

Living Image®

Release Notes Version 4.8.2

1 Contents

1	Contents	1
2	Purpose	1
3	Improvements	1
4	Bug Fixes	2
5	Known Issues	2
6	Changes to the Product Documentation	3

2 Purpose

This document briefly describes improvements, and relevant bug fixes in Living Image® (LI) 4.8.2, along with the changes to the hardware and software manuals.

LI 4.8.2 supports:

- IVIS Lumina III
- IVIS Lumina LT
- IVIS Lumina XRMS
- IVIS Spectrum
- IVIS Spectrum CT
- IVIS Spectrum 2
- IVIS Spectrum CT 2

3 Improvements

This release includes enhancements such as:

- All branding has been updated to Revvity Inc.

- Support for the new IVIS Spectrum 2 and IVIS SpectrumCT 2 instruments.
- Camera chiller utility. A service-based function to maintain the camera temperature setpoint when LI is not running.
- New agents: redFluc, nanoluc, and luc 2
- IVIS Spectrum 2, IVIS Spectrum CT 2, Spectrum, Spectrum CT support the SmartTray™ and will support segmentation. See the known issues section.

4 Bug Fixes

The release includes these major bug fixes:

- National Instruments card is now detected by the software on a retry basis before initialization, which allows Spectrum Instruments to use LI4.8.2.
- LI no longer crashes when acquiring a sequence of multiple images and when the photograph was not reused.
- Surface Topology Tool palette does not disappear now.
- Only the 5-mouse tray checkbox OR the MIS checkbox is available when the other is in checked.
- A correction to the UI in the ROI Tools palette when loading a CT dataset that prevented the ability to load/save 3D ROI.
- A warning when the number of keyframes exceeds the time scale limit in 3D animation.
- The Recent Files feature has been fixed.
- A crash on exiting LI for analysis workstations has been fixed.

Several minor bugs have also been addressed but are not listed here.

5 Known Issues

There is one known critical severity issue:

- The SmartTray™ segmentation functionality does not work in all cases.

There are several major severity issues open and are scheduled to be fixed in 4.8.3 or a future version, such as:

- Tool palette disappears when opening some DLIT images.
- Lens distortion correction is not applied to a Structure Only image.
- Issues with auto exposure not correctly setting Binning and/or F-Stop and exposure.
- On a Lumina XRMS, the scintillator arm will collide with the stage if multiple X-ray images are taken without a corresponding new photographic image. The

workaround is to unclick the 'reuse photo' button to force a photograph and x-ray image at each line in a sequence.

There are several medium severity software issues open, such as:

- When you change the text size in user preferences, the units field overlaps with the color scale field.
- The initial position of the mouse organ atlas does not match the optical/CT data. It is easily repositioned manually.
- The fluorescence signal does not always align with the optical image using the image math functions.

There are several low severities instrument issues open, such as:

- (Lumina S5, Lumina X5) The animal height at FOV D is capped at 1.98cm.
- (Lumina X5) Occasionally if the scintillator arm moves past the operating position during initiation, the initialization fails and reports the incorrect motor has failed. The workaround: manually move the scintillator arm to a location within the operating range.
- Living Image 4.8.2 has not been tested with Spectrum CT using the Dexela PCIE card. The last version tested with this instrument is LI 4.7.3.

6 Changes to the Product Documentation:

- Combined the software (LI 4.8.2) and hardware manuals.
- Added relevant software and hardware changes, as noted throughout these release notes.
- Rebranded technical notes