



Technical Specifications

IncuCyte® S3 Live-Cell Analysis System for Neuroscience

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 |
|-----------------------------------|---------|------------|------------|
| | Orange | 513–568 nm | 577–641 nm |
| | NIR | 648–674 nm | 684–771 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 Conditions: 0° to 33°C
5% to 90% RH Non-Condensing

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

