warehouse to store data.
- Role-based access controls to limit access to the Snowflake database with separate Cytiva administrator roles.
- Option for customers to integrate with their own environments to transfer data as required.
- Multiple options to on-board new customers.
- Possibility to access audit logs via Snowflake.
- Storing eData in XML document format and relational database format in Snowflake with secure download option.
- Query data bridge relational database.
- Retrieve data bridge documents linked to database records.
- Segregate accessible eData between customers.
- Manage access to data bridge on product and eData type level.

- Extract eData from systems of record, transform to target format, load into the system.
- Access to customer Support Team for inquiries.
- Access to data bridge documentation including FAQ, help resources, user manuals, etc.

Page intentionally left blank



**Give feedback on this document**

Visit [cytiva.com/techdocfeedback](https://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.

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

© 2024 Cytiva

Data bridge © 2024 Cytiva

Any use of Data bridge is subject to Cytiva Terms of Service for Cloud Products.

For local office contact information, visit [cytiva.com/contact](https://cytiva.com/contact)

29747241 AA V:1 04/2024