



Image Studio™ Software 5.x

Image Acquisition and Import

- **Application-Driven Ribbon Interface (Figure 1)**
 - Each analysis type has a corresponding ribbon with tools for analysis and formatting that are only relevant for that application

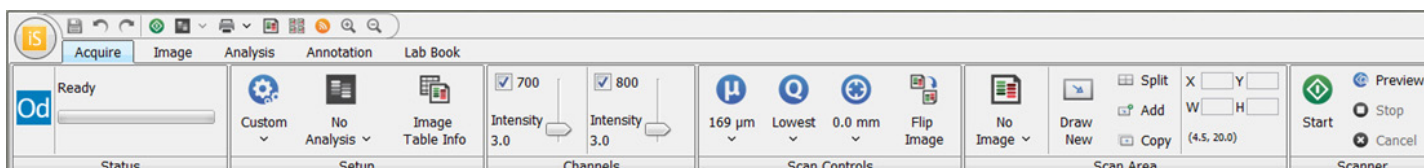


Figure 1. Image Studio is a ribbon-based application that displays analysis and formatting tools for user selection and implementation.

- **Ability to Scan Multiple Areas in a Single Scan (for Odyssey® Classic and Odyssey CLx Imaging Systems)**
 - With an enhanced acquisition interface, the user can divide a single scan into multiple scan areas
- **Ability to Import TIFF, PNG, or JPEG Files**

Data Analysis

- **Definition of Lanes and Bands**
 - Simply define the sample area and the total number of lanes
- **Placement of Shapes by Simple Point-and-Click Method**
 - Click on a band to see a rectangle placed around it automatically
- **Rotation of Objects**
 - If a band or blot are not in the desired orientation, objects can be rotated as a group or individually
- **Data Analysis Tools**
 - The user has full control of the data display and can rearrange each column through sorting and filtering
 - Create customized charts
- **Annotation Controls**
 - Simple to change fonts and controls, including rotational control of text and arrows
- **Linking of Look-Up-Tables**
 - View different scans under identical display conditions
- **Supports Many Types of Analysis, Including:**
 - Western Blots
 - In-Cell Western™ Assays
 - Multi-Well Plates
 - Coomassie Gel Documentation
 - MPX™ Western Blots
 - EMSAs
 - Tissue Imaging
 - Grid Arrays
 - DNA Gels
 - Small Animal Imaging

Image and Data Export and Printing

- **Image Export Tools**
 - Export experimental images and tables for further analysis in several downstream software packages, such as Adobe® Photoshop, Microsoft® PowerPoint, and GraphPad Prism® Software
 - Image and accompanying information can be easily transferred from one computer to another for data analysis
 - New sizing and resolution options are now available, as well as expanded file formats
- **Data Export Tools**
 - Launch an external spreadsheet program and automatically import selected rows of data into the spreadsheet program
- **Customizable Lab Notebooks (Figure 2)**
 - Ability to create a direct-to-PDF export of data with separate templates for each analysis type
 - PDF format is fully customizable to meet different documentation requirements



Figure 2. Create electronic or hard-copy lab reports customized to meet specific needs and documentation requirements, such as GLP or ISO.

Data Storage

- **Simplified File Access**
 - Easier, more convenient access to user data

Minimum Requirements

- **Odyssey® Classic Infrared Imaging System:**
 - SN <3000: Embedded software, version 2.1.15
 - SN >3000: Embedded software, version 3.0.X or higher
- **Odyssey CLx Infrared Imaging System: Embedded software, version 1.0.9 or higher**
- **Odyssey Fc Dual-Mode Imaging System: Embedded software, version 1.0.14 or higher**
- **Aerius Imaging System: Embedded software, version 2.0.14**
- **Odyssey Sa/Sa Express Imaging System: Embedded software, version 2.0.13**
- **C-DiGit® Blot Scanner: Embedded software, version 1.1.3 or higher**
- **Pearl® Imaging Systems: Embedded software, version 1.2.6 or higher**
- **Pearl Trilogy Imaging System: Embedded software, version 1.3.x**
- **Computer Operating Systems:**
 - Windows Systems: 4GB RAM • Windows® 7, 8, 8.1, or 10
 - Macintosh Computer: 4GB RAM • OS 10.9, 10.10, 10.11, and 10.12

Get Image Studio Today at www.licor.com/imagestudio

LI-COR is an ISO 9001:2015 registered company. ©2020 LI-COR, Inc. Specifications subject to change. LI-COR, In-Cell Western, MPX, Pearl, C-DiGit, Image Studio, and Odyssey are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. Macintosh is a trademark of Apple Inc. All other trademarks belong to their respective owners. For patent information, visit www.licor.com/patents



4647 Superior Street • Lincoln, NE 68504 • Phone: +1-402-467-0700 • Toll free: 800-645-4267 • Fax: +1-402-467-0819 • biosales@licor.com

Regional Offices

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

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

LI-COR Distributor Network: www.licor.com/bio/distributors • www.licor.com/bio