



Image Studio 3.1 Imaging Software

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

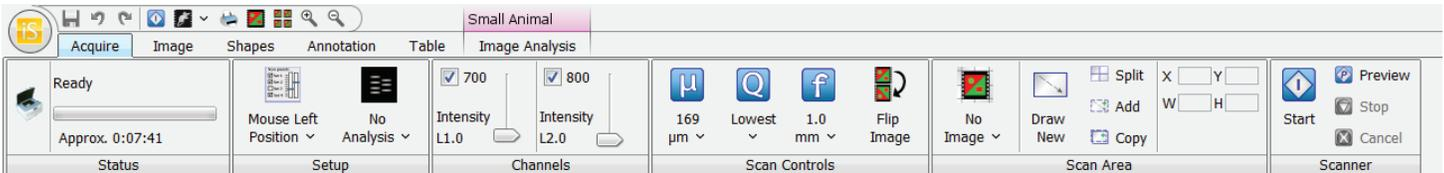


Figure 1. Image Studio is a ribbon-based application that displays analysis and formatting tools for user selection and implementation.

- **Ability to scan multiple areas in a single scan**
 - With an enhanced acquisition interface, the user can divide a single scan into multiple scan areas
- **Ability to import .tif, .jpg and .png 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
- **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 & 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

Aerius Imaging System: Embedded software, version 2.0.14

Odyssey Sa Imaging System: Embedded software, version 2.0.13

Pearl® Imaging System: Embedded software, version 1.2.6

Computer Operating Systems:

- | | |
|---|--|
| <ul style="list-style-type: none"> ▫ Windows® Systems <ul style="list-style-type: none"> ◆ 4GB RAM ◆ Windows® 7, XP, or Windows Vista® Operating System | <ul style="list-style-type: none"> ▫ Macintosh® Computer <ul style="list-style-type: none"> ◆ 2GB RAM ◆ OS 10.6 or above |
|---|--|

LI-COR is an ISO 9001 registered company. ©2013 LI-COR, Inc. Specifications subject to change.
LI-COR, In-Cell Western, MPX, and Odyssey 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 | PO Box 4000 | Lincoln, Nebraska 68504 USA

LI-COR Biosciences North America: 800-645-4267 | 402-467-0700

Fax: 402-467-0819 | Technical Support: 800-645-4260

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

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

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

www.licor.com/distributors | www.licor.com/bio