

Odyssey® CLx, Odyssey Fc, and Pearl® Imagers

Note: You must have Administrative privileges for the following steps.

Connect to Image Studio

1. Start Image Studio and create or open a Work Area.
2. If a connection is established, the buttons on the Acquire ribbon will become active.
3. If a network connection is not established, wait a few minutes then click the application button in the upper left-hand corner of the window.
4. Hover over Instrument and click Connect.

Computer Network Settings

Windows® 7 and 10

1. Go to Control Panel > Network and Sharing Center. Alternatively, go to Control Panel > Network & Internet > Network & Sharing Center.
2. Under “View your active networks,” click the Local Area Connection associated with the network card that is connected to the Odyssey and choose Properties.
3. Ensure that Internet Protocol Version 6 (TCP/IPv6) is checked.
4. Click Properties. Select “Obtain an IPv6 address automatically.”

Mac®

1. Click the Apple® icon in the upper left-hand corner of the screen and choose System Preferences.
2. In the Internet and Wireless section, click Network.
3. Click the Network object on the left associated with the network card that is connected to the instrument. Normally, the Network object is Ethernet.
4. In the Configure IPv4 section, choose Using DHCP.
5. Click Apply.

Troubleshooting

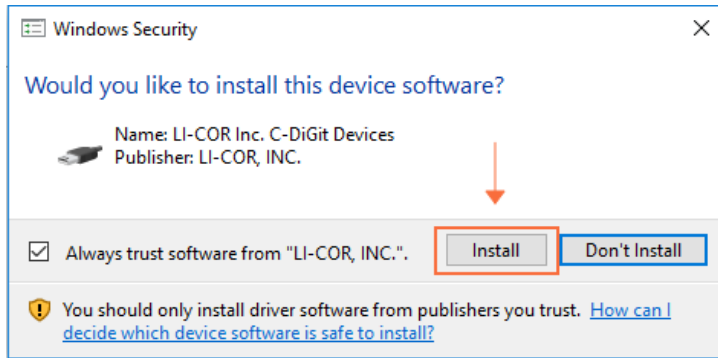
- If required, disable all firewall and anti-virus protection on the network card connected to the instrument.
- Do not block IPv6 traffic on the network card connected to the instrument.
- Leave Port 50000 open.
- Leave Port 5353 open for Bonjour mDNSResponder to Auto Discover the instrument.
- Do not block Bonjour Services.
- Do not block the Image Studio executable in a firewall.
- Do not block java running from "<Image Studio Install Folder>/jre" in a firewall.

C-DiGit® Blot Scanner

Ensure you have the correct permissions to install software on the computer you want to use.

Note: If you are using Windows 10, a separate driver must be installed to connect your C-DiGit Blot Scanner to Image Studio Software via USB.

1. Ensure your C-DiGit Blot Scanner is powered off and not connected to your computer.
2. Ensure Image Studio Software is closed.
3. [Download the driver.](#)
4. Once the installer has downloaded, click the installer to start the Setup Wizard.
5. Click Next and proceed through the Setup Wizard.
6. If the Windows Security dialog appears during the install, click Install or the driver will not install.



7. After the driver has been installed, connect the USB cable from the C-DiGit® Blot Scanner to the computer.
8. Turn on the C-DiGit Blot Scanner.
9. Start Image Studio™ Software, choose a Work Area, and select C-DiGit Blot Scanner in the Instrument Selection dialog.

Odyssey® Sa and Aerius Imagers

For an Odyssey Sa or Aerius that is directly connected to the computer, follow the instructions below. For an Odyssey Sa or Aerius that is connected to a network, please refer to its support page or contact Biotechnology Support (biohelp@licor.com).

Note: You must have Administrative privileges for the following steps.

Step 1. Configure Network Card of Computer

Windows® 7 and 10

1. Go to Control Panel > Network and Sharing Center. Alternatively, go to Control Panel > Network & Internet > Network & Sharing Center.
2. Under “View your active networks,” click the Local Area Connection associated with the network card that is connected to the Odyssey and choose Properties.
3. Select Internet Protocol Version 4 (TCP/IPv4) and click Properties.
4. Check “Use the following IP address” and enter “172.24.41.181.”
5. If not automatically filled in, enter “255.255.0.0” for the Subnet mask.

Mac®

1. Click the Apple® icon in the upper left-hand corner of the screen and choose System Preferences.
2. In the Internet and Wireless section, click Network.
3. Click the Network object on the left associated with the network card that is connected to the instrument. Normally, the Network object is Ethernet.
4. In the Configure IPv4 section, choose Manually and enter “172.24.41.181.”
5. If not automatically filled in, enter “255.255.0.0” for the Subnet mask.
6. Click Apply.

Step 2. Connect to Image Studio

1. Start Image Studio and create or open a Work Area.
2. Enter a username and password.
3. If a network connection is not established, click the application button in the upper left-hand corner of the window.
4. Hover over Instrument. Click Management then click Scanner List.
5. If there are no instruments listed, click Auto Discover.
6. If nothing appears in the list, click Add and enter the information manually.

Note: “Name” can be anything meaningful to the user. “Host name” is the IP address. Do not include numbers after the “/”. Additionally, check the IP address of the instrument by pressing the Next button on the front panel of the instrument 4-5 times. The default IP address is 172.24.41.180.

7. Click OK then OK again to close.

8. Go back to Application > Instrument > Connect. When the login dialog appears, select the instrument and enter the username and password.

Troubleshooting

- If required, disable all firewall and anti-virus protection on the network card connected to the instrument.
- Leave Port 80 open.
- Do not block the Image Studio™ Software executable in a firewall.
- Do not block java running from "<Image Studio Install Folder>/jre" in a firewall.

Odyssey® Classic Imager

For an Odyssey Classic that is directly connected to the computer, follow the instructions below.

For an Odyssey Classic that is connected to a network, please refer to its support page or contact Biotechnology Support (biohelp@licor.com).

Note: You must have Administrative privileges for the following steps.

Step 1. Configure Network Card of Computer

Windows® 7 and 10

1. Go to Control Panel > Network and Sharing Center. Alternatively, go to Control Panel > Network & Internet > Network & Sharing Center.
2. Under "View your active networks," click the Local Area Connection associated with the network card that is connected to the Odyssey and choose Properties.
3. Select Internet Protocol Version 4 (TCP/IPv4) and click Properties.
4. Check "Use the following IP address" and enter "172.24.41.181."
5. If not automatically filled in, enter "255.255.0.0" for the Subnet mask.

Mac®

1. Click the Apple® icon in the upper left-hand corner of the screen and choose System Preferences.
2. In the Internet and Wireless section, click Network.
3. Click the Network object on the left associated with the network card that is connected to the instrument. Normally, the Network object is Ethernet.
4. In the Configure IPv4 section, choose Manually and enter "172.24.41.181."
5. If not automatically filled in, enter "255.255.0.0" for the Subnet mask.
6. Click Apply.

Step 2. Connect to Image Studio

1. Start Image Studio and create or open a Work Area.
2. Enter a username and password.
3. If a network connection is not established, click the application button in the upper left-hand corner of the window.
4. Hover over Instrument. Click Management then click Scanner List.
5. If there are no instruments listed, click Auto Discover.
6. If nothing appears in the list, click Add and enter the information manually.

Note: "Name" can be anything meaningful to the user. "Host name" is the IP address. Do not include numbers after the "/". Additionally, check the IP address of the instrument by pressing the Next button on the front panel of the instrument 4-5 times. The default IP address is 172.24.41.180.

7. Click OK then OK again to close.
8. Go back to Application > Instrument > Connect. When the login dialog appears, select the instrument and enter the username and password.

Troubleshooting

- If required, disable all firewall and anti-virus protection on the network card connected to the instrument.
- Leave Port 80 open.
- Do not block the Image Studio™ Software executable in a firewall.
- Do not block java running from "<Image Studio Install Folder>/jre" in a firewall.



© 2020 LI-COR, Inc. LI-COR, C-DiGit, Odyssey, Pearl, and Image Studio are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com