



Living Image™

Release Notes

Living Image Analysis software for Macintosh (macOS).

Version 4.8.2

Purpose

The Living Image 4.8.2 analysis software for Macintosh computers provides Revvity Branding, supports new Apple Silicon-based computers, includes bug fixes, and uses a new installer for expanded macOS release support.

Analysis PC Requirements

Living Image 4.8.2 analysis software for Macintosh computers runs on the following macOS versions:

- macOS 10.15 Catalina (Intel chipset)
- macOS 11 Big Sur (Intel chipset)
- macOS 12 Monterey (Intel or Apple chipset)
- macOS 13 Ventura (Intel or Apple chipset)
- macOS 14 Sonoma (Intel or Apple chipset)

4GB RAM or higher is recommended for 3D and IVIS Spectrum CT data analysis.

For Apple computers running Intel chipsets, an ATI Radeon video card or certain Intel Iris Graphics chipsets are required for multi-modality support. See the section on Video Card Requirements.

For computers running Apple chipsets, the built-in graphics are adequate, no additional cards are needed.



Living Image™

Video Card Requirements

For Apple computers running older OS versions and Intel chipsets, to use 3D multi-modality analysis, a graphics card with the following minimal requirements must be met. If the graphics card does not meet these requirements, the 3D multi-modality tools will not appear in the tool palate.

Specification	Description
OpenGL Version	OpenGL 2.0 and above
OpenGL Extension	GL-EXT-Texture3D
Graphics card memory	Recommended: 1GB (Dedicated)

New Features/Functions

New installer with Revvity branding and improved support for current macOS versions.

Revvity Branding. All user-facing software items reflect the new company name, colors, and logo.

Support for Apple Silicon. OpenGL is deprecated and replaced with QT classes.

Support for macOS versions Catalina through Sonoma

Bugs Fixed

- Unable to Maximize 3D Plot Window
- Several other minor or trivial bugs



Living Image™

Known Issues

The following are known issues planned for future versions of Living Image Analysis Software for Macintosh

- Selected window stays in front even after surface reconstruction
- Aborting while saving a movie in Kinetic datasets can result in a hang
- Option to choose the installation path in the installer is not functional
- Mac installer lacks build number information
- Application crash in macOS11 Big Sur on "Save as a Movie" in kinetic datasets

Installation

Please refer to the installation and Licensing guide for instructions.