



Cytiva data bridge

User Manual

Table of Contents

1	Introduction	3
2	Log in to Snowflake	4
3	View databases and tables in Snowflake	6
4	Run a query in Snowflake	7
5	Example SQL queries	9
6	Dashboards	10

1 Introduction

This user manual describes the features available in Cytiva data bridge and provides instructions for managing, utilising, and accessing data.

Data bridge delivers an infrastructure where data can be shared with customers through a cloud database solution. The initial release will focus on raw materials data, although the platform is designed to be scalable to incorporate additional data types in the future.

Scope

The initial release of data bridge uses the Snowflake user interface. The scope of this user manual is to help the user get started with the Snowflake interface. For detailed information on using Snowflake, see their documentation: <https://docs.snowflake.com/>.

Definitions/abbreviations

Abbreviation	Definition
SQL	Structured Query Language. For more information, see the Snowflake SQL command reference at https://docs.snowflake.com/en/sql-reference/constructs .

2 Log in to Snowflake

Step	Action
------	--------

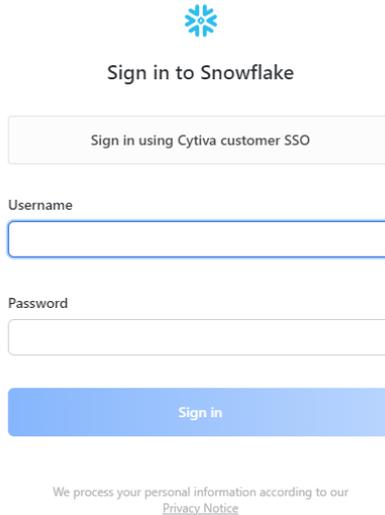
1	Log in at https://www.cytiva.com
---	---

The screenshot shows the Cytiva website's login interface. At the top, there is a navigation bar with the Cytiva logo on the left, a search bar in the center, and links for 'Quotes' and 'Cart' on the right. Below the navigation bar, there are links for 'Products', 'Applications', 'Service & Support', 'Resources', and 'Our Company'. A 'Track Order' link and a highlighted 'Register/Log in' button are also present. The main content area is titled 'Log In' and contains a form with the following elements:

- A 'Username' label followed by a text input field.
- A 'Password' label followed by a text input field and a 'Show' icon.
- A green 'Log In' button.
- A link for 'Forgot Password?'.
- A text prompt 'Not registered with us yet?' followed by a 'Register Now' button.

Step Action

- 2 Copy-paste the URL for your given Snowflake account into the same browser. Example URL:
customername_databridge.snowflakecomputing.com



The screenshot shows the Snowflake sign-in interface. At the top center is the Snowflake logo, a blue snowflake icon. Below it is the text "Sign in to Snowflake". Underneath is a light gray button labeled "Sign in using Cytiva customer SSO". Below this button are two input fields: "Username" and "Password". The "Username" field is currently selected with a blue border. At the bottom of the form is a blue button labeled "Sign in". Below the button, there is a small line of text: "We process your personal information according to our [Privacy Notice](#)".

Result:

This will automatically authenticate using Cytiva Single sign-on (SSO).

The process is the same if a user has access to other Snowflake accounts. Once the user has signed in with Cytiva.com, they should be able to copy-paste the Snowflake account URL into the browser and authenticate with Cytiva.com SSO.

3 View databases and tables in Snowflake

You can use Snowflake to view the eData databases, their columns and the data they contain. For more information, see the Snowflake documentation: <https://docs.snowflake.com/>.

BATCH_ID	DATA_PARTY_EMAIL	GENERATION_DATE	GENERATION_TIME	COMMENTS
20240314_003	null	2024-03-14 18:03:05	null	This product is manufactured using 30% simethicone
21	20240314_003	null	2024-03-14 18:03:05	This product is manufactured using 30% simethicone
22	20240314_003	null	2024-03-14 18:03:05	This product is manufactured using 30% simethicone

Step Action

- 1 Click **Data** → **Databases** in the left-hand menu.
- 2 Click a database to view its contents.
In the above screenshot, **EDATA_STAGING** is a table within the **EDATA_STAGING** database. This information is important when you run a query. For more information, see [Chapter 4 Run a query in Snowflake, on page 7](#).
- 3 Click **Data Preview** to see a table displaying a limited number of data rows (4) from the selected database.

Now that you know the names of the databases, their tables, their columns, and their data, you can run a query in Snowflake. See [Chapter 4 Run a query in Snowflake, on page 7](#).

4 Run a query in Snowflake

You can use Snowflake to access eData via SQL queries.



Step Action

- 1 In the left-hand menu, click **Projects** → **Worksheets**
- 2 In the top right corner, click **+** → **SQL Worksheet**

Result:

An empty SQL worksheet opens:



- 3 Click **No Database selected** to select a database.
- 4 Enter your SQL query in the text area.

Example query: `select batch_id from ecoa_stage where lot like ('%AJ307%') and mfg_date like ('%2023-08-22%') and specification = 'Satisfactory'`

Note:

For more information on SQL syntax, see the *Snowflake SQL command reference*: <https://docs.snowflake.com/en/sql-reference-commands>.

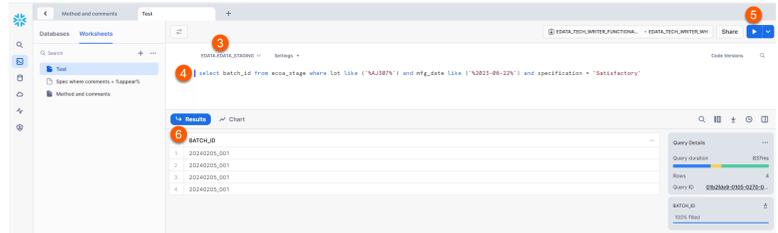
Note:

The **from** keyword must be followed by the name of a **table** in your selected **database**. In the example query, that table is **ECOA_STAGE**, found in the **EDATA_STAGING** database.

See [Chapter 3 View databases and tables in Snowflake](#), on page 6 for more information.

Step	Action
------	--------

The table name is not case sensitive when writing a query.



5 Click the arrow button in the top right corner to run the SQL query.

6 The results will be shown in a table below the SQL query. Click **Chart** to view a chart visualization of the result data.

For more information on visualizing data, see the Snowflake documentation: <https://docs.snowflake.com/en/user-guide/ui-snowsight-visualizations>.

5 Example SQL queries

These examples are given to help users get started with writing their own queries. For more information, see the Snowflake documentation: <https://docs.snowflake.com/en/guides-overview-queries> and <https://docs.snowflake.com/en/sql-reference/constructs>.

select specification **from** ecoa_stage **where** method = 'Endotoxin' **and** comments **like** ('%Cell growth%') **order by** lot_date **asc**

Note: The **select** keyword selects the data column(s) which will be visible in the results. In this case, the **specification** column.

Tip: Use **select *** to select all data columns.

Tip: To search for partial phrases within text strings, use the keyword **like**, followed by the search phrase enclosed as such: (**%search_phrase%**). In this example, we are looking for **comments** that include the phrase "Cell growth".

Tip: Use **order by column asc** to sort the data by ascending values. Use **desc** to sort by descending values.

You can also order the displayed result data by the values in a different data column. In this case, we display the **specification** data, but we order the results by the **lot_date**.

select batch_id **from** ecoa_stage **where** lot **like** ('%AJ307%') **and** mfg_date = '2023-08-22' **and** specification = 'Satisfactory'

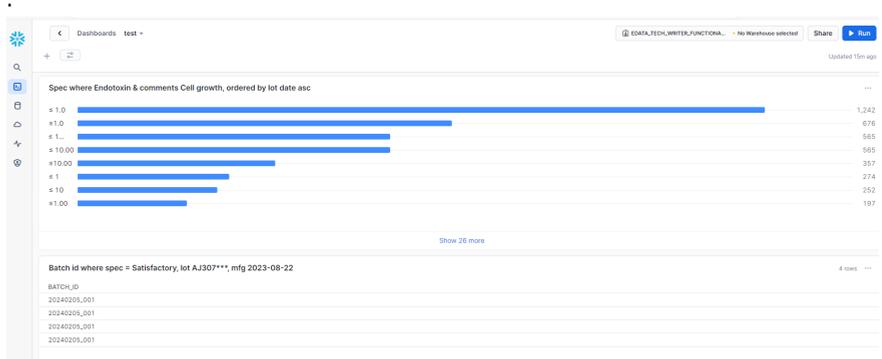
Note: In the above example, we display the batch ID for lots that fit the below criteria:

- Lot ID begins with AJ307. (The **like** keyword will look for "AJ307" anywhere in the column data, but in this example data set, we know that AJ307 only occurs at the start of the lot ID.)
- Manufactured on 2023-08-22.
- The **specification** is defined as "Satisfactory."

6 Dashboards

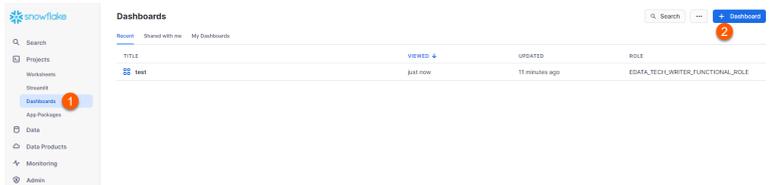
In Snowflake, you can create custom dashboards to quickly view your data queries. For more information, see the Snowflake guide on dashboards and visualizations:

- <https://docs.snowflake.com/en/user-guide/ui-snowsight-dashboards>
- <https://docs.snowflake.com/en/user-guide/ui-snowsight-visualizations>



Step Action

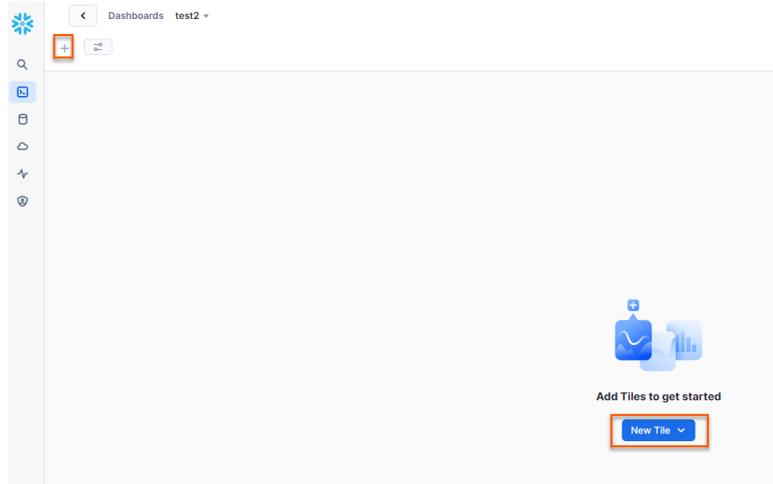
- 1 Click **Projects** → **Dashboards** in the left-hand menu.



- 2 Click the **+ Dashboard** button in the upper right corner.
- 3 In the following pop-up, name your dashboard and click **Create Dashboard**.

Step Action

- 4 Click the **+** button in the top left corner, or **New Tile**, to add an SQL query worksheet to the dashboard.





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.

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

29744780 AA V:3 05/2024