revvity

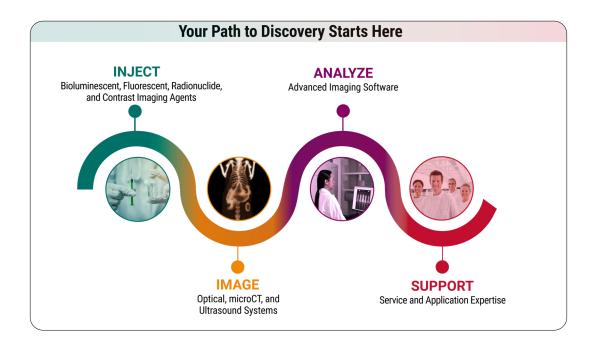
Visualize everything biology has to show you.



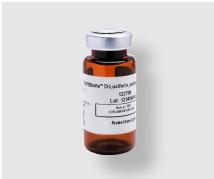
In vivo imaging solutions

Whether your research involves better understanding molecular pathways of disease, tracking disease progression, or evaluating therapeutic effectiveness of drug candidates, your cutting-edge research commands high-sensitivity and reliable *in vivo* imaging data.

Revvity is here to help you achieve your research goals with our leading molecular imaging platforms and diverse range of bioluminescent and fluorescent reagents. From single mode 2D optical and 3D tomography to multimode integrated systems, we have the tools you need to help you get the answers you are searching for to move your research forward.

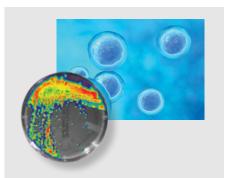


Reagents



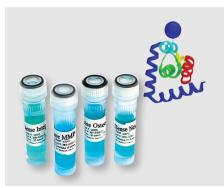
IVISbrite™ substrates

- D-Luciferin potassium salt
- D-Luciferin in RediJect™ Solution
- D-Luciferin Ultra in RediJect™ Solution
- Coelenterazine h in RediJect™ Solution



IVISbrite™ cells and bacteria

- Tumor cell lines with enhanced Red-Fluc vector
- Dual optical tumor cell lines with Red-Fluc and Green Fluorescent Protein (GFP)
- Lentiviral particles
- Bacterial strains labeled with lux operon gene



IVISense™ fluorescent probes*+

- Broad range of fluorescent probes for imaging biological targets, processes, and protease activity
 - Activatable
 - Targeted
 - Vascular
- Fluorescent panels prepackaged and targeted for your research



IVISense™ fluorescent labels and nanoparticles*

- Dyes and labels
- Nanoparticles



VesselVue® Microbubble **Contrast Agent**

P-Formulation (polydisperse)

⁺ Volume discount on selected reagents
* Select products available in sample sizes. Contact your sales representative for details.

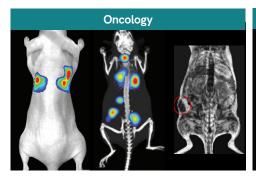
In vivo imaging instruments

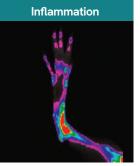
	Optical	MicroCT	Ultrasound	
	2D	2D and 3D	3D	3D
Alvaaa	rewrty	Auroau	Arvae	resorts -
IVIS® Lumina LT	IVIS Lumina S5	IVIS Spectrum 2	Quantum GX3	Vega®
Entry level BLI and FLI	High-throughput BLI and FLI	2D/3D BLI and FLI	High-resolution, low-dose	Automated, hands-free,
Leaving County	Ajma	Tevrty (microCT	high-throughput ultrasound system
IVIS Lumina III	IVIS Lumina X5	IVIS SpectrumCT 2		
IVIS Lumina III XRMS BLI and FLI with integrated x-ray	High-throughput BLI and FLI with integrated x-ray	2D/3D BLI and FLI with integrated microCT		

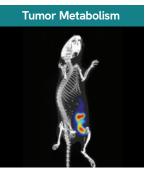
Comparison of IVIS optical imaging systems

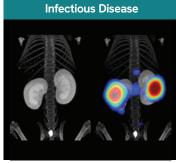
	Lumina LT	Lumina III	Lumina XRMS	Lumina S5	Lumina X5	Spectrum 2	Spectrum CT 2
Animal capacity range	1-5	1-5	1-3	1-10	1-10	1-10	1-10
2D Bioluminescence	√	√	√	√	√	√	√
2D Fluorescence	√	√	√	√	√	√	√
Near infrared fluorescence	√	√	√	√	√	√	√
Spectral unmixing (CPS)	Upgrade	√	√	√	√	√	√
Modes of fluorescent excitation	epi-illumination	epi-illumination	epi-illumination	epi-illumination	epi-illumination	epi-illumination and trans-illumination	epi-illumination and trans-illumination
2D X-ray imaging			√		√		
2D High resolution x-ray imaging					√		
3D Bioluminescence						√	√
3D Fluorescence						√	√
microCT							√

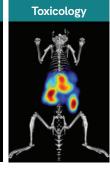


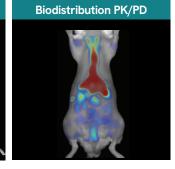


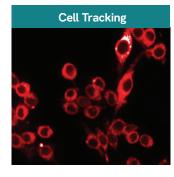


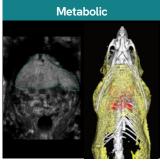




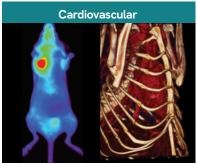


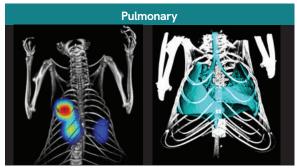












For more information on our in vivo imaging solutions visit www.revvity.com

