



# Technical Specifications

## IncuCyte® S3 Live-Cell Analysis System

**Operating Systems:** 64-bit Windows 7 / Windows 8 /  
Windows 8.1 / Windows 10

**Objectives:** 4X / 10X / 20X

**Image Resolution:** 4X: 2.82 um/pixel  
10X: 1.24 um/pixel  
20X: 0.62 um/pixel

**Exported Image Formats:** JPEG, PNG, TIFF,  
MetaMorph Multi Dimensional File v2.0,  
Meta Imaging Series TIFF (Single / Multi-Plane) v1

**Exported Movie Formats:** WMV / AVI / MPEG-4

Fluorescence Excitation/Emission:	Channel	Excitation	Emission
	Green	440–480 nm	504–544 nm
	Red	565–605 nm	625–705 nm

**Internal Data Storage:** 16.4 TByte RAID Hard Drive  
48 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 22.7"  
14.0 cm X 43.2 cm X 57.7 cm

**Controller Weight:** 41 lbs / 18.6 Kg

**Microscope Dimensions (H X W X D):** 12.6" X 17.7" X 18.8"  
32.0 cm X 45.0 cm X 47.8 cm

**Microscope Weight:** 52 lbs / 23.6 Kg

**Controller Operating Environmental** 0° to 33°C  
**Conditions:** 5% to 90% RH Non-Condensing

**Microscope Operating Environmental** 0° to 42°C  
**Conditions:** 5% to 95% RH Non-Condensing

