



# IncuCyte® ZOOM live-cell analysis system

## Technical Specifications

OPERATING SYSTEMS Windows XP / Windows Vista / Windows 7

OBJECTIVES 4X / 10X / 20X

IMAGE RESOLUTION **4X:** 3.05 um/pixel  
**10X:** 1.22 um/pixel  
**20X:** 0.61 um/pixel

EXPORTED IMAGE FORMATS JPEG, PNG, TIFF, RAW

EXPORTED MOVIE FORMATS WMV / AVI

FLUORESCENCE EXCITATION/EMISSION DUAL COLOR MODULE 4459	Channel	Excitation	Emission
	Green	440–480 nm	504–544 nm
	Red	565–605 nm	625–705 nm

INTERNAL DATA STORAGE 9 TByte RAID Hard Drive  
 12 GByte RAM

NETWORK CONNECTION Ethernet 10/100/1000 Mbps

POWER 100-240 VAC 50/60Hz  
 3.5 Amps max @ 120V  
 1.75 Amps max @240V

CONTROLLER DIMENSIONS (H X W X D) 5.5" X 17.0" X 21.5"  
 14.0 cm X 43.2 cm X 54.6 cm

CONTROLLER WEIGHT 38 lbs / 17.2 Kg

GANTRY DIMENSIONS (H X W X D) 12.4" X 17.7" X 18.5"  
 31.5 cm X 45.0 cm X 47.0 cm

GANTRY WEIGHT 44 lbs / 20 Kg

CONTROLLER OPERATING 0° to 33°C  
 ENVIRONMENTAL CONDITIONS 5% to 90% RH Non-Condensing

MICROSCOPE OPERATING 0° to 42°C  
 ENVIRONMENTAL CONDITIONS 5% to 95% RH Non-Condensing