

Thermo Scientific™ DeviceLink™ Connect LAN (Ethernet) Commissioning/Setup

Commissioning of the DeviceLink Connect can be done both via Wi-Fi, as described in the Getting Started section of the user manual, or via a direct LAN (ethernet) connection between the DeviceLink Connect mote and a computer. The following guide describes the commissioning process via LAN.

Components Needed:

- (1x) DeviceLink Connect mote (provided with your DeviceLink Connect kit)
- (1x) PoE wall adapter and PoE cable (provided with your DeviceLink Connect kit)
- (2x) Ethernet cables
- (1x) Computer

Hardware Setup:



- The DeviceLink Connect needs to have an active network connection on the "POE LAN" port, either via a PoE network jack, or via the PoE power injector "IN" port (as shown above)
- The DLC port labeled "LAN" is for ethernet commissioning **only**. It is not for providing a long-term network connection to the device

Commissioning Procedure:

1. On your computer, navigate to your network settings and configure your ethernet/LAN port with the following static IP:
 - a. IP: 192.168.5.10
 - b. Subnet mask: 255.255.255.0
 - c. Gateway and DNS servers should be blank

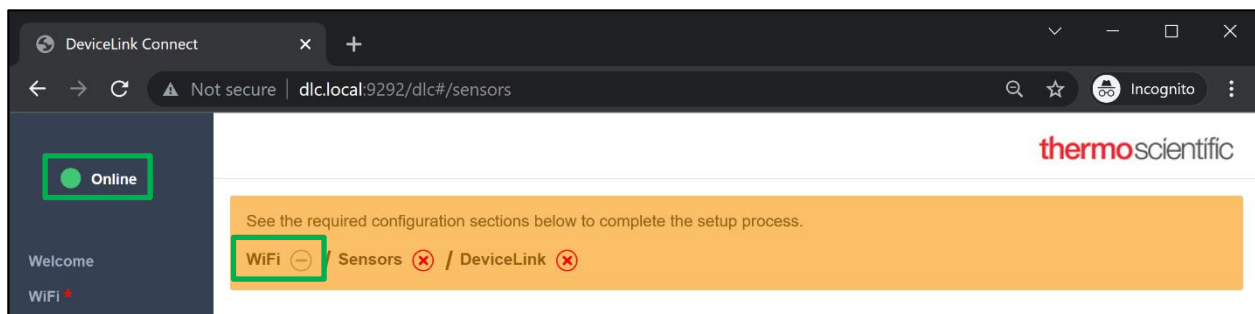
For more information on configuring a static IP, refer to the [Configuring your ethernet \(LAN\) port](#) below.

2. Open a web browser and navigate to: <http://192.168.5.5:9292/dlc>

3. You will be prompted to login. The first time you login use the following credentials:
 - a. **Username:** admin
 - b. **Password:** changeThisPassword

After logging in you will be prompted to change the password and re-login with the updated password.

4. Upon loading, the commissioner will display a status header bar at the top of the screen. If the DeviceLink Connect has an active network connection via the POE LAN port, as described above, and is able to connect to the Thermo Connect (cloud) network, "WiFi" will show a *dash* symbol. You will also see an *Online* indication in the left navigation menu. If there is no connection established, the status bar will be a *red X*.



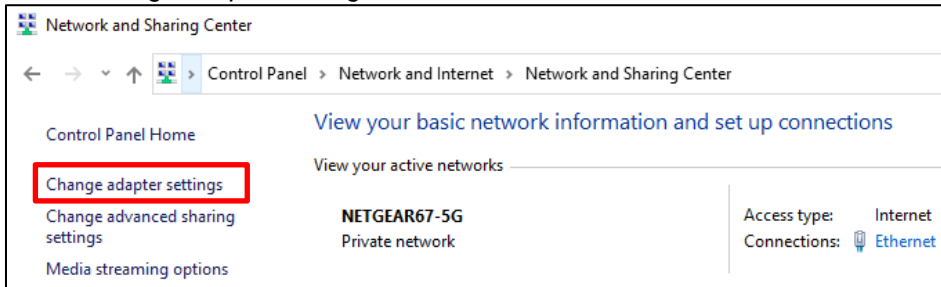
If the device is offline (no connection):

- a. Verify the cable connections are as indicated in the above hardware setup
 - b. Consult your network/IT administrator as the DeviceLink Connect may need to be whitelisted on your network. Ports may also need to be opened. Refer to the [Ports and Whitelisting](#) guide below for more information.
5. Continue the commissioning of your device per the DeviceLink Connect Quick Start Guide starting with the setup of your **Sensors**.

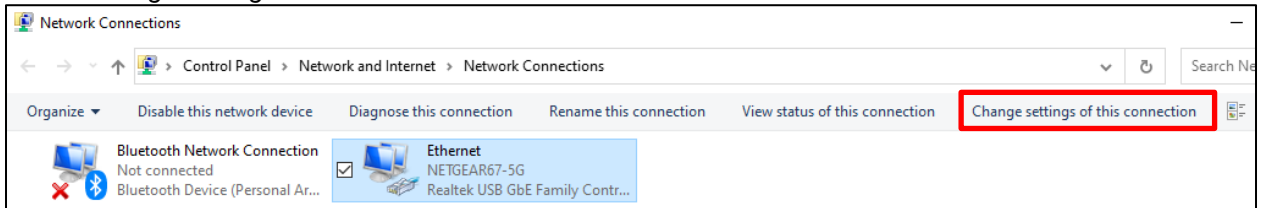
Configuring your ethernet (LAN) Port

Windows PC:

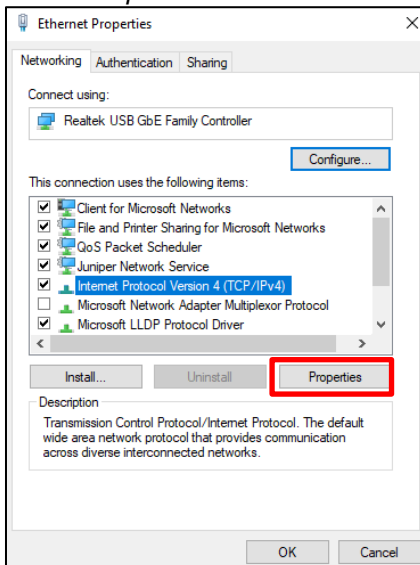
1. Navigate to the Control Panel
2. Select **Network and Sharing Center**
3. Select *Change adapter settings*



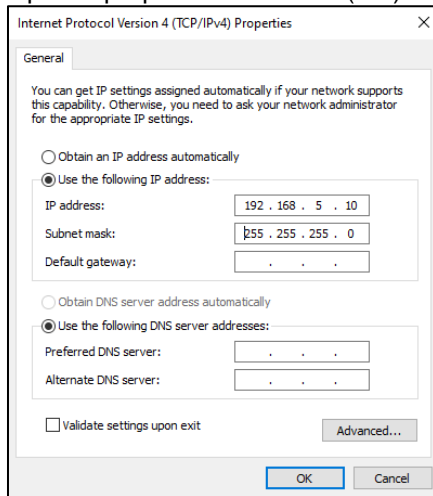
4. Select your ethernet adapter from the list
5. Select *Change settings of this connection*



6. Select the TCP/IPv4 from the connection list
7. Select *Properties*

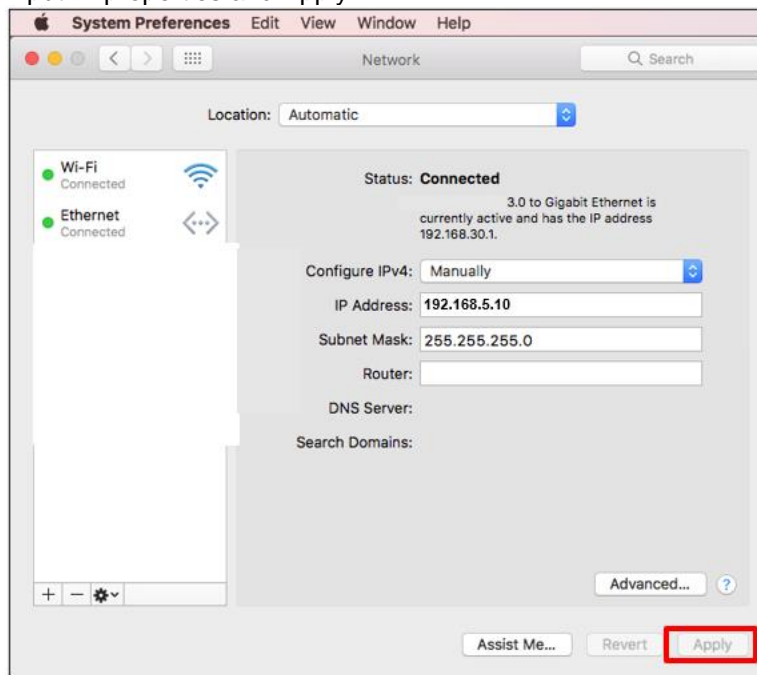


8. Input IP properties and save (OK).



MAC OS:

1. Navigate to **System Preferences** (either from your dock or by clicking the Apple menu at the top of your screen).
2. Select **Network**
3. Select your Ethernet connection from the left list menu
4. Change the Configure IPv4 to "Manually"
5. Input IP properties and Apply.



Ports and Whitelisting:

Port **80** (HTTP) or Port **443** (HTTPS) with full access to **api.traxxekg.com** must be open/configured.

Static IP addresses for api.traxxekg.com: 34.209.13.39, 35.155.85.162, or 52.27.148.178

An explicit HTTP proxy is configurable via the DeviceLink Connect Commissioner application (Cloud menu).