

# Cloud Connect Utility

## HELP

Publication Number MAN0017164

Revision A.0

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**NOTICE TO PURCHASER: DISCLAIMER OF LICENSE:** Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Life Technologies Corporation, either expressly, or by estoppel.

**Revision history:** Pub. No. MAN0017164

Revision	Date	Description
A.0	18 August 2017	Help system. Instructions for installation and use of the Cloud Connect Utility.

**Corporate entity:** Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

**Trademarks:** All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2017 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

- **About the Cloud Connect Utility ..... 4**
  - Download and Install the Cloud Connect Utility ..... 4
  - Set up the Cloud Connect Utility connection ..... 4
  - Upload files using the Cloud Connect Utility ..... 5
  - Manage your connection ..... 5
  - Quit the application ..... 6
  - Delete the connection ..... 6
  - Install updates ..... 6
- Documentation and support ..... 7**
  - Customer and technical support ..... 7
  - Limited product warranty ..... 7




# About the Cloud Connect Utility

This help guide provides instruction on the installation, initial connection set up, and use of the Thermo Fisher Scientific Cloud Connect Utility to automatically transfer files of any type to your Thermo Fisher Connect account.



Use the **Table of Contents** in the left navigation pane to find your topic of interest, or enter your search term(s) in the **Search** field above.

## Download and Install the Cloud Connect Utility


1. Sign in to Thermo Fisher Connect using your username and password. If not already registered, click **Sign Up Now**.
2. Download the utility to your computer.
  - a. Click  (Apps) to open the **AppConnect** screen.
  - b. Scroll to the **Utilities** section, then in the **Cloud Connect Utility** tile for your operating system, click **Download**. The file will automatically begin downloading.
    - ThermoFisherCCUv2.exe (*Windows*)
    - ThermoFisherCCUv2.dmg (*Mac*)
3. Double-click the .exe or .dmg file to launch the installer.
4. Enter your Thermo Fisher Connect username and password, then click **Sign In**.

## Set up the Cloud Connect Utility connection

After installing the software, you must configure the utility to link with your Thermo Fisher Connect account and designate source and destination folders for file transfers. The utility monitors the source folder for new or updated files and uploads them to the destination folder automatically.

1. Select your local folder.
  - a. Click  **Select your local folder**.
  - b. Navigate to your folder of choice, or create a new folder.
  - c. Click **Select Folder**.
2. Select your destination folder in the Thermo Fisher Connect DataConnect screen. Alternatively, create a new destination folder.
  - a. Click  **Select a DataConnect folder**.



- b. Navigate to your folder of choice, or click  **Create new folder**. Enter a folder name, then click ✓.
- c. Click **Select**, then click **Next**.





3. Enter a **Connection name**, then click **Next**.

The utility will automatically launch and begin uploading files. The connection will appear as a tile in the **PCs & smartphones** tab of your Thermo Fisher Connect account **InstrumentConnect** screen with the name you gave the connection.

## Upload files using the Cloud Connect Utility

Files placed in your source folder are automatically uploaded to your destination folder in **DataConnect** as long as you have an active connection. To reestablish your connection and upload files to your destination folder in Thermo Fisher Connect:

**Note:** On first setup the utility launches automatically and starts uploading files that are in your source folder.

- (Windows)
  - a. Double-click the desktop shortcut or click **Start ▶ All Programs ▶ Thermo Fisher Scientific ▶ Cloud Connect Utility** to launch the utility.
  - b. Transfer files for upload into your source folder.  
Upload progress is indicated in the  **Upload activity** screen.
  - c. Click  **DataConnect folder** to verify your upload.
- (Mac)
  - a. Double-click the **Cloud Connect Utility** icon in the **Dock**, on the **Launchpad**, or in your **Applications** folder.
  - b. Transfer files for upload into your source folder.  
Upload progress is indicated in the  **Upload activity** screen.
  - c. Click  **DataConnect folder** to verify your upload.

## Manage your connection


To make any change to your connection, first delete the existing connection, then set up a new connection with a new configuration.

Click the  **Settings** tab to view:

- Username and Password
- Connection name
- Source folder
- DataConnect folder
- Delete connection



## Quit the application

Click the  **Quit application** tab, then click **Yes, exit applicaiton**.

**Note:** If you quit the application, file uploads that are in progress will be cancelled.



To restart the Cloud Connect Utility:

- (Windows) Double-click the desktop shortcut or click **Start ▶ All Programs ▶ Thermo Fisher Scientific ▶ Cloud Connect Utility**.
- (Mac) Double-click the **Cloud Connect Utility** icon in the **Dock**, on the **Launchpad**, or in your **Applications** folder.

## Delete the connection


To make any change to your connection, first delete the existing connection, then set up a new connection with a new configuration.

**Note:** File uploads that are in progress will be cancelled, and the connection with Thermo Fisher Connect will be deleted.

- In **Cloud Connect Utility**, click  **Settings ▶ ↔ Delete connection**, then click **Yes, delete connection**.
- or*
- In **InstrumentConnect ▶ PCs & smartphones**, select the **Cloud Connect Utility** tile to be deleted, click  **Disconnect**, then click **Confirm**.

## Install updates

When an update to the Cloud Connect Utility is available an **Update** button will appear in the left navigation menu.

Click  **CCU update available**.

- Click **Yes** to quit and relaunch the application immediately. The update will be installed.
- Click **No** to continue without installing the update.

**Note:** The application will be updated the next time you quit and relaunch the Cloud Connect Utility.

# Documentation and support

## Customer and technical support

Visit <http://thermofisher.com/support> for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

**Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at <http://www.thermofisher.com/support>.

