

# QuantStudio™ 6 Pro and 7 Pro Real-Time PCR Systems

Pub. No. MAN0018160 Rev. A.0

## Let's connect

To review our security white paper and detailed information about the features and benefits of our Connect platform, click [here](#).

**Questions?** We're here to help. Send your questions to [TFC.Device.Connectivity.Support@thermofisher.com](mailto:TFC.Device.Connectivity.Support@thermofisher.com). You must complete and return this checklist before installation and training can be scheduled.

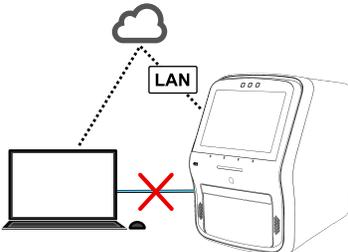
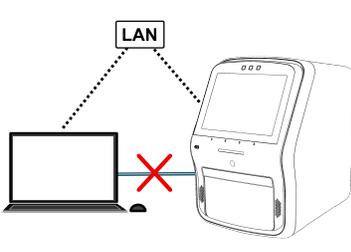
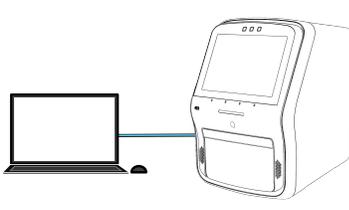
**IMPORTANT!** The instrument must be networked and connected to the internet in order to enable the Hands Free and Smart Help features.

- Step 1: Instrument administrator: sign into your Connect account and select your connection type
- Step 2: Network or IT specialist: complete networking requirements
- Step 3: Return the completed checklist

## Step 1: Instrument administrator: sign into your Connect account and select your connection type

Start by registering or signing into your Connect account at [thermofisher.com/connect](http://thermofisher.com/connect).

Have the username and password available during installation and training. The first username that is used on the instrument or linked to the instrument on Connect automatically becomes the instrument administrator. This administrator can add more users to the instrument with or without administrative privileges.

<p>Select one:</p> <p><input type="radio"/> Connect platform connection (recommended)</p> <p><input type="radio"/> Local area network (LAN) connection</p> <p><input type="radio"/> Direct connection</p> <p><b>Additional information about connection options and networking</b></p>	<p>Connect platform connection wired <i>or</i> wireless</p>  <p><b>Best</b></p>	<p>Local area network (LAN) connection wired <i>or</i> wireless</p>  <p><b>Better</b></p>	<p>Direct connection wired</p>  <p><b>Good</b></p>
--	---	--	--

If you cannot complete these requirements, please let us know why:

Signature indicates requirement has been met unless noted above.

Instrument administrator name:

Instrument administrator title:

Instrument administrator email address:

Instrument administrator telephone number:

Instrument administrator signature and date:

## Step 2: Network or IT specialist: complete networking requirements

Instrument information					
MAC address: <input type="text"/>	Room number: <input type="text"/>				
Serial number: <input type="text"/>	Port number: <input type="text"/>				
Computer information					
MAC address: <input type="text"/>	Room number: <input type="text"/>				
Service tag number: <input type="text"/>	Port number: <input type="text"/>				
Networking requirements					
<p><b>Firewall exception requirements:</b> The system should be configured behind a firewall. If outbound traffic is limited, these firewall exceptions are required to support system features:</p> <table border="0"> <tr> <td style="vertical-align: top;"> <p><b>Ports</b></p> <ul style="list-style-type: none"> <li>outbound 443</li> <li>outbound 80</li> </ul> <p><b>Remote services</b></p> <p><a href="https://apps.lifetechnologies.com/app-gateway">https://apps.lifetechnologies.com/app-gateway</a>  <a href="https://apps.thermofisher.com/g2i-gw/">https://apps.thermofisher.com/g2i-gw/</a>  <a href="https://drm.appliedbiosystems.com">https://drm.appliedbiosystems.com</a>  <a href="https://china.apps.thermofisher.com/app-gateway">https://china.apps.thermofisher.com/app-gateway</a> (China only)</p> </td> <td style="vertical-align: top;"> <p><b>File storage</b></p> <p><a href="https://dhap-cloudstorage.s3.amazonaws.com">https://dhap-cloudstorage.s3.amazonaws.com</a></p> <p><b>Firmware downloads</b></p> <p><a href="http://downloads.lifetechnologies.com/">http://downloads.lifetechnologies.com/</a>  <a href="http://downloads.thermofisher.com">http://downloads.thermofisher.com</a></p> </td> </tr> </table>		<p><b>Ports</b></p> <ul style="list-style-type: none"> <li>outbound 443</li> <li>outbound 80</li> </ul> <p><b>Remote services</b></p> <p><a href="https://apps.lifetechnologies.com/app-gateway">https://apps.lifetechnologies.com/app-gateway</a>  <a href="https://apps.thermofisher.com/g2i-gw/">https://apps.thermofisher.com/g2i-gw/</a>  <a href="https://drm.appliedbiosystems.com">https://drm.appliedbiosystems.com</a>  <a href="https://china.apps.thermofisher.com/app-gateway">https://china.apps.thermofisher.com/app-gateway</a> (China only)</p>	<p><b>File storage</b></p> <p><a href="https://dhap-cloudstorage.s3.amazonaws.com">https://dhap-cloudstorage.s3.amazonaws.com</a></p> <p><b>Firmware downloads</b></p> <p><a href="http://downloads.lifetechnologies.com/">http://downloads.lifetechnologies.com/</a>  <a href="http://downloads.thermofisher.com">http://downloads.thermofisher.com</a></p>		
<p><b>Ports</b></p> <ul style="list-style-type: none"> <li>outbound 443</li> <li>outbound 80</li> </ul> <p><b>Remote services</b></p> <p><a href="https://apps.lifetechnologies.com/app-gateway">https://apps.lifetechnologies.com/app-gateway</a>  <a href="https://apps.thermofisher.com/g2i-gw/">https://apps.thermofisher.com/g2i-gw/</a>  <a href="https://drm.appliedbiosystems.com">https://drm.appliedbiosystems.com</a>  <a href="https://china.apps.thermofisher.com/app-gateway">https://china.apps.thermofisher.com/app-gateway</a> (China only)</p>	<p><b>File storage</b></p> <p><a href="https://dhap-cloudstorage.s3.amazonaws.com">https://dhap-cloudstorage.s3.amazonaws.com</a></p> <p><b>Firmware downloads</b></p> <p><a href="http://downloads.lifetechnologies.com/">http://downloads.lifetechnologies.com/</a>  <a href="http://downloads.thermofisher.com">http://downloads.thermofisher.com</a></p>				
<p><b>Open port requirements:</b> If the system will use the <b>LAN connection option</b>, then these ports must be allowed on the network:</p> <ul style="list-style-type: none"> <li><b>5353:</b> Port 5353 is used for instrument discovery and uses multicast address 224.0.0.251.</li> <li><b>7000:</b> Port 7000 outbound calls are used by the to send commands to the system.</li> </ul>					
<p>If you cannot complete these requirements, please let us know why:</p> <div style="border: 1px solid black; height: 40px; width: 100%;"></div>					
<p>Signature indicates requirements have been met unless noted above.</p> <table border="0"> <tr> <td style="vertical-align: top;"> <p>Network or IT specialist name: <input type="text"/></p> <p>Network or IT specialist title: <input type="text"/></p> <p>Network or IT specialist email address: <input type="text"/></p> <p>Network or IT specialist telephone number: <input type="text"/></p> </td> <td style="vertical-align: top;"> <p>Network or IT specialist signature and date: <div style="border: 1px solid black; height: 100px; width: 100%;"></div></p> </td> </tr> <tr> <td></td> <td style="text-align: center;"> <input type="text"/> <input type="text"/> <input type="text"/> </td> </tr> </table>		<p>Network or IT specialist name: <input type="text"/></p> <p>Network or IT specialist title: <input type="text"/></p> <p>Network or IT specialist email address: <input type="text"/></p> <p>Network or IT specialist telephone number: <input type="text"/></p>	<p>Network or IT specialist signature and date: <div style="border: 1px solid black; height: 100px; width: 100%;"></div></p>		<input type="text"/> <input type="text"/> <input type="text"/>
<p>Network or IT specialist name: <input type="text"/></p> <p>Network or IT specialist title: <input type="text"/></p> <p>Network or IT specialist email address: <input type="text"/></p> <p>Network or IT specialist telephone number: <input type="text"/></p>	<p>Network or IT specialist signature and date: <div style="border: 1px solid black; height: 100px; width: 100%;"></div></p>				
	<input type="text"/> <input type="text"/> <input type="text"/>				

## Step 3: Return the completed checklist

To return the checklist:

- North America: Email the completed checklist to [TFC.Device.Connectivity.Support@thermofisher.com](mailto:TFC.Device.Connectivity.Support@thermofisher.com).
- All other regions: Unless asked to do otherwise, email the completed checklist to the Thermo Fisher Scientific representative who sent it to you.

## Additional information about connection options and networking

- **Connect platform connection (recommended):** Allows you to access all available features on our Connect platform and IoT-enabled functionality on the instrument such as Smart Help and voice commands and remote run monitoring on smart phone; includes support for Macintosh™ and Windows™ operating systems; does not require a Thermo Fisher Scientific-supplied computer, and allows you to connect multiple computers to the instrument via the Connect platform.
- **Local area network (LAN) connection:** Enables data sharing and run monitoring within the network; includes support for Macintosh™ and Windows™ operating systems; does not require a Thermo Fisher Scientific-supplied computer, and allows you to connect multiple computers to the instrument via the network.
- **Direct connection:** Supports a conventional system setup with a dedicated Thermo Fisher Scientific-supplied computer. No networking capability required.

For more information about service remote monitoring features, see the **Remote Monitoring and Diagnostics Service FAQs**.



Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256  
For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://thermofisher.com/symbols-definition).

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

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

**Revision history:** Pub. No. MAN0018160

Revision	Date	Description
A.0	18 April 2019	New document.

**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.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.