

Cloud connectivity with the NanoDrop Ultra UV-Vis Spectrophotometers and Fluorometers

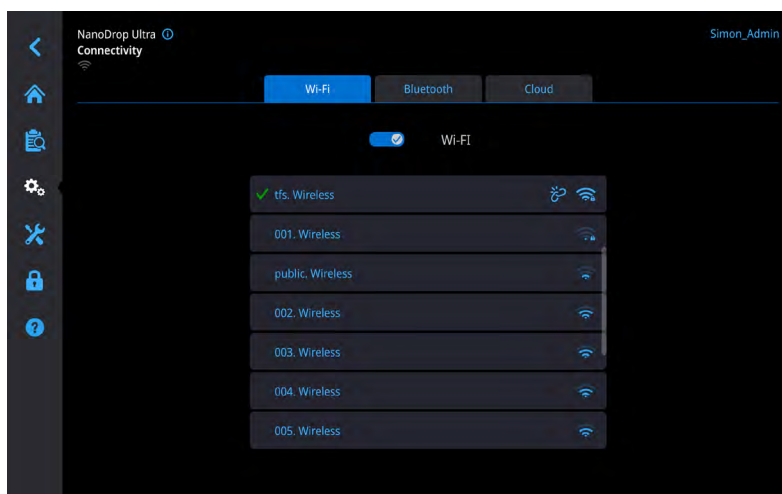
Introduction

In the modern laboratory landscape, seamless integration of analytical instruments with cloud-based data management systems is paramount for enhancing data accessibility, collaboration, and storage efficiency. The Thermo Scientific™ NanoDrop™ Ultra UV-Vis Spectrophotometers and Fluorometers not only deliver high precision in measuring nucleic acid and protein concentrations but also offer seamless data management and accessibility through integrated cloud connectivity. By leveraging cloud connectivity, researchers can ensure secure, real-time access to their spectrophotometric data from popular cloud platforms like Microsoft OneDrive, Google Drive, and Thermo Fisher Connect, thereby facilitating better data management and collaborative research efforts. This document will outline the steps required to configure the NanoDrop Ultra for cloud integration, the prerequisites for establishing a connection, and best practices for ensuring data integrity and security during the exportation process.

Prerequisites

Network connection

1. Establishing a connection between a cloud platform and the NanoDrop Ultra spectrophotometers and fluorometers first involves proper network setup to ensure both systems can communicate effectively. This can be achieved through either Wi-Fi or direct ethernet connection. The NanoDrop Ultra USB Wi-Fi and Bluetooth Dongle must be inserted if Wi-Fi is to be used for the network connection. Additionally, Wi-Fi must be enabled: open settings > **Connectivity** > enable the Wi-Fi toggle.

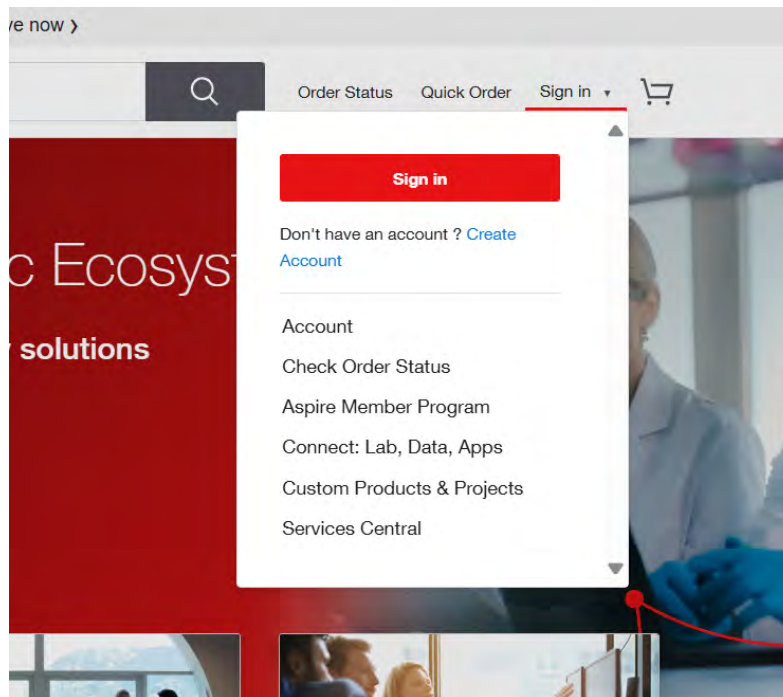


Prerequisites (continued)

Create a Thermo Fisher Scientific account

Connection to the various available cloud platforms is executed through the Thermo Fisher Cloud Connect interface. This application requires a Thermo Fisher Scientific account.

1. From an internet connected web browser, browse to www.thermofisher.com/cloud.
2. Select **Sign in** in which will display a dropdown menu.



3. Select **Create Account** and complete the additional steps to create a free account.

Create a Thermo Fisher Cloud Connect PIN

1. After signing into your Thermo Fisher Scientific account in your preferred browser, select **My Account**.
2. Select **Connect: Lab, Data, Apps** which will take you to the Thermo Fisher Connect home page.
3. Open the settings by selecting ☰ at the top left of the screen.
4. Open the **Instruments** dropdown, then select **Connected Instruments**.
5. Select **Change PIN** at the upper right of the Instrument dashboard.
6. Enter a unique PIN number and save this information as you will need it later, then select **Confirm**.

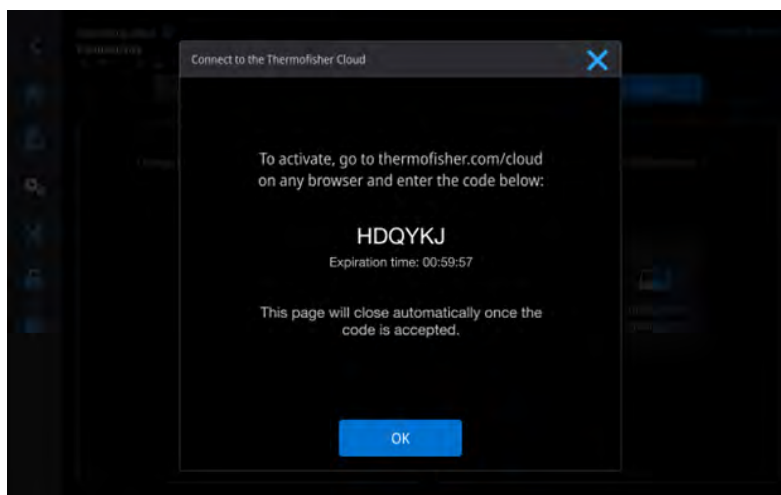
Cloud connectivity

Thermo Fisher Connect

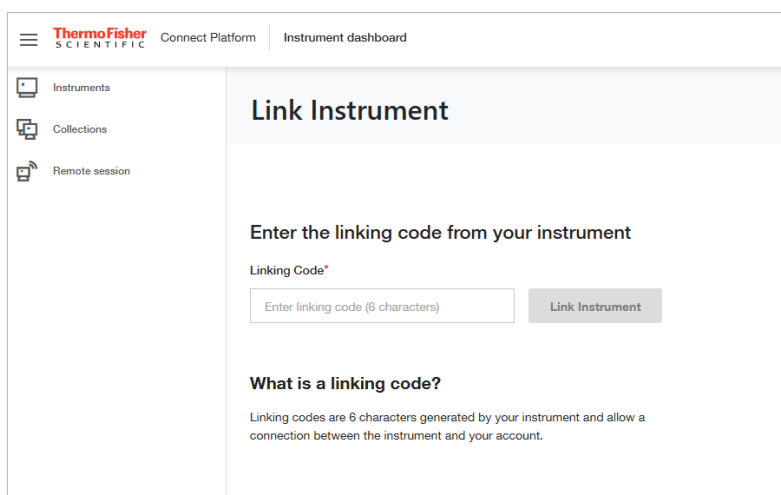
1. From the NanoDrop Ultra home screen, open the settings options by selecting ⚙️ on the navigation bar.
2. Select **Connectivity**, then open the **Cloud** tab.
3. If you are located within China, select **China** as the region; otherwise ensure **Others** is selected.

Cloud connectivity (continued)

4. Select **Dashboard** on **Internet Browser** to display a unique linking code.



5. From the Thermo Fisher Connect home page, open the settings by selecting ☰ at the top left of the screen.
6. Open the **Instruments** dropdown, then select **Connected Instruments**.
7. Select **Link instrument**, then enter the unique linking code provided by the NanoDrop Ultra.
8. Select **Link Instrument** to confirm.



9. The connected Thermo Fisher Scientific account will now appear in the **Cloud** tab within the settings on the instrument with the active status and will now appear as an export option.

Google Drive and Microsoft OneDrive

The connection to the Thermo Fisher Connect platform must be established before connecting to other cloud platforms, see the above section for more information.

1. From the Thermo Fisher Connect home page, open the settings by selecting ☰ at the top left of the screen.
2. Open the **Instruments** dropdown, then select **Connected Instruments**.
3. Use the instrument settings ⋮ button below the now connected NanoDrop Ultra and select **Storage Settings**.

Instrument dashboard

Sort by:

Select sort key
Name A-Z



Firmware Version: 1.0.0.1

OS Version: 1.0.0.3

Software Version: 1.0.0.33

Package Version: 1.0.0.21

Social Number

Online

Calendar, Star, and More icons

4. Select **Add account**.
5. Select **Add account** under the desired cloud platform.

Add account [Close]

Add one Google or OneDrive account each to move data from your Instrument_NUI2400038

Google Drive
Max. 1 account
Share your instrument files to your Google drive account.
[Add account](#)

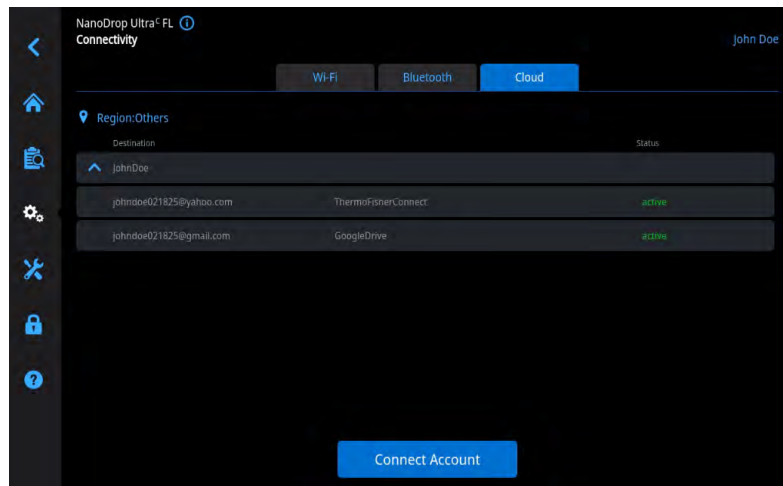
One Drive
Max. 1 account
Share your instrument files to your One drive account.
[Add account](#)

6. Grant thermofisher.com access to the cloud drive.
 - Ensure the option **See, edit, create, and delete all of your Google Drive files** is checked off for Google Drive connection; if it is unchecked the files will not be saved to the drive.

See, edit, create, and delete all of your Google Drive files. [Learn more](#)

7. In the **Cloud** tab within the instrument settings, select the down arrow followed by the PIN to view the newly added cloud drive(s). The newly added cloud drive should be listed as active and will now appear as an export option.

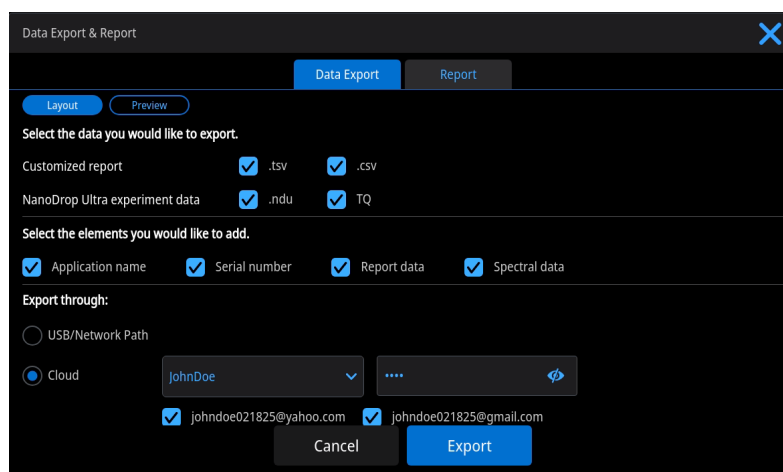
Cloud connectivity (continued)



Export to the cloud

Exportation of data can occur at the end of an experiment or by opening a previously performed experiment.

1. From the Data Export & Report window select the Data Export tab and ensure you are on the **Layout** subtab.
2. Select the desired export file type(s) by checking the box beside each option.
3. Select the desired report elements you would like to appear in the report by checking the box beside each option.
4. Select the **Cloud** radio button.
5. Select from the list of available connected Thermo Fisher Scientific accounts by using the provided dropdown menu, then entering that account's unique PIN.
6. Select the desired cloud platform(s) to export to by checking the box beside each option.
7. Select **Export**.



**Cloud connectivity
(continued)**

The data file(s) will now appear in the respective cloud drive(s).

