



Modalities combined. Intelligence driven. Efficiency reimaged.

Streamline *in vivo* imaging analysis with
Living Image Synergy AI multimodality software.

Living Image™ Synergy AI software applies AI-powered algorithms that help accelerate *in vivo* analysis and reduce variability, while easily integrating optical, microCT, and ultrasound data into a unified imaging workflow.



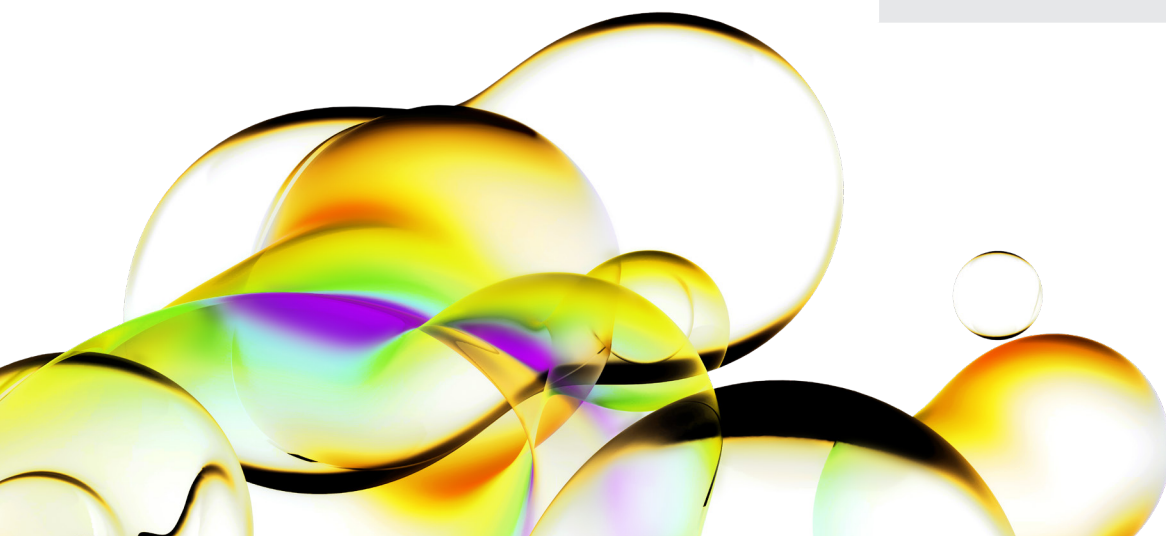
Features:

- Multiple AI-assisted organs, tumors, and whole-body segmentation models
- One analysis platform for multiple compatible Revvity modalities
- Automated bioluminescence kinetic curves
- Streamlined batch processing of studies
- Intuitive co-registration capability

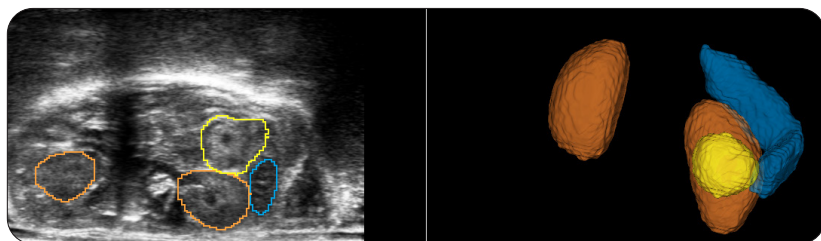
What's new in version 1.1.0

- Additional AI organ segmentation models
- Fluorescence imaging visualization, segmentation, and measurement support
- Editor effect sculpting tool for easier manipulation of 3D segmentations in 3D space

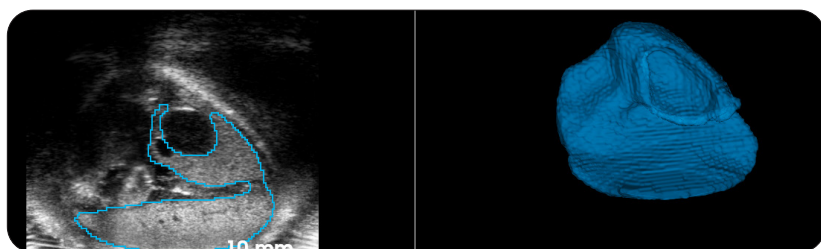
Imaging system / modality	AI segmentation model
Vega™ ultrasound	▪ Spleen ▪ Kidney ▪ Liver ▪ Tumors (orthotopic pancreatic, and subcutaneous tumors)
Quantum™ GX3 uCT	▪ Lung (total volume and air volume)
IVIS™ optical platform	▪ Head ▪ Thorax ▪ Abdomen ▪ Body contour



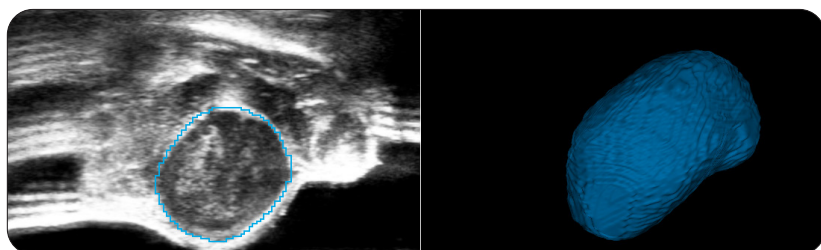
Examples of AI segmentation models using Living Image Synergy AI software.



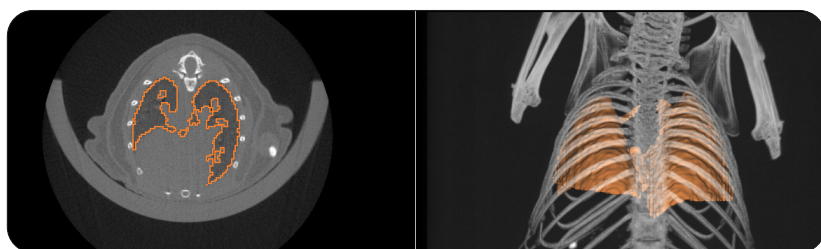
KPC 4662 tumor cells implanted into the pancreas of C57BL/6 mouse. Tumor (yellow), kidneys (orange), and spleen (blue) are shown in 2D axial cross-section (left) and 3D rendering (right). Imaged using the Vega ultrasound system.



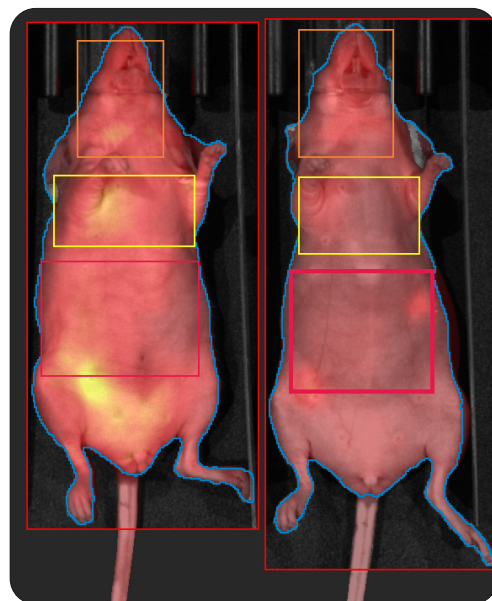
Total liver volume segmentation. Shown in 2D axial cross-section (left) and 3D rendering (right). Imaged using the Vega ultrasound system.



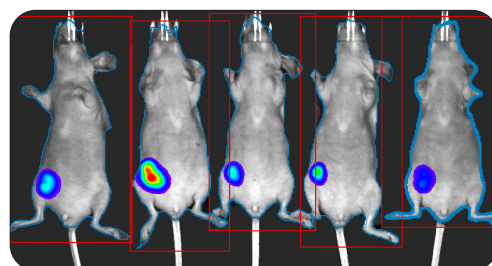
Triple negative breast cancer (TNBC) tumor cells implanted into the mammary fat pad of a female Balb/c mouse. Shown in 2D axial cross-section (left) and 3D rendering (right). Imaged using the Vega ultrasound system.



Lung air volume segmentation. Shown in 2D axial cross-section (left) and 3D rendering (right). Imaged using the Quantum GX3 microCT system.



Fluorescence imaging showing whole body contour (blue), head (orange), thoracic (yellow), and abdominal (red) segmentation regions. Imaged using the IVIS Spectrum 2 system.



Bioluminescence imaging of tumor-bearing mice with whole-body contour segmentation. Imaged using the IVIS Spectrum 2 system.

revvity