



Image Studio™ 4.x.x Installation Instructions for the C-DiGit® Blot Scanner



LI-COR®

Image Studio 4.x.x Installation Instructions

Mac® OS X® 10.8 (Mountain Lion) or 10.9 (Mavericks):

Installing via CD:

- 1) Place installation CD into drive.
- 2) If installation does not start immediately, browse to the CD and double-click on MacImageStudioDigit_4.x.x.dmg.
- 3) Click **Agree** to accept the license agreement.
- 4) To install, drag the Image Studio icon into the Applications folder (right).



Installing via download:

- 1) Go to your Mac System Preferences and click **Security & Privacy** to open the Security & Privacy window.
- 2) Click the lock in the bottom left corner of the window to allow editing, if necessary.
- 3) Under the **General** Tab, click **Mac App Store and identified developers** under the **Allow applications downloaded from** option.
- 4) Download, and double-click on the MacImageStudioDigit_4.x.x.dmg file on your computer.
 - a. Click **Agree** to accept the license agreement.
 - b. To install, drag the Image Studio icon into the Applications folder.
 - c. Open the Image Studio Application by double-clicking on the Image Studio icon.
 - d. If a window appears that says "ImageStudio.app is an application downloaded from the internet. Are you sure you want to open it?" click **Open**.
- 5) If you have purchased additional application keys, you will need to import those keys again to utilize their functionality.

Windows 7, 8, or 8.1 Operating Systems:

IMPORTANT: The installer must have administrator privileges on the computer to install Image Studio 4.x.x. Log on to the computer as Administrator, or log on as a user who has administrative permissions. Computer requirements: Minimum 4 GB of RAM, Win 7, 8, or 8.1 Operating System.

Installing via CD:

- 1) Place installation CD into drive.
- 2) If installation does not start immediately, browse to the CD and double-click on Win_Image StudioDigit_4.x.x.exe.
- 3) Click **Next** and follow instructions in the Setup Wizard.

Installing via download:

- 1) Download and double-click on the Win_Image StudioDigit_4.x.x.exe file on your computer.
- 2) Click **Next** and follow the instructions in the Setup Wizard.

Image Studio Setup Wizard

- 1) Open Win Image Studio Installer 4.x.x on your computer.

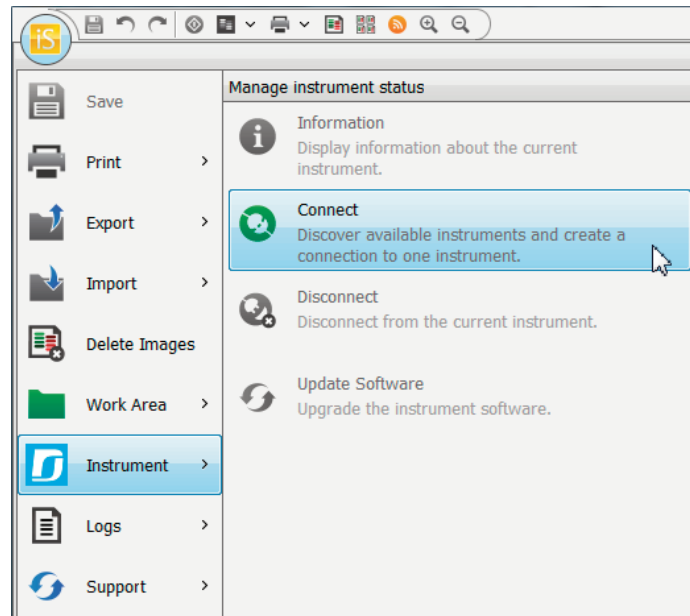


- 2) Follow instructions in the Setup Wizard.

NOTE: After the Image Studio installation process is completed, installation of the communication program (mDNSResponder) will begin automatically. If upgrading from a previous version of Image Studio, this installation may not be needed.

- 3) Please refer to instructions included with your additional application keys, if applicable, for new installations.
- 4) If you are upgrading from a previous version of Image Studio software on a Mac, you will need to reinstall additional application keys.

Instrument Connection



The first time you connect, power the C-DiGit Blot Scanner on and wait several minutes to allow the instrument to establish a network address.

- 1) Click the **Image Studio** icon in the upper left-hand corner of the Image Studio window.
- 2) Hover over Instrument and click **Connect**.

The software searches for and discovers the C-DiGit Blot Scanner automatically.

NOTE: If you experience connection problems, a connection diagnostic will launch to help you diagnose the problem.

LI-COR is an ISO 9001 registered company. ©2014 LI-COR, Inc. LI-COR, Odyssey, 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. For patent information, visit www.licor.com/patents.



4647 Superior Street • P.O. Box 4000 • Lincoln, Nebraska 68504 USA
Phone: 402-467-0700 • Toll free: 800-645-4267 • Fax: 402-467-0819 • www.licor.com/bio

LI-COR GmbH Germany, Serving Europe, Middle East, and Africa: +49 (0) 6172 17 17 771

LI-COR UK Ltd. UK, Serving UK, Ireland, and Scandinavia: +44 (0) 1223 422104

All other countries, contact LI-COR Biosciences or a local LI-COR distributor:

<http://www.licor.com/distributors>