

Living Image 4.8.2 - ReadMe.txt

1. WHAT IS NEW IN LIVING IMAGE 4.8.2?

Living Image 4.8.2 serves as an update for Living Image 4.8.0 on all Revvity branded instruments.

On instruments listed below LI 4.8.2 has been fully tested:

Lumina III

Lumina LT

Lumina XRMS

Spectrum

Spectrum CT

Spectrum 2

SpectrumCT 2

All user facing branding has been updated to Revvity Inc.

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

2. SYSTEM REQUIREMENTS

For 4.8.2, we still need to install and run as admin (using the local IVISUSER account should cover this).

For optimum performance, the Living Image 4.8.2 software requires hardware and software that meet or exceed the following requirements:

Windows 10 64-bit

2GHz Quad Core (i5, i7 equivalent or better) processor

8GB RAM is the absolute minimum for IVIS Lumina instruments.

16 GB RAM is the absolute minimum for all Spectrum class instruments.

32 GB RAM is the absolute minimum for all Spectrum class instruments.

3D multi-Modality tools require that the graphics processing unit (GPU) meets the minimum specifications shown below.

If the appropriate license is not installed or the GPU does not meet these specifications, the 3D multi-Modality tools will not appear in the tool palette.

OpenGL Version Requirement* OpenGL 2.0 and above

OpenGL Extension Requirement* GL-EXT-Texture3D

Graphics Card Memory: Recommended: 1GB (Dedicated)

Consumer Graphics Cards (Desktop/ Mobile, Windows/Mac)

Supported:

- NVIDIA® GeForce® 8 Series and above (8, 9, 100, 200, 300 and 400 series)
- Intel HD 3000 and above and Intel Iris/Iris Pro Graphics (Mac)

Recommended:

- Desktop - NVIDIA GeForce GT 240 and above
- Mobile - NVIDIA GeForce GT 230M and above

Workstation Graphics Cards (Desktop/Mobile, Windows/Mac) Supported:

- NVIDIA® Quadro® NVS Series and Above (NVS & FX series)

- Recommended:

- Desktop - Quadro FX 1800 and above
- Mobile - Quadro FX 880M and above

*If these specifications are not met, the 3D multi-Modality tools will not appear in the tool palette.

3. INSTALLATION INSTRUCTIONS

Living Image 4.8.2 offers different installation packages depending on the computer operating system where Living Image is being installed. Choose from Windows 64-bit and MAC OS options.

3.1. Windows 64-bit version

- Recommended for data acquisition on all IVIS instruments excluding IVIS Spectrum CT.

- Recommended for analysis on Windows 10 64-bit operating systems.
- Supported for acquisition on all IVIS instruments excluding the older Spectrum CT with the PCIE Dexela x-ray detector set.

NOTE: make sure license server is whitelisted by your IT Department. Server information is:

Domain name: activation.caliperls.com

IP address: 165.88.8.121

Port: 443

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.