

2D OPTICAL BENCHTOP SYSTEMS

2D AND 3D OPTICAL TOMOGRAPHY SYSTEMS

IVIS SYSTEM COMPARISON	IVIS LUMINA LT	IVIS LUMINA III	IVIS LUMINA XRMS	IVIS LUMINA S5	IVIS LUMINA X5	IVIS SPECTRUMBL	IVIS SPECTRUM	IVIS SPECTRUMCT
Detector Type	0.5" Back-thinned, back-illuminated Grade 1 CCD	0.5" Back-thinned, back-illuminated Grade 1 CCD	0.5" Back-thinned, back-illuminated Grade 1 CCD	1" Back-thinned, back-illuminated Grade 1 CCD	1" Back-thinned, back-illuminated Grade 1 CCD	1" Back-thinned, back-illuminated Grade 1 CCD	1" Back-thinned, back-illuminated Grade 1 CCD	1" Back-thinned, back-illuminated Grade 1 CCD
Camera Temp	(-90 °C)	(-90 °C)	(-90 °C)	(-90 °C)	(-90 °C)	(-90 °C)	(-90 °C)	(-90 °C)
CCD Size	1.3 cm X 1.3 cm	1.3 cm X 1.3 cm	1.3 cm X 1.3 cm	2.7 cm X 2.7 cm	2.7 cm X 2.7 cm	2.7 cm X 2.7 cm	2.7 cm X 2.7 cm	2.7 cm X 2.7 cm
CCD Pixel Size (um)	13	13	13	13.5	13.5	13.5	13.5	13.5
Optical FOV (cm)	5 - 12 cm (2.6 cm with ZFOV attachment)	5 - 12 cm (2.6 cm with ZFOV attachment)	5 - 12 cm (2.6 cm with ZFOV attachment)	10-20 cm	10-20 cm	4 - 22.5 cm	4 - 22.5 cm	4 - 22.5 cm
Optical Resolution (Microns)	50 (35 with ZFOV)	50 (35 with ZFOV)	50 (35 with ZFOV)	50	50	20	20	20
Imaging Pixels	1024 x 1024	1024 x 1024	1024 x 1024	2048 x 2048	2048 x 2048	2048 x 2048	2048 x 2048	2048 x 2048
Wavelength Range (nm)	415-875	410 - 860	410 - 860	410 - 860	410 - 860		415 - 850	415 - 850
Number of Emission Filters/ Bandwidth	4/Variable	7/30 nm	7/30 nm	7/30 nm	7/30 nm		18/20 nm	18/20 nm
Emission Filters	515-575 nm, 575-650 nm, 695-770 nm and 810-875 nm	520, 570, 620, 670, 710, 790, 845	520, 570, 620, 670, 710, 790, 845	520, 570, 620, 670, 710, 790, 845	520, 570, 620, 670, 710, 790, 845		500, 520, 540, 560, 580, 600, 620, 640, 660, 680, 700, 720, 740, 760, 780, 800, 820, 840	500, 520, 540, 560, 580, 600, 620, 640, 660, 680, 700, 720, 740, 760, 780, 800, 820, 840
Number of Excitation Filters/ Bandwidth	10/30 nm	19/20 nm	19/20 nm	19/20 nm	19/20 nm		10/30 nm	10/30 nm
Excitation Filters	430, 465, 500, 535, 570, 605, 640, 675, 710, 745	420, 440, 460, 480, 500, 520, 540, 560, 580, 600, 620, 640, 660, 680, 700, 720, 740, 760, 780	420, 440, 460, 480, 500, 520, 540, 560, 580, 600, 620, 640, 660, 680, 700, 720, 740, 760, 780	420, 440, 460, 480, 500, 520, 540, 560, 580, 600, 620, 640, 660, 680, 700, 720, 740, 760, 780	420, 440, 460, 480, 500, 520, 540, 560, 580, 600, 620, 640, 660, 680, 700, 720, 740, 760, 780		430, 465, 500, 535, 570, 605, 640, 675, 710, 745	430, 465, 500, 535, 570, 605, 640, 675, 710, 745
Excitation Light Source	Extended NIR Range 150 W Tungsten EKE-ER	Extended NIR Range 150 W Tungsten EKE-ER	Extended NIR Range 150 W Tungsten EKE-ER	Extended NIR Range 150 W Tungsten EKE-ER	Extended NIR Range 150 W Tungsten EKE-ER		Extended NIR Range 150 W Tungsten EKE	Extended NIR Range 150 W Tungsten EKE
Heated Chamber	20 - 40 °C	20 - 40 °C	20 - 40 °C	20 - 40 °C	20 - 40 °C	20 - 40 °C	20 - 40 °C	20 - 40 °C
Accessory Line	Zoom lens, MVI-2, expansion lens, isolation chamber, ECG, injection system, anesthesia, pull out tray, calibration tools, phantom mice	Zoom lens, MVI-2, expansion lens, isolation chamber, ECG, injection system, anesthesia, pull out tray, calibration tools, phantom mice	Zoom Lens, MVI-2, Isolation chamber, ECG, injection system, anesthesia, pull out tray, calibration tools, phantom mice	Quick connect anesthesia system, 5 mouse tray, posing station, RFID and P-chip support, MVI-2, Isolation chamber, ECG, injection system, anesthesia, calibration tools, phantom mice	Quick connect anesthesia system, 5 mouse tray, posing station, RFID and P-chip support, MVI-2, Isolation chamber, ECG, injection system, anesthesia, calibration tools, phantom mice	Isolation chamber, injection system, anesthesia, calibration tools, phantom mice, Multimodality Software and Mouse Imaging Shuttle	10 mouse manifold, injection system, Isolation chamber, anesthesia, calibration tools, phantom mice, Multimodality Software and Mouse Imaging Shuttle	Injection system, anesthesia, calibration tools, phantom mice, Multimodality Software

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



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2014-2017, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.