

# Flexible and sensitive 2D optical imaging in a benchtop format.

### Key features

- Large field of view enabling imaging of up to 10 mice\*
- Optional Smart animal handling accessories increasing throughput and streamlining imaging workflow
- Spectral unmixing for sensitive mulitspectral imaging for monitoring multiple biological events
- Full fluorescence tunability through the NIR spectrum
- High sensitivity 2D bioluminescence imaging
- Software wizard for simplified experimental workflow

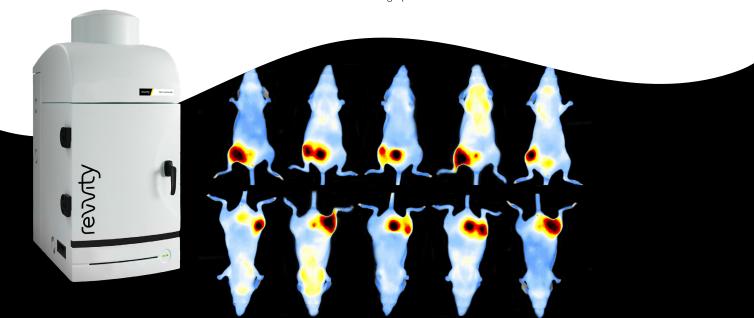
The IVIS® Lumina S5 combines best in class high throughput 2D *in vivo* bioluminescence and fluorescence imaging into a compact system that fits easily on a benchtop. With an expanded field of view for 2D optical imaging plus our unique line of accessories to accelerate setup, it has never been easier or faster to get robust data – and answers – on anatomical and molecular aspects of disease. The IVIS Lumina S5 also includes state of the art spectral unmixing features for sensitive multispectral imaging to monitor multiple biological events in the same animal.

# Improve productivity with high-throughput optical

The IVIS Lumina S5 integrates a one inch CCD camera into our benchtop instrument providing a high-throughput  $22.5 \times 22.5 \text{ cm}$  FOV sufficient for imaging 5 animals at a time with bioluminescence and fluorescence (Figure 1) or up to 10 mice with optional manifold.

**IVIS® Lumina S5** 

\*Using optional manifold



As with other IVIS Lumina systems, the IVIS Lumina S5 is equipped with 26 filters tunable to image fluorescent sources that emit from green to near-infrared. Novel illumination technology using our extended NIR range Tungsten lamp, increases fluorescent transmission through 900 nm. Moreover, the IVIS Lumina S5 incorporates our patented Compute Pure Spectrum (CPS) algorithm for spectral library generation software tools to ensure accurate autofluorescence removal, unmixing and fluorophore quantitation.

Standard on all IVIS instruments, absolute calibration affords consistent and reproducible results independent of magnification, filter selection, or acquisition settings from one instrument to any IVIS instrument within an organization or around the world.

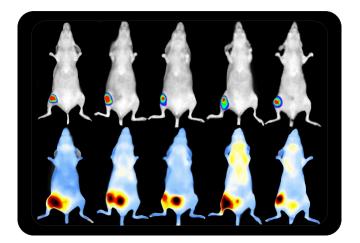


Figure 1: IVISbrite™ 4T1 Red F-luc bioluminescent tumor cells injected subcutaneously into flank and imaged on the IVIS Lumina S5. (Top) Bioluminescent signal acquisition, (bottom) fluorescence detection of IVISense™ Integrin Receptor 750 probe accumulation at tumor site.

#### Simplified workflow

Not only does the IVIS Lumina S5 offer higher throughput imaging via the one inch CCD, but it is also compatible with a set of Smart animal handling accessories (purchased separately) designed with throughput and safety in mind.

Smart loading trays (Figure 2) enable users to pose animals on the benchtop before placing the tray into the IVIS. Using fiducials built into the tray, the software can automatically recognize and draw ROIs providing automated animal identification.

Animal trays are designed with ease of use and user safety in mind. No nose cones are required thus minimizing cleanup. All tray parts have been tested and are resistant to repeated use with common laboratory disinfectants. Furthermore when used with the anesthesia unit (RAS-4), strong vacuum capabilities minimize excess gas from escaping thus preventing exposure of users to anesthetic gas.

Finally, Living Image® software brings IVIS technology to life by facilitating an intuitive workflow for *in vivo* optical analysis and data organization. The software will support input of unique animal IDs when using chip technologies and readers from third party vendors thus streamlining labeling, setup and subsequent export of data for analysis



Figure 2: Smart Animal Management accessories were designed with safety, ease of use and speed of acquisition in mind. A quick connect anesthesia port and 5 mouse/2 rat manifold come standard with the IVIS Lumina S5 (A1). The high throughput five mouse tray (purchased separately) connects seamlessly to the supplied quick connect port (a2). The ergonomic tray design does not require nose cones; new baffles securely isolate signal for data integrity; and fiducials present in the tray allow automatic subject recognition during image acquisition (B). The tray can connect either inside the IVIS (C) or it can be used for prepositioning animals (D) on the benchtop when used in conjunction with the benchtop posing station (sold separately).

www.revvity.com 2

# Improve productivity with high-throughput optical

Excitation Filter Ranges (nm)	Emission Filter Ranges (nm)	Common Dyes / Agents / Reporters
410-430	500-540	IVISense™ targeted, vascular, & activatable probes
430-450	550-590	IVISense dyes
450-470	600-640	IVISense self-quenching dyes
470-490	650-690	IVISense cell labeling dyes
490-510	690-730	Alexa Fluor® 600-750
510-530	770-810	Су5-Су7.5
530-550	825-865	DsRed, Doxorubicin**
550-570		mCherry**
570-590		tdTomato**
590-610		GFP*
610-630		FITC*
630-650		ICG
650-670		
670-690		
690-710		
710-730		
730-750		
750-770		
770-790		

<sup>\*</sup> Best used with in vitro, ex vivo and surface imaging techniques

www.revvity.com 3

<sup>\*\*</sup> Enhanced quantification with Spectral Unmixing

#### Inside the IVIS Lumina S5

- Back-illuminated, thermoelectric cooled 2.7 x 2.7 cm grade
   1 CCD provides high quantum efficiency over the entire visible to near-infrared spectrum
- · Light-tight imaging chamber
- 19 excitation filters and eight emission filters support CPS spectral unmixing
- LED lamps for photographic images
- Heated stage to maintain optimum body temperature
- Motor controlled stage, filter wheels, lens position, and f-stop

#### **Optional Accessories**

- Smart animal handling tray to accelerate setup, labeling and analysis with benchtop posing accessory
- Heavy duty workbench with integrated keyboard tray and monitor stand.
- $MVI-2^{TM}$  for 360 degree surface mapping

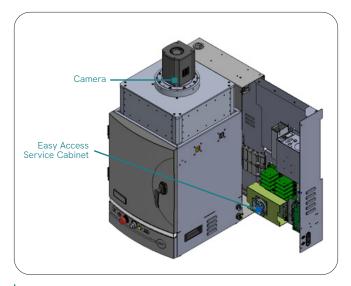


Figure 3. Diagram of external features of IVIS Lumina S5.

#### IVIS Lumina platform system accessories - Expand your IVIS Lumina Instrument with features when you need them!



Smart high throughput imaging kit CLS148874



MVI-2 Module for 360 degree surface mapping CLS143807



RAS-4 Rodent anesthesia system CLS146737

www.revvity.com

# The IVIS Lumina platform offers a selection of instruments tailored to your *in vivo* imaging needs

Features	IVIS Lumina LT	IVIS Lumina III	IVIS Lumina XRMS	IVIS Lumina S5	IVIS Lumina X5
Capacity	Up to 5 mice*	Up to 5 mice*	Up to 3 mice	Up to 10 mice**	Up to 10 mice**
1.3 x 1.3 cm CCD - 90°C	√	√	√		
2.7 x 2.7 cm CCD - 90°C				$\checkmark$	$\checkmark$
2D Bioluminescence	√	√	√	√	√
2D Fluorescence	$\checkmark$	$\checkmark$	$\checkmark$	$\checkmark$	$\checkmark$
Enhanced fluorescence		√	√	√	√
Extended range 150 W Tungsten EKE	√	√	$\checkmark$	√	$\checkmark$
Narrow bandpass excitation filters Supporting CPS Spectral Unmixing		<b>√</b>	√	√	<b>√</b>
Standard X-ray Package			$\checkmark$		
High-Resolution X-ray					√
Mouse tray with automatic subject recognition (optional)				$\checkmark$	$\checkmark$

#### Table 2. IVIS Lumina S5 Specifications

Imaging system components:     Specifications       Camera sensor     Back-illuminated thermoelectric, cooled Grade 1 CCD       Ccd size     2.7 x 2.7 cm       Ccd operating temperature     -90°C       Imaging pixels     2048 x 2048       Quantum efficiency     >85% at 500 nm, >80% at 650 nm       Pixel size     13.5 microns       Optical field of view (fov)     10 x 10, 15 x 15, 22.5 x 22.5 cm       Lens     f/1 - f/8       Minimum image pixel resolution     50 microns       Minimum read noise (e)     2e RMS for Bin 1       Dark current     < 0.0008e/pix/sec       Illumination source     Extended NIR Range 150W Tungsten EKE       Excitation fluorescence filters, number/ range     19 hard-coated narrow band pass filters       Emission fluorescence filters, number/range     7       Imaging system space requirement     48 x 69 x 122 cm (W x D x H)       Imaging chamber interior dimension     37 x 37 x 43 cm (W x D x H)       Power requirements     3.0 A max @ 100-127 Vac, 50-60 Hz (± 10%)       1.5 A max @ 200-240 Vac, 50-60 Hz (± 10%)       Stage temperature     20-40 °C       Computer (minimum specifications)     Windows* 10.4 GB RAM, pixidia Quadro 600, 250 GB and 1 TB HD 24"		
Ccd size2.7 x 2.7 cmCcd operating temperature-90°CImaging pixels2048 x 2048Quantum efficiency>85% at 500 nm, >80% at 650 nmPixel size13.5 micronsOptical field of view (fov)10 x 10, 15 x 15, 22.5 x 22.5 cmLensf/1 - f/8Minimum image pixel resolution50 micronsMinimum read noise (e)2e RMS for Bin 1Dark current< 0.0008e/pix/secIllumination sourceExtended NIR Range 150W Tungsten EKEExcitation fluorescence filters, number/ range19 hard-coated narrow band pass filtersEmission fluorescence filters, number/range7Imaging system space requirement48 x 69 x 122 cm (W x D x H)Imaging chamber interior dimension37 x 37 x 43 cm (W x D x H)Power requirements3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)Stage temperature20-40 °C	Imaging system components:	Specifications
Ccd operating temperature  -90°C  Imaging pixels  2048 x 2048  Quantum efficiency  >85% at 500 nm, >80% at 650 nm  Pixel size  13.5 microns  Optical field of view (fov)  10 x 10, 15 x 15, 22.5 x 22.5 cm  Lens  f/1 - f/8  Minimum image pixel resolution  50 microns  Minimum read noise (e)  2e RMS for Bin 1  Dark current  < 0.0008e/pix/sec  Illumination source  Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  Fmission fluorescence filters, number/range  Tmaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature	Camera sensor	Back-illuminated thermoelectric, cooled Grade 1 CCD
Imaging pixels  Quantum efficiency  >85% at 500 nm, >80% at 650 nm  Pixel size  13.5 microns  Optical field of view (fov)  10 x 10, 15 x 15, 22.5 x 22.5 cm  Lens  f/1 - f/8  Minimum image pixel resolution  50 microns  Minimum read noise (e')  2e' RMS for Bin 1  Dark current  < 0.0008e'/pix/sec  Illumination source  Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%)  1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Ccd size	2.7 x 2.7 cm
Quantum efficiency >85% at 500 nm, >80% at 650 nm  Pixel size 13.5 microns  Optical field of view (fov) 10 x 10, 15 x 15, 22.5 x 22.5 cm  Lens f/1 - f/8  Minimum image pixel resolution 50 microns  Minimum read noise (e) 2e RMS for Bin 1  Dark current < 0.0008e /pix/sec  Illumination source Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range 19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range 7  Imaging system space requirement 48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension 37 x 37 x 43 cm (W x D x H)  Power requirements 3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature 20-40 °C	Ccd operating temperature	-90°C
Pixel size  13.5 microns  Optical field of view (fov)  10 x 10, 15 x 15, 22.5 x 22.5 cm  Lens  f/1 - f/8  Minimum image pixel resolution  50 microns  Minimum read noise (e')  2e' RMS for Bin 1  Dark current  < 0.0008e'/pix/sec  Illumination source  Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Imaging pixels	2048 x 2048
Optical field of view (fov)  10 x 10, 15 x 15, 22.5 x 22.5 cm  Lens  f/1 - f/8  Minimum image pixel resolution  50 microns  Minimum read noise (e')  2e' RMS for Bin 1  Dark current  < 0.0008e'/pix/sec  Illumination source  Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Quantum efficiency	>85% at 500 nm, >80% at 650 nm
Lens f/1 - f/8  Minimum image pixel resolution 50 microns  Minimum read noise (e') 2e' RMS for Bin 1  Dark current < 0.0008e'/pix/sec  Illumination source Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range 19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range 7  Imaging system space requirement 48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension 37 x 37 x 43 cm (W x D x H)  Power requirements 3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature 20-40 °C	Pixel size	13.5 microns
Minimum image pixel resolution  50 microns  Minimum read noise (e')  2e' RMS for Bin 1  < 0.0008e'/pix/sec  Illumination source  Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Optical field of view (fov)	10 x 10, 15 x 15, 22.5 x 22.5 cm
Minimum read noise (e')  2e' RMS for Bin 1  2e' RMS for Bin 1  2e' RMS for Bin 1  200008e'/pix/sec  Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Lens	f/1 - f/8
Dark current < 0.0008e*/pix/sec  Illumination source Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range 19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range 7  Imaging system space requirement 48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension 37 x 37 x 43 cm (W x D x H)  Power requirements 3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature 20-40 °C	Minimum image pixel resolution	50 microns
Extended NIR Range 150W Tungsten EKE  Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Minimum read noise (e <sup>-</sup> )	2e <sup>-</sup> RMS for Bin 1
Excitation fluorescence filters, number/ range  19 hard-coated narrow band pass filters  Emission fluorescence filters, number/range  7  Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Dark current	< 0.0008e <sup>-</sup> /pix/sec
Emission fluorescence filters, number/range  7 Imaging system space requirement  48 x 69 x 122 cm (W x D x H) Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  Power requirements  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Illumination source	Extended NIR Range 150W Tungsten EKE
Imaging system space requirement  48 x 69 x 122 cm (W x D x H)  Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Excitation fluorescence filters, number/ range	19 hard-coated narrow band pass filters
Imaging chamber interior dimension  37 x 37 x 43 cm (W x D x H)  3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature  20-40 °C	Emission fluorescence filters, number/range	7
Power requirements 3.0A max @ 100-127 Vac, 50-60 Hz (± 10%) 1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature 20-40 °C	Imaging system space requirement	48 x 69 x 122 cm (W x D x H)
1.5A max @ 200-240 Vac, 50-60 Hz (± 10%)  Stage temperature 20-40 °C	Imaging chamber interior dimension	37 x 37 x 43 cm (W x D x H)
	Power requirements	
Computer (minimum specifications) Windows® 10, 4 GB RAM, nVidia Quadro 600, 250 GB and 1 TR HD, 24"	Stage temperature	20-40 °C
wildescreen LED Monitor	Computer (minimum specifications)	Windows® 10, 4 GB RAM, nVidia Quadro 600, 250 GB and 1 TB HD, 24" widescreen LED Monitor
Living image® software Included with IVIS purchase	Living image® software	Included with IVIS purchase

www.revvity.com

<sup>\*</sup> Using expansion lens \*\*Using optional manifold kit

#### In Vivo Imaging Solutions

#### **MICRO-CT Ultrasound Optical** Reagents **IVIS® Lumina Series III** IVIS® Lumina S5 & X5 **Vega®** IVISbrite™ **IVIS® Spectrum 2 Series Quantum GX3** • 2D optical imaging • 2D & 3D optical Automated, • 2D optical imaging • High-resolution, Bioluminescent imaging low-dose microCT hands-free substrates, cells, and • Imaging up to 5 mice • Imaging of up to 10 lentiviral particles mice using optional • Imaging of up to 10 • Cardiac & High-throughput Optional integrated manifold mice using optional respiratory gating 3 mice imaging x-ray IVISense™ manifold Optional integrated Scan times in · Fluorescent probes, high-resolution • Fully automated, < 1 minute labels, and dyes x-ray one-click • Whole body field of co-registration with **VesselVue®** Optional Smart view **IVIS SpectrumCT** accessories to • Multiple 3D imaging Microbubble Seamlessly streamline imaging modes contrast agent for workflow co-register 3D vascular ultrasound Elastography optical and hi-res MVI-2 for (tissue stiffness) imaging microCT data automated - B-mode (soft tissue 360 degree imaging • Two powerful modes imaging) of fluorescence • 4D B-mode/M-mode excitation—epi- and (cardiac imaging) trans- illumination Acoustic angiography (microvessel networks)



