

LI-COR Aerius™ Automated Infrared Imaging System

Advanced Automation for Protein Discovery

Aerius introduces a new level of automation for In-Cell Western assays and an expanding list of other infrared plate-based assays. Aerius provides fully automated data analysis of In-Cell Westerns for true, one-touch, walk-away automation.

Aerius combines the infrared advantages of the Odyssey® Infrared Imaging System—extreme sensitivity and quantitative analysis of proteins—with a fast and efficient microplate reader. In addition to processing up to 30 microplates per run for large-scale analysis, the Aerius scans single membranes or microplates for validating antibodies.



Aerius is available with an optional Bio-Tek® Bio-Stack™ Automated Microplate Stacking System. LI-COR also offers COM software for integration with other robotic systems.



Scanning, spot finding, intensity measurements, and percent response calculations for In-Cell Westerns are fully automatic on the Aerius System.

SCANNING FEATURES

- Two-color scanning for simultaneous detection of two protein targets or the use of one channel for normalization.
- 20 – 500 µm resolution.
- Scans standard microplates or membranes using the Aerius Membrane Carrier.
- Integration of scanning and automatic analysis for In-Cell Western assays.
- Simple operation. Scanning and In-Cell Western analysis are reduced to choosing an express run template.

AUTOMATION FEATURES

User choice of fully automated or interactive analysis of scanned images.

Bio-Tek® Bio-Stack™ Automated Microplate Stacking System

- Aerius Express run software integrates with the plate stacker to serially scan up to 30 microplates per run.
- Express run templates allow you to save configurations for an entire stack of plates.
- During a run, each plate is scanned, analyzed with In-Cell Western analysis software, and data output in reports or data files.
- Automatic output to a network location in delimited text format for spreadsheets, laboratory information systems, or analysis software.

COM Interface

- A programming, language-independent interface to Aerius useful for small-scale laboratory automation.

SPOT FINDING FOR MICROPLATES

- Automatic: Well locations are determined by analyzing the image and placing an array of markers over the wells.
- Manual: A user-defined template determines initial location of well markers.

LANE FINDING

- Flexible lane tools make it easy to mark both straight and curved lanes.
- Bands are found automatically and accurately as lanes are marked.
- Band finding can be verified by viewing intensity curves for each lane.
- Band markers can be re-sized and moved on the image.

BAND SIZING

- The image is automatically calibrated for molecular weight after standards are identified.
- Molecular weights are quickly and accurately corrected on images with smiles.

- Molecular weight standards need only be run using one of the two colors, reducing reagent costs.
- All bands in sample lanes are automatically sized after molecular weight standards are identified.

QUANTIFICATION

- A variety of shape tools make it easy to mark and analyze areas on the image.
- A Details View with curve data and an enlarged image can be opened to verify accurate placement of markers.
- Quantification data can be reviewed in table format for rapid comparison.

CUSTOM REPORT GENERATION

- Customizable reports for In-Cell Westerns, band sizing, or quantification.
- Export macros can start programs such as Microsoft® Excel and export the data directly into a new document.

SECURITY

- Only authorized users can start scans or analyze images.
- Access permission can be changed only by users with Administrator accounts.

AERIUS HARDWARE SPECIFICATIONS

- Laser Lifetime: 60,000 hours typical.
- 700 Channel Laser Source: Solid-state diode laser at 685 nm.
- 800 Channel Laser Source: Solid-state diode laser at 785 nm.
- Detectors: Silicon avalanche photodiodes.
- Scanning Speed: 6 minutes per microplate with 100 μ m resolution.
- Power Requirements: 100-127 VAC or 200-240 VAC.
- Dimensions:
Instrument Only: 36H \times 45W \times 57D cm (14 \times 18 \times 22 inches).
Instrument with Stacker: 85H \times 110W \times 45D cm (33 \times 44 \times 18 inches)
- Bar Code Reader symbology: Code 39, preferred. Compatible with Code 128 and other symbologies. Contact LI-COR for more information.

Specifications subject to change.

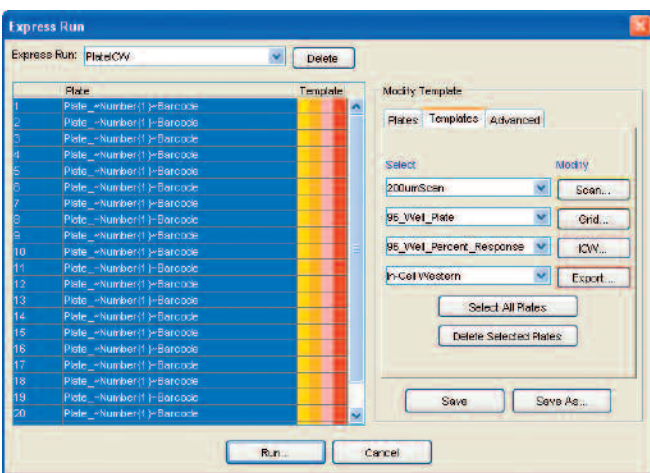
AERIUS SYSTEM CONFIGURATIONS

Aerius Automated Infrared Imaging System

- Includes the Aerius Instrument, computer with LCD monitor, Aerius Application Software, and Aerius Accessory Pack which includes a membrane carrier.

Aerius Automated Infrared Imaging System—With Plate Stacker

- Includes the Aerius Instrument, Bio-Stack Microplate Stacker, Bio-Stack Plate Stacker install kit (includes stand, alignment tools and cables), computer with LCD monitor, Aerius Application Software, and Aerius Accessory Pack which includes a membrane carrier.



For repetitive runs, scanning is as simple as choosing an express run template that defines scanning and analysis for up to 30 microplates.

LI-COR[®]

Biosciences

4308 Progressive Ave. • P.O. Box 4000 • Lincoln, Nebraska 68504 USA
North America: 800-645-4267 • International: 402-467-0700 • Fax: 402-467-0819
LI-COR GmbH (Germany, Austria, Switzerland): +49 (0) 6172 17 17 771
LI-COR UK Ltd.: +44 (0) 1223 422104 • www.licor.com/aerius

LI-COR is an ISO 9001 registered company. © 2004 LI-COR Inc. LI-COR, Odyssey, and Aerius are trademarks or registered trademarks of LI-COR Inc. Microsoft is a registered trademark of Microsoft Corporation. Bio-Stack and Bio-Tek are trademarks or registered trademarks of Bio-Tek Instruments, Inc.