

Microscopy Specifications for NIR Imaging

Olympus Deconvolution Microscope

Specifications based on Olympus IX81 Inverted Microscope

Quantity	Catalog Number	Description
1	IX81F5	IX81F3-5; IX81 FRAME W/ MOT Z & 6-POS. NOSEPIECE, REQS PRISM
1	U-I881	PRISM UNIT IX2-PRS1; FOR LEFT SIDEPORT, BI/SIDE/0/100%
1	H5-UL361	IX2-ILL100-1-5; ILL.PILLAR TILT NON-ROT.100W HOUSING DOVETAIL
1	H9-U115	LIGHT BALANCE DAYLIGHT FILTER 45MM HIGH QUALITY INTERFERENCE
1	H9-B323	FILTER, GREEN 43IF550, 45MM INTERFERENCE, IN MOUNT
1	H5-UL123	U-LH100L-3-5; 12V/100W HALOGEN LAMPHOUSE, LONG CORD
1	H5-UL705	U-RMT-1-5; EXTENSION CABLE FOR HAL LAMPHOUSE, 2M
2	H8-C406	12V 100W HALOGEN BULB FOR BHS OS 64623, 7724, AXIOVERT, EVA
1	H3-U245	STRAIGHT IX BINOCULAR U-B190CT WITH CENTERING TELESCOPE
2	H2-U1005	WHN10X-1-2; 10X EYEPIECE FOR BX,IX,HI EYEPNT,RETSHELF, FN22
1	Z-81225	IMMOIL-F30CC; IMMERSION OIL 30CC, LOW AUTO-FL
5	C-0100	"PROFESSIONAL" LENS PAPER 4" X 6" BOOK (100 SHEETS)
1	OMT-010	MT-010; LOGO CLOTH DUST COVER 11X25X26",ANTI-STATIC BX / IX

Objectives

Quantity	Catalog Number	Description
1	H1-U2C524	UPLFLN10X2PH; UPLAN FL 10X OBJ PH1, NA0.3, WD 10mm
1	H1-U2C375	LUCPLFLN20XPH; LWD UPLAN FL PH1, NA.45,WD6.4-7.6MM, W/CC
1	H1-U2C377	LUCPLFL40XPH; PH2 LWD 0.6NA WD2.9-4.2 W/C-COLLAR
1	H1-U2B825	BX UPLAN S-APO 20X 0.75NA WD 0.65MM
1	H1-U2B828	UPLSAPO40X2; U PLAN S APO 40X OBJ, NA 0.95, WD 0.18MM
1	H1-U2B832	UPLSAPO60XO; UPLAN SAPO 60X OIL OBJ. NA 1.35, WD 0.15MM

Condenser and Contrast Optics

Quantity	Catalog Number	Description
1	H6-U211	IX2-LWUCD; LONG WD DIC/PHASE CONDENSER NA 0.55 WD 27MM
1	U-P106	U-DICTS; TRANSMITTED DIC PRISM SLIDER, SHIFT TYPE
1	U-P230	POLARIZER FOR IX-LWUCD AND FOR IX-ULWCD
1	U-P236	IX2-AN; ANALYZER SLIDER FOR IX2
1	U2-CD412	IX2-DIC20; NOMARSKI PRISM FOR 20X, 30MM, IX2-LWUCD
1	U2-CD414	IX2-DIC40; NOMARSKI PRISM FOR 40X, 38MM, IX2-LWUCD
1	U2-CD416	IX2-DIC60; NOMARSKI PRISM FOR 60X, 38MM, IX2-LWUCD
1	U-CD405	PH1 PHASE ANNULUS FOR IX-LWUCD
1	U-CD406	PH2 PHASE ANNULUS FOR IX-LWUCD

Stage and Accessories

Quantity	Catalog Number	Description
1	H4-U201	PLAIN STAGE FOR IX2/81/71/51 IX2-SP

1	H4-U205	ATTACHABLE MECHANICAL STAGE 85x130mm FOR IX-SP 4-U200
1	AB4489	SLIDE HOLDER FOR 1"x3" SLIDES IX-HOS-CK40MVR/CK30/CKX40 STGE
1	H4-U293	IX2-GCP-SP; GLASS STAGE PLATE W/ COLOR BANDS AND OILING HOLE
1	H4-U923	PETRI DISH HOLDER IX-HOP FOR 65mm, 54mm, 35mm DIA.

Fluorescence Components

Quantity	Catalog Number	Description
1	H5-UR401	IX2-RFA; STRAIGHT FL ILLUMINATOR W/FS
1	LB-XL	SUTTER 15,000 HR LIGHT SOURCE
1	LLG-5XL	LIQUID LIGHT GUIDE FOR LB-XL 2M LONG, 5mm DIAMETER
1	LG-Y71-LF	LIGHT GUIDE LARGE FORMAT ADAPTER
1	IQ25-XL	25mm SMART SHUTTER W/ HOUSING TO FIT IN LAMBDA XL
1	LB-SC	LAMBDA SC CONTROL UNIT FOR SMART SHUTTER

Fluorescence Filters

Quantity	Catalog Number	Description
1	OSF-0007Z	BRIGHTLINE DAPI HI CONTRAST FILTER SET,EX387/11,409DM,ZPIX
1	U-N49002	ET-GFP; MAGNETRON SPUTTER COAT EX:488/20,DC495LP,EM:525/50
1	U-N49005	ET-DSRED(TRITC/CY3)EXET545/30X BS T570LP EM ET620/60M
1	U-N49006	ET-CY5; MAGNETRON SPUTTER COAT EX:620/60; DC660LP; EM:700/75
1	U-N49007	ET-CY7; MAGNETRON SPUTTER COAT EX:710/75 DC:760LP EM:810/90

Deep Cooled High QE NIR-CCD & Camera Mount

Quantity	Catalog Number	Description
1	OH-ORCA-R2	HAMAMATSU COOLED MONOCHROME INTERLINE FIREWIRE CCD CAMERA
1	U-V111	U-TV1X-2 CCD CAMERA ADAPTER 1X, REQUIRES CAMERA MOUNT
1	U-V311	U-CMAD3; C-MOUNT CAMERA ADAPTER

Olympus Motorized Components

Quantity	Catalog Number	Description
1	U-I115	IX2-UCB-2-5; UNIVERSAL CONTROL UNIT FOR IX2,INCL. BSW ROHS
1	U-IF181	U-ZPCB-2-5; Z-DRIVE CONTROL BOARD F/ U-UCB,BX61/62TRF
1	U-R914	HANDSET U-HSTR2 FOR BX2 MOTORIZED COMPONENTS
1	U-I501	IX2-SHA-5; MOTORIZED SHUTTER FOR USE ON TRANSMITTED LIGHT
1	H5-UR412A	IX2-RFACA-1-5; MOTORIZED FLUOR 6-POSITION CUBE
1	U-B126	SERIAL DIRECT CABLE,DB9 FEMALE TO FEMALE,F/ IX81&BX61, 15-FT.

Olympus Slidebook for Habanero™ Software

Quantity	Catalog Number	Description
1	O3I-SLBKPC	2656; SLIDEBOOK V.4.2 FOR PC HRDWR CONTROL & ANAL. SOFTWARE
1	O3I-DCON	SLIDEBOOK DECONVOLUTION (WIN OR MAC)
1	O3I-3D	SLIDEBOOK ADV. 3D RENDERING MODULE (WIN OR MAC)



Biosciences
www.licor.com