

Connect your NanoDrop™ One Instrument to the Thermo Fisher Cloud


Pub. No. MAN0017041 Rev. A.0

Link your NanoDrop™ One instrument with Thermo Fisher Cloud

Linking your NanoDrop™ One instrument profile with the Thermo Fisher Cloud allows you to:

- View your instrument status from the Cloud.
- Transfer run data from the instrument to your Cloud storage.

Link an existing profile with Thermo Fisher Cloud

1. Touch  **Settings**.
2. Open the **Networking** tab.
3. Touch **Cloud Connect**.
4. Enter your username and password, then touch **Connect**.
5. Touch **OK**.

A message confirming your successfully registered your instrument is displayed.

Your instrument profile is linked with the Thermo Fisher Cloud.

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

thermofisher.com/support | thermofisher.com/askaquestion

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

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.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.