







Table 1. IVIS Lumina platform comparison.

Features	Lumina LT	Lumina III	Lumina XRMS	Lumina S5	Lumina X5
Capacity	3 mice	3 mice	3 mice	5 mice	5 mice
1.3 x 1.3 cm CCD -90 °C	√	√	√		
2.7 x 2.7 cm CCD -90 °C				√	√
2D Bioluminescence	√	√	√	√	√
2D Epifluorescence Imaging	√	√	√	√	√
Extended Range 150 W Tungsten EKE	√	√	√	√	√
Narrow Bandpass Excitation Filters Supporting CPS Spectral Unmixing		√	√	√	√
Standard X-ray Package			√		
Microfocus X-ray Source – High-Resolution X-ray					√
Animal ID Support (optional)				√	√
Mouse Tray with Automatic Subject Recognition (optional)				√	√

Table 2. IVIS Lumina S5 Specifications.

Imaging System Components:	Specifications
Camera Sensor	Back-illuminated, AR coated 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 <sup>-</sup> ) (Typical)	2e <sup>-</sup> RMS for Bin 1
Dark Current	< 0.0008e <sup>-</sup> /pix/sec
Illumination Source	Extended NIR Range 150 W Tungsten EKE
Excitation Fluorescence Filters, Number/Range	19/410 - 790 nm
Emission Fluorescence Filters, Number/ Range	7/500 - 865 nm
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
Computer (Minimum specifications)	Windows® 10, 4 GB RAM, nVidia Quadro 600, 250 GB and 1 TB HD, 24" widescreen LED Monitor
Living Image® Software	1 acquisition copy and 4 analysis copies of Living Image® software version 4.0 and higher

For more information, please visit our website at [www.perkinelmer.com/invivo](http://www.perkinelmer.com/invivo)

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

67658 (S5)

PKI

DRAFT