



Image Studio™ Software Installation with the Pearl® Imager

LI-COR

Image Studio Software 5.x.x Installation Instructions

Mac OS X 10.9, 10.10, 10.11, or 10.12:

Installing via CD

- 1) Place installation CD into drive.
- 2) If installation does not start immediately, browse to the CD and double-click the installer file.
- 3) Click **Agree** to accept the license agreement.
- 4) To install, drag the Image Studio™ Software icon into the Applications folder (below).



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 the installer file on your computer.
- 5) Click **Agree** to accept the license agreement.
- 6) To install, drag the Image Studio™ Software icon into the Applications folder.
- 7) Open the Image Studio Application by double-clicking the Image Studio™ Software icon.
- 8) 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**.

Windows® 7, 8, 8.1, or 10 Operating Systems:

Note: The installer must have administrator privileges on the computer to install Image Studio™ Software 5.x.x. Log on to the computer as an Administrator, or log on as a user who has administrative permissions. Computer requirements: Minimum 4 GB of RAM and Windows 7, 8, 8.1, or 10 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 Studio Full_5.x.x.exe.
- 3) Click **Next** and follow the instructions in the Setup Wizard.

Installing via download

- 1) Download and double-click the Win_Image Studio Full_5.x.x.exe file on your computer.
- 2) Click **Next** and follow the instructions in the Setup Wizard.

Image Studio Software Setup Wizard:

- 1) Open Win Image Studio Installer 5.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.

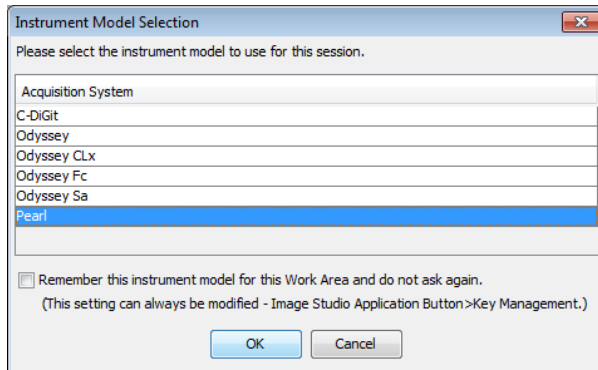
Note: If you are upgrading from a previous version of Image Studio™ Software on a Mac, you will need to reinstall additional application keys.

Connect to the Instrument:

The first time you connect:

- 1) Power the Pearl Imaging System on and wait several minutes to allow the instrument to establish a network address.
- 2) Start Image Studio™ Software. Instruments will be detected automatically.
- 3) The **Instrument Model Selection** dialog will open. Select Pearl and click **OK**.

Image Studio™ Software will open your **Work Area** and search for the Pearl Imaging System .



Note: You can save this as your default instrument for the current Work Area by checking 'Remember this instrument model for this Work Area and do not ask again' box.

- 4) If the software discovers multiple instruments, it will prompt you to select the instrument to be used. Select the Serial Number of the desired instrument, and click **OK**.

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



© 2020 LI-COR, Inc. LI-COR, 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