

Living Image 4.8.2 mac - ReadMe.txt

1. WHAT IS NEW IN LIVING IMAGE 4.8.2 MAC RELEASE?

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

For the full overview of Living Image 4.8.2 improvements and bug fixes, please refer to the Release Notes.

2. SYSTEM 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.

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)

INSTALLATION INSTRUCTIONS

Please, refer to Installation & Licensing Guide for step-by-step installation instructions.

4. GETTING HELP

Revvity, Inc.
68 Elm Street
Telephone: +1.508.435.9500
www.revvity.com

Technical Support
www.revvity.com/contact-us
Telephone: +1.800.762.4000
Fax: +1.203.944.4904

Copyright
© 2019 - 2024 Revvity, Inc. and its parent, affiliated, and subsidiary companies. All rights reserved, including but not limited to those rights to reproduce this publication or parts thereof.