

# Microscopy Specifications for NIR Imaging

---

## Zeiss Deconvolution Microscope

### Zeiss Specs

Quantity	Catalog Number	Description
1	423622	Shutter High Speed f/ Imager
4	424931	Reflector Module FL EC P&C
1	4880019901	Filterset 01 shift free
1	48801	FL Filter Set 10, Ex BP450-490 Shift Free
1	4880209901	Filterset 20 shift free
1	4108121999188	41037 LI-COR F/IR Dye 800
1	4326069901	Power Supply N XBO 75
1	0000001007981	Lamphousing N XBO 75 w/o Ignt Devc
1	0000001007976	Collector Lens HBO/XBO f/Lamp Z+N
1	380053987	Burner Xenon XB075 W/2 Ofr
1	423631	Adjusting Aid HBO/XBO f/ Axio Imager
1	4300049901	Axio Imager M1 Stand Mot
1	430051	Z-drive Operation Flat Rgt Imgr/Obsrvr
1	424505	Nosepiece 6x H/DIC Mot M27 f/Imager
1	4239019901	Transmitted Light Mot f/Imager
1	4236219901	Shutter for Transmitted Light f/Imager
1	42830119901	Filter Wheel Excitation 2x Mot f/Imager
1	487931	Neutral Density Filter Set D f/Imager
1	425512	Binoc Phototube Ergo 20/23 100:0/0:100
2	455043	Eyepiece W PL 10x/23 Focusing
2	444801	Folding Eyecup
1	4320019901	Mechanical stage 75x50/240 R
1	432308	Specimen Holder 76x26, immersion
1	428305	Filter Carrier 3x for filter d=32 Imager
1	467852	Conversion Filter CB 3, D=32
1	434304	Dust Protection Set L

### Halogen Illuminator

Quantity	Catalog Number	Description
1	423	Lamphousing 12V/100W w/Collector
2	380079954	Bulb HAL 12V/100W Rectangular Filament

### Epi-Fluorescence Equipment

Quantity	Catalog Number	Description
1	423603	FL/HD Light Train Man f/ Imager A,M

## Camera Adapter

Quantity	Catalog Number	Description
1	426114	C-Mount Adapter 60N C 1" 1.0x
1	4265099901	AxioCam MRm Rev. 3 with FireWire
1	41013003	AV Rel 4.6 Software

## Zeiss Optical Equipment

Quantity	Catalog Number	Description
1	42064099	Plan-Apochromat 10x/0.45 M27 D=2.0
1	4206509901	Plan-Apochromat 20x/0.8 M27 D=0.55
1	42076298	Plan-Apochromat 40x/1.3 Oil DIC UVVis-IR
1	42079099	Plan-Apochromat 100x/1.4 Oil D=0.17 M27
1	4242019901	Condsr Achr Apl 0.9 H/D/Ph/DIC Mot

## Zeiss Camera Adapter

Quantity	Catalog Number	Description
1	0000001222047	AV4 Mod MC/Z/T
1	0000001235922	AV4 Mod 3D Deconvolution
1	4108809559	Zeiss 30A Deluxe W/S SOLD W/AV
1	4108809566	ViewSonic 22" Widescreen LCD Monitor

## Chroma Filters

Quantity	Catalog Number	Description
1	HQ830/50M	
1	HQ760/40X	
1	790DCXXR	

**LI-COR**

5 05101005

[www.licor.com](http://www.licor.com)