



Microscopy Specifications for NIR Imaging

Many labs are interested in obtaining a microscope for optimal detection of near-infrared (NIR) dyes. At LI-COR, we use microscopes compatible 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 NIR optimization discussions with your microscope manufacturer. Information is provided for the Olympus IX83 Inverted Microscope and the Olympus IX73 Inverted Microscope.

Olympus IX83 Inverted Microscope

Quantity	Catalog Number	Description
1	IX83P2F-1	IX83P2ZF-1-2; MICROSCOPE FRAME WITH 2 DECKS
1	3-U148	U-TBI90; TILTING BINOCULAR HEAD FOR GX/IX/IX2, 35-85 DEGREES
1	2-U1007	WHN10X-1-7; 10X EYEPIECE FOR BX, IX HI EYEPNT, RETSHELF, FN:22
1	2-U100H6	WHN10X-H-1-7; 10X EYEPIECE, FOC. BX, IX HI EPNT, RET SHELF, FN:22
1	U-I120	IX3-CBH; CONTROL BOX FOR IX83
1	U-R944	U-MCZ-1-2; FOCUS CONTROL MODULE FOR BX63/IX83
1	U-B30200	U-IFCBL200; CONNECTION CABLE FOR IX3/BX3/STM7 - 200CM
1	U-B30015	U-IFCBL15; CONNECTION CABLE FOR IX3/BX3 - 15CM
3	UYCP-11	UYCP-11; US STYLE 3-PRONG POWER CORD
1	Z-81226	IMMERSION OIL 30CC, LOW AUTO-FL SET W/SDS
1	OMT-010	MT-010; LOGO CLOTH DUSTCOVER 11x25x26", ANTI-STATIC BX, IX

Manual Stage and Inserts

Quantity	Catalog Number	Description
1	4-U320	IX3-SVR; RIGHT HANDLE STAGE
1	4-U324	IX3-HOW-2; WELL PLATE HOLDER
1	4-U321	IX3-H035D; 35MM DISH HOLDER
1	4-U323	IX3-HOS; SLIDE HOLDER

Objectives

Quantity	Catalog Number	Description
1	1-U2C524	UPLFLN10X2PH; U PLAN FL 10X PHASE OBJ.; PH1, NA 0.3, WD 10MM
1	1-U2C375	LUCPLFLN20XPH; LWD UPLAN FLUOR 20X PH OBJ, Ph1, NA 0.45, WD 6.6-7.8MM, W/CC
1	1-U2C377	LUCPLFLN40XPH; LWD UPLAN FLUOR 40X PH OBJ, Ph2, NA 0.6, WD 3.0-4.2MM, W/CC
1	1-U2B825	UPLSAPO20X; U PLAN S-APO 20X OBJECTIVE, NA 0.75, WD 0.6MM
1	1-U2B828	UPLSAPO40X2; U PLAN S-APO 40X OBJ, NA 0.95, WD 0.18MM, W/CC
1	1-U2B832	UPLSAPO60X0; U PLAN S-APO 60X OIL OBJ, NA1.35, WD 0.15MM, BFP1

Transmitted Light Pathway and Contrast Optics

Quantity	Catalog Number	Description
1	5-UL1237	U-LH100L-3-7;12V/100W HALOGEN LAMPHOUSE, 0.8M LONG CORD, WEEE
1	5-UL705	1 U-RMT-1-5; EXTENSION CABLE FOR HAL LAMPHOUSE, 2M
1	6-U211	IX2-LWUCD; LONG WD DIC/PHASE CONDENSER, NA0.55, WD27MM
2	8-C406	JC12V100WHAL-L; 12V, 100W HALOGEN BULB, BHS, AH2/3, U-LH, AX-LH
1	9-B323	43IF550W45; 45MM GREEN INTERFERENCE 550NM FILTER
1	9-U115	45-LBD-IF; 45MM LIGHT BALANCING DAYLIGHT FILTER
1	U-CD405	IX-PH1; PH1 PHASE ANNULUS FOR IX2-LWUCD, 30MM
1	U-CD406	IX-PH2; PH2 PHASE ANNULUS FOR IX2-LWUCD, 30MM
1	U2-CD412	IX2-DIC20; NOMARSKI PRISM FOR 20X OBJECTIVE, 30MM, IX2-LWUCD
1	U2-CD414	IX2-DIC40; NOMARSKI PRISM FOR 40X OBJECTIVE, 38MM, IX2-LWUCD
1	U2-CD416	IX2-DIC60; NOMARSKI PRISM FOR 60X OBJECTIVE, 38MM, IX2-LWUCD
1	U-M636I	IX3-FDICT; ANALYZER CUBE FOR DIC W/IX3-RFACS
1	U-P106	U-DICTS; TRANSMITTED DIC PRISM SLIDER, SHIFT TYPE

Fluorescence Pathway

Quantity	Catalog Number	Description
1	5-UR404	1 IX3-RFAL; L-SHAPED ILLUMINATOR
1	OSI-LB-XL	LB-XL; LAMBDA XL LT SOURCE ONLY, REQ LLG
1	OSI-LLG/5XL	LLG/5XL; LLG FOR LB-XL, 2M LONG, 5MM D, W LENS, TUBE
1	OSI-LG-Y73LF	LG-Y73-LF; LIQUID LIGHT GUIDE ADAPTER F/ IX73/83, MOUNTS TO RFA
1	OSI-IQ25-XL	IQ25-XL; 25MM SMARTSHUTTER FOR LAMBDA XL
1	OSI-LB-SC	LB-SC; SMART SHUTTER CONTROLLER, USB CONTROL

Fluorescence Filters

Quantity	Catalog Number	Description
1	OCT-49000BX3	49000-BX3; ET-DAPI 350/50X, BS400, 460/50M
1	OCT-49002BX3	49002-BX3; ET-EGFP/FITC/CY2 470/40X, BS495, 525/50M
1	OCT-49004BX3	49004-BX3; ET-CY3/TRITC 545/25X, BS565, 605/70M
1	OCT-49006BX3	49006-BX3; ET-CY5 620/60X, BS660, 700/75M
1	OCT-49007BX3	49007-BX3; ET-CY7 710/75X, BS760, 810/90M

Cameras and Adapters

Quantity	Catalog Number	Description
1	OH-FLASH4V3	C13440-20CU-KIT; Flash4.0 V3 sCMOS, USB 3.0, 40 FPS
1	OH-CAMRA-1007	CAMRA-1007-000-KIT; Firebird CamLink Board, Full Height Bracket, 2 SDR-SDR Cables
1	U-V360	IX3-TV1XTL; T-MOUNT ADAPTR FOR IX83/73 & LARGE FORMT CAMERAS
1	U-V411	U-CMT; C MOUNT ADAPTER FOR IX OR T-MOUNT

Olympus cellSens Software

Quantity	Catalog Number	Description
1	CS-S-CCD	CS-S-CCD-V1; HIGH END CAMERA
1	CS-DI-V1	CS-DI-V1; CELLSSENS DIMENSION - VERSION 1
1	CS-S-DE	CS-S-DE-V1; 3D DECONVOLUTION

Computer and Accessories

Quantity	Catalog Number	Description
1	D-M27FPW3	3000012208656.1; DELL ULTRASHARP 27-INCH U2717D
1	HP-Z440WIN10-1	Z8W31UC; Z440 Tower, Xeon E5-1620v3 3.5GHz, 32GB DDR4-2400, 512GB SATA SSD, 2x1TB SATA HDD, Quadro K620 2GB

Olympus IX73 Inverted Microscope

Quantity	Catalog Number	Description
1	IX73P2F	IX73P2F; MICROSCOPE FRAME WITH 2 DECKS
1	U-R380	IX3-D6RES; 6-POSITION IX NOSEPIECE CODED, DIC
1	3-U148	U-TBI90; TILTING BINOCULAR HEAD FOR GX/IX/IX2, 35-85 DEGREES
1	2-U1007	WHN10X-1-7; 10X EYEPIECE FOR BX, IX HI EYEPNT, RET SHELF, FN:22
1	2-U100H6	WHN10X-H-1-7; 10X EYEPIECE, FOC. BX, IX HI EPNT, RET SHELF, FN:22
1	UYCP-11	UYCP-11; US STYLE 3-PRONG POWER CORD
1	Z-81226	IMMERSION OIL 30CC, LOW AUTO-FL SET W/SDS
1	OMT-010	MT-010; LOGO CLOTH DUSTCOVER 11x25x26", ANTI-STATIC BX, IX

Manual Stage and Inserts

Quantity	Catalog Number	Description
1	4-U320	IX3-SVR; RIGHT HANDLE STAGE
1	4-U324	IX3-HOW-2; WELL PLATE HOLDER
1	4-U321	IX3-HO35D; 35MM DISH HOLDER
1	4-U323	IX3-HOS; SLIDE HOLDER

Objectives

Quantity	Catalog Number	Description
1	1-U2C524	UPLFLN10X2PH; U PLAN FL 10X PHASE OBJ.; PH1, NA 0.3, WD 10MM
1	1-U2C375	LUCPLFLN20XPH; LWD UPLAN FLUOR 20X PH OBJ, Ph1, NA 0.45, WD 6.6-7.8MM, W/CC
1	1-U2C377	LUCPLFLN40XPH; LWD UPLAN FLUOR 40X PH OBJ, Ph2, NA 0.6, WD 3.0-4.2MM, W/CC
1	1-U2B825	UPLSAPO20X; U PLAN S-APO 20X OBJECTIVE, NA 0.75, WD 0.6MM
1	1-U2B828	UPLSAPO40X2; U PLAN S-APO 40X OBJ, NA 0.95, WD 0.18MM, W/CC
1	1-U2B832	UPLSAPO60X0; U PLAN S-APO 60X OIL OBJ, NA 1.35, WD 0.15MM, BFP1

Transmitted Light Pathway and Contrast Optics

Quantity	Catalog Number	Description
1	5-UL1237	U-LH100L-3-7; 12V/100W HALOGEN LAMPHOUSE, 0.8M LONG CORD, WEEE
1	5-UL705	U-RMT-1-5; EXTENSION CABLE FOR HAL LAMPHOUSE, 2M
1	5-UT410	TH4-100-1-5; 12V/100W POWER SUPPLY, 100V INPUT, ROHS
1	6-U211	IX2-LWUCD; LONG WD DIC/PHASE CONDENSER, NA0.55, WD27MM
2	8-C406	JC12V100WHAL-L; 12V, 100W HALOGEN BULB, BHS, AH2/3, U-LH, AX-LH
1	9-B323	43IF550W45; 45MM GREEN INTERFERENCE 550NM FILTER
1	9-U115	45-LBD-IF; 45MM LIGHT BALANCING DAYLIGHT FILTER
1	U-CD405	IX-PH1; PH1 PHASE ANNULUS FOR IX2-LWUCD, 30MM
1	U-CD406	IX-PH2; PH2 PHASE ANNULUS FOR IX2-LWUCD, 30MM
1	U2-CD412	IX2-DIC20; NOMARSKI PRISM FOR 20X OBJECTIVE, 30MM, IX2-LWUCD
1	U2-CD414	IX2-DIC40; NOMARSKI PRISM FOR 40X OBJECTIVE, 38MM, IX2-LWUCD
1	U2-CD416	IX2-DIC60; NOMARSKI PRISM FOR 60X OBJECTIVE, 38MM, IX2-LWUCD
1	U-M636I	IX3-FDICT; ANALYZER CUBE FOR DIC W/IX3-RFACS
1	U-P106	U-DICTS; TRANSMITTED DIC PRISM SLIDER, SHIFT TYPE

Fluorescence Pathway

Quantity	Catalog Number	Description
1	5-UR404	IX3-RFAL; L-SHAPED ILLUMINATOR
1	OSI-LB-XL	LB-XL; LAMBDA XL LT SOURCE ONLY, REQ LLG
1	OSI-LLG/5XL	LLG/5XL; LLG FOR LB-XL, 2M LONG, 5MM D, W LENS, TUBE
1	OSI-LG-Y73LF	LG-Y73-LF; LIQUID LIGHT GUIDE ADAPTER F/ IX73/83, MOUNTS TO RFA
1	OSI-IQ25-XL	IQ25-XL; 25MM SMARTSHUTTER FOR LAMBDA XL
1	OSI-LB-SC	LB-SC; SMART SHUTTER CONTROLLER, USB CONTROL

Fluorescence Filters

Quantity	Catalog Number	Description
1	OCT-49000BX3	49000-BX3; ET-DAPI 350/50X, BS400, 460/50M
1	OCT-49002BX3	49002-BX3; ET-EGFP/FITC/CY2 470/40X, BS495, 525/50M
1	OCT-49004BX3	49004-BX3; ET-CY3/TRITC 545/25X, BS565, 605/70M
1	OCT-49006BX3	49006-BX3; ET-CY5 620/60X, BS660, 700/75M
1	OCT-49007BX3	49007-BX3; ET-CY7 710/75X, BS760, 810/90M

Cameras and Adapters

Quantity	Catalog Number	Description
1	OH-FLASH4V3	C13440-20CU-KIT; Flash4.0 V3 sCMOS, USB 3.0, 40 FPS
1	OH-CAMRA-1007	CAMRA-1007-000-KIT; Firebird CamLink Board, Full Height Bracket, 2 SDR-SDR Cables
1	U-V360	IX3-TV1XTL; T-MOUNT ADAPTR FOR IX83/73 & LARGE FORMT CAMERAS
1	U-V411	U-CMT; C MOUNT ADAPTER FOR IX OR T-MOUNT

Olympus cellSens Software

Quantity	Catalog Number	Description
1	CS-S-CCD	CS-S-CCD-V1; HIGH END CAMERA
1	CS-DI-V1	CS-DI-V1; CELLSSENS DIMENSION - VERSION 1
1	CS-S-DE	CS-S-DE-V1; 3D DECONVOLUTION

Computer and Accessories

Quantity	Catalog Number	Description
1	D-M27FPW3	3000012208656.1; DELL ULTRASHARP 27-INCH U2717D
1	HP-Z440WIN10-1	Z8W31UC; Z440 Tower, Xeon E5-1620v3 3.5GHz, 32GB DDR4-2400, 512GB SATA SSD, 2x1TB SATA HDD, Quadro K620 2GB

Contact us: biosales@licor.com
licor.com/microscopy

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com

LI-COR Distributor Network

www.licor.com/bio/distributors

Regional Offices

LI-COR Biosciences GmbH

Siemensstraße 25A
61352 Bad Homburg • Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.

St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com

LI-COR and IRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

©2018 LI-COR, Inc.



982-17462 04/18