

Microscopy Specifications for NIR Imaging

Many labs are interested in obtaining a microscope for optimal detection of infrared dyes. At LI-COR, we use microscopes with the following specifications in order to optimally image IRDye® 800CW and IRDye 680RD. These specifications are meant to provide you with a starting point in your infrared optimization discussions with your microscope manufacturer.

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 lgnt 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

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
1	HQA30/50M
1	HQ760/40X
1	790DCXXR

Contact us: biosales@licor.com
www.licor.com/microscopy
982-14682 8/14

