

Essen BioScience IncuCyte™ Specifications

	IncuCyte™	IncuCyte™ HD	IncuCyte™ FLR	IncuCyte™ EX
Operating Systems	Microsoft Windows® XP Microsoft Windows® Vista Windows® 7	Microsoft Windows® XP Microsoft Windows® Vista Windows® 7	Microsoft Windows® XP Microsoft Windows® Vista Windows® 7	Microsoft Windows® XP Microsoft Windows® Vista Windows® 7
Image Resolution	0.93µm 1280 x 1024 Pixels	0.93µm 1280 x 1024 Pixels	10X Objective : 1.49µm or 20X Objective : 0.74µm 1280 x 1024 Pixels	0.93µm 1280 x 1024 Pixels
Objective	20X	20X	20X or 10X	20X
Native Image Format	JPEG	JPEG	Phase : JPEG Fluorescence : 16-bit PNG	JPEG
Exported Image Format	JPEG, TIFF	JPEG, TIFF	JPEG, PNG, TIFF, RAW	JPEG, TIFF
Exported Movie Format	WMV, AVI, Metamorph®	WMV, AVI, Metamorph®	WMV, AVI, Metamorph®	WMV, AVI, Metamorph®
High Definition (HD) Optics	No	Yes	Yes	Optional
Image Size (typical)	150 Kbytes	400 Kbytes	Phase : 400 Kbytes Fluorescence : 1.1 Mbytes	150 Kbytes with HD: 400 Kbytes
Fluorescence Excitation/Emission	n/a	n/a	450-490 nm/500-530 nm	n/a
Internal Image Storage	500 Gbyte Hard Disk 2 Gbyte RAM	500 Gbyte Hard Disk 2 Gbyte RAM	>900 Gbyte Hard Disk 2 Gbyte RAM	>900 Gbyte Hard Disk 2 Gbyte RAM
Network Connection	Ethernet 10/100 Mbps RJ45 Connector	Ethernet 10/100 Mbps RJ45 Connector	Ethernet 10/100 Mbps RJ45 Connector	Ethernet 10/100 Mbps RJ45 Connector
Power	100-240 VAC 47-63Hz 2.0 Amps max @ 120V 1.0 Amps max @ 230V	100-240 VAC 47-63Hz 2.0 Amps max @ 120V 1.0 Amps max @ 230V	100-240 VAC 47-63Hz 2.0 Amps max @ 120V 1.0 Amps max @ 230V	100-240 VAC 47-63Hz 2.0 Amps max @ 120V 1.0 Amps max @ 230V
Controller Dimensions (H x W x D)	3.75" x 17.25" x 20.5" 9.5cm x 44cm x 52cm	3.75" x 17.25" x 20.5" 9.5cm x 44cm x 52cm	3.75" x 17.25" x 20.5" 9.5cm x 44cm x 52cm	3.75" x 17.25" x 20.5" 9.5cm x 44cm x 52cm
Microscope Dimensions (H x W x D)	8.8" x 17.7" x 18.3" 224mm x 450mm x 465mm	8.8" x 17.7" x 18.3" 224mm x 450mm x 465mm	9.9" x 17.7" x 18.3" 251mm x 450mm x 465mm	8.8" x 17.7" x 18.3" 224mm x 450mm x 465mm
Controller Weight	30 lbs 13.6 kg	30 lbs 13.6 kg	33 lbs 15.0 kg	33 lbs 15.0 kg
Microscope Weight	32.5 lbs 14.8 kg	32.5 lbs 14.8 kg	36 lbs 16.3 kg	37 lbs 16.8 kg
Operating Environmental Controller	0°C to 33°C 5% to 90% RH Non-Condensing	0°C to 33°C 5% to 90% RH Non-Condensing	0°C to 33°C 5% to 90% RH Non-Condensing	0°C to 33°C 5% to 90% RH Non-Condensing
Operating Environmental Microscope	0°C to 42°C 5% to 95% RH Non-Condensing	0°C to 42°C 5% to 95% RH Non-Condensing	0°C to 42°C 5% to 95% RH Non-Condensing	0°C to 33°C 5% to 90% RH Non-Condensing

Windows® is a registered trademark of Microsoft Corporation
MetaMorph® is a registered trademark of Molecular Devices

8000-0010-C



Essen BioScience Inc.
300 West Morgan Road.
Ann Arbor, MI 48108
www.essenbioscience.com