

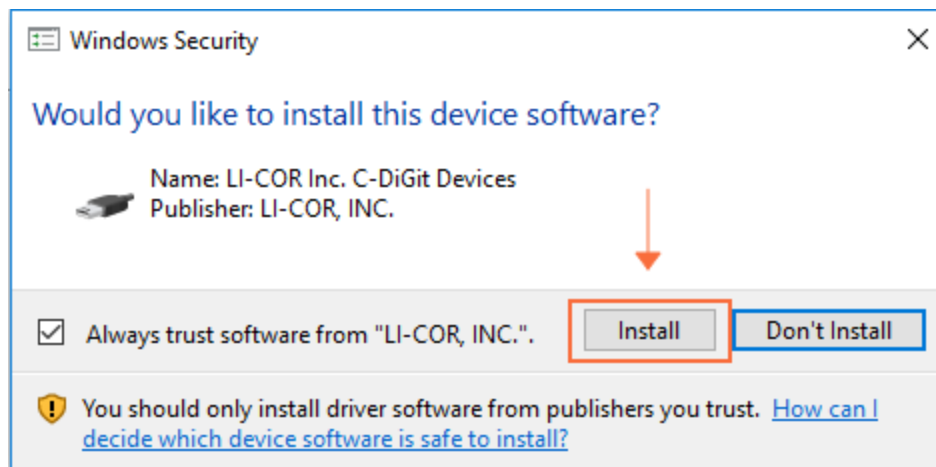
# Install the C-DiGit® Blot Scanner Windows 10 Driver

**LI-COR**

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.

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

1. | Ensure your C-DiGit Blot Scanner is powered off and is 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, choose a Work Area, and select C-DiGit Blot Scanner in the Instrument Selection dialog.

You should now be able to use the C-DiGit with the USB Cable on Windows 10.

© 2017 LI-COR, Inc. LI-COR, C-DiGit, 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 is an ISO 9001 registered company.



Global Headquarters: 4647 Superior Street • Lincoln, NE 68504  
Toll free: 800-645-4267 • Fax: +1-402-467-0819 • [biosales@licor.com](mailto:biosales@licor.com)

LI-COR GmbH, Germany, Serving Europe, Africa, and the Middle East.  
LI-COR Biosciences GmbH • Siemensstraße 25A • 61352 Bad Homburg • Germany  
Phone: +49 (0) 6172 17 17 771 • Fax: +49 (0) 6172 17 17 799 • [bio-eu@licor.com](mailto:bio-eu@licor.com)

LI-COR Ltd., United Kingdom, Serving Denmark, Finland, Iceland, Ireland, Norway, Sweden, and UK  
LI-COR Biosciences UK Ltd. • St. John's Innovation Centre • Cowley Road • Cambridge • CB4 0WS  
United Kingdom • Phone: +44 (0) 1223 422104 • Fax: +44 (0) 1223 422105 • [bio-eu@licor.com](mailto:bio-eu@licor.com)

LI-COR Biotechnology Distributor Network: [www.licor.com/bio/distributors](http://www.licor.com/bio/distributors)

[www.licor.com/bio](http://www.licor.com/bio)