

# LI-COR Biotechnology Pricelist



 Germany

**LI-COR**<sup>®</sup>

# Content

	Page
<b>INSTRUMENTS &amp; ACCESSORIES</b>	
Odyssey <sup>®</sup> M	3
Odyssey <sup>®</sup> DLx	4
Odyssey <sup>®</sup> XF, D-DiGit <sup>™</sup> Gel Scanner	5
C-DiGit <sup>®</sup> Blot Scanner & WesternSure <sup>®</sup> Reagents	6
Pearl <sup>®</sup> Imager Family- Near-Infrared <i>in vivo</i> Imaging Systems	7
Pearl <sup>®</sup> Accessories	8
MPX <sup>™</sup> (Multiplexer) Blotting System	9
<b>COMPUTER SYSTEMS &amp; IMAGER TRAINING</b>	
Training for Odyssey Family Instruments & Pearl	9
<b>ANALYSIS SOFTWARE &amp; LICENSES</b>	
Application and Data Analysis Software for Odyssey Family Instruments & Pearl	9
<b>IRDye<sup>®</sup> CONJUGATES</b>	
IRDye Secondary Antibodies	10
<b>IRDye<sup>®</sup> AND VRDye<sup>™</sup> LABELING REAGENTS</b>	
IRDye NHS Esters and Maleimide	12
VRDye Secondary Antibodies	12
IRDye and VRDye Labeling Kits	13
Click Chemistry	13
<b>NORMALISATION REAGENTS, PROTEIN MARKER, PRIMARY ANTIBODIES etc.</b>	
Protein Marker	14
Streptavidin	14
IRDye Quencher	14
EMSA Reagents	14
Cell & Protein Stains	15
Normalisation Reagents, Primary Antibodies	15
<b>BrightSite<sup>™</sup> SMALL ANIMAL IMAGING AGENTS</b>	
IRDye Cancer Probes	15
IRDye Structural & Control Probes	15
<b>KITS, BUFFERS &amp; MORE</b>	
Western Kits, Stripping Buffer, Sample Buffer	16
Intercept Blocking Buffer	16
Intercept Antibody Diluents, Blot Storage	17
<b>INCUBATION BOXES</b>	17
<b>WESTERN &amp; IN-CELL WESTERN KITS</b>	18
<b>SERVICE CONTRACTS / EXTENDED WARRANTIES</b>	19
<b>SHIPPING CHARGES</b>	20
<b>CONTACT INFORMATION</b>	21
<b>STANDARD TERMS &amp; CONDITIONS OF SALE</b>	21

## Odyssey Family Instruments & Accessories

Part No.	Description	€
3340-BA 3350-BA	<b>Odyssey M (Fluorescence)- Data Integrity Bundle, Basic</b> <b>Odyssey M (Chemiluminescence/Fluorescence)- Data Integrity Bundle, Basic</b>	please inquire please inquire
contains:	Odyssey M Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training	
3340-STA 3350-STA	<b>Odyssey M (Fluorescence)- Data Integrity Bundle, Standard</b> <b>Odyssey M (Chemiluminescence/Fluorescence)- Data Integrity Bundle, Standard</b>	please inquire
contains:	Odyssey M Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Computer <small>Intel Core i5-10500 Processor, 4.5 GHz, 16 GB DDR4 RAM, 2 TB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day</small> Monitor 24" TFT Widescreen Monitor	
3340-PRE 3350-PRE	<b>Odyssey M (Fluorescence)- Data Integrity Bundle, Premium</b> <b>Odyssey M (Chemiluminescence/Fluorescence)- Data Integrity Bundle, Premium</b>	please inquire
contains:	Odyssey M Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Standard Extended Warranty Computer <small>Intel Core i5-10500 Processor, 4.5 GHz, 16 GB DDR4 RAM, 2 TB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day</small> Monitor 24" TFT Widescreen Monitor	

### ***Odyssey M Accessories***

926-18971	Microscope Slide Holder	1.631 €
926-18972	Microplate Insert	1.631 €
926-70003	Silicone Mat (19 x 26 cm)	104 €

## Odyssey Family Instruments & Accessories

Part No.	Description	€
<b>9142-BA</b>	<b>Odyssey DLx - Data Integrity Bundle, Basic</b>	<b>please inquire</b>
contains:	Odyssey DLx Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training	
<b>9142-STA</b>	<b>Odyssey DLx - Data Integrity Bundle, Standard</b>	<b>please inquire</b>
contains:	Odyssey DLx Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Computer Intel Core i5-8500 Processor, 4.1 GHz, 8 GB DDR4 RAM, 256 GB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day Monitor 24" TFT Widescreen Monitor	
<b>9142-PRE</b>	<b>Odyssey DLx - Data Integrity Bundle, Premium</b>	<b>please inquire</b>
contains:	Odyssey DLx Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Standard Extended Warranty Computer Intel Core i5-8500 Processor, 4.1 GHz, 8 GB DDR4 RAM, 256 GB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day Monitor 24" TFT Widescreen Monitor	

### ***Odyssey DLx/CLx Accessories***

926-72000	1 MP Alignment Guide (for single plate)	60 €
926-72200	6 MP Alignment Guide (for multiple plates)	69 €
926-70000	Silicone Mat (13 x 13 cm)	59 €
926-70001	Silicone Mat (22 x 22cm)	139 €
926-71000	4" Soft Roller	72 €
926-71804	Odyssey Pen 4-Pack, black fine-liner for membranes	48 €
929-97901	Mouse Nose Cone	107 €
929-97902	Rat Nose Cone	107 €

## Odyssey Family Instruments & Accessories, D-DiGit

Part No.	Description	€
<b>2802-BA</b>	<b>Odyssey XF - Data Integrity Bundle, Basic</b>	<b>please inquire</b>
contains:	Odyssey XF Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training	
<b>2802-STA</b>	<b>Odyssey XF - Data Integrity Bundle, Standard</b>	<b>please inquire</b>
contains:	Odyssey XF Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Computer Intel Core i5-8500 Processor, 4.1 GHz, 8 GB DDR4 RAM, 256 GB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day Monitor 24" TFT Widescreen Monitor	
<b>2802-PRE</b>	<b>Odyssey XF - Data Integrity Bundle, Premium</b>	<b>please inquire</b>
contains:	Odyssey XF Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Standard Extended Warranty Computer Intel Core i5-8500 Processor, 4.1 GHz, 8 GB DDR4 RAM, 256 GB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day Monitor 24" TFT Widescreen Monitor	
<b>Odyssey XF/Fc Accessories</b>		
926-40020	Odyssey Fc Imaging Tray, 5 pack	81 €
9957-046	Desiccant Replacement Kit	410 €
<b>3500-00</b>	<b>D-DiGit Gel Scanner</b>	<b>4.476 €</b>
3500-80	D-DiGit One Year Extended Warranty	343 €
3500-81	D-DiGit Two Years Extended Warranty <i>Only available at the initial D-DiGit purchase</i>	616 €
<b>3600-10</b>	<b>DiGit Duo</b>	<b>11.341 €</b>
	contains: C-DiGit Blot Scanner, D-DiGit Gel Scanner, Laptop PC	

## C-DiGit Blot Scanner

Part No.	Description	€
<b>3600-00</b>	<b>C-DiGit Blot Scanner</b>	7.488 €
contains:	C-DiGit Instrument C-DiGit Image Studio Analysis Software (1 user license) Quick Start Guide Resource CD (Tutorial, User Guide)	
3600-500	C-DiGit Image Studio Analysis Software - 1 user license	2.000 €
3600-500U	C-DiGit Image Studio Analysis Software - Upgrade Package	1.075 €
3600-501	C-DiGit Image Studio Analysis Software - additional 1 user license	157 €
3600-505	C-DiGit Image Studio Analysis Software - additional 5 user license	427 €
3600-510	C-DiGit Image Studio Analysis Software - additional 25 user license	790 €
3600-80	C-DiGit One Year Extended Warranty	525 €
3600-81	C-DiGit Two Years Extended Warranty	1.020 €
	<i>Both only available at the initial C-DiGit purchase</i>	
616-06141LCDG	Laptop PC for the C-DiGit	1.687 €

## WesternSure Chemiluminescent Reagents

Part No.	Description	€
<b>HRP Conjugates</b>		
926-80010	WesternSure Goat anti-Mouse HRP, 1.0 ml	152 €
926-80011	WesternSure Goat anti-Rabbit HRP, 1.0 ml	152 €
<b>Chemiluminescence Substrates</b>		
926-95000	WesternSure PREMIUM Chemiluminescent Substrate, 100 ml kit	464 €
926-95010	WesternSure PREMIUM Chemiluminescent Substrate, 20 ml kit	133 €
<b>Protein Marker</b>		
926-91000	WesternSure Pen	115 €
926-98000	WesternSure Pre-Stained Chemiluminescent Protein Ladder	254 €

## Pearl Instruments

Part No.	Description	€
<b>9430-01</b> contains:	<b>Pearl Trilogy Standard</b> <b>Pearl Trilogy Small Animal Imaging System</b> Pearl Imager (Rotameter) Pearl Docking Bed Components: Pearl Docking Bed and Mouse Bed Spares Kit: O-Rings (Pack of 25), O-Ring Cord Stock (10'), Nose Cone Plug (6), Nose Cones, (3), Neoprene Cord, Test Probe Crown Pearl Image Studio Analysis Software (1 User License) Pearl Accessory Pack Components: Tech Help Card, Network Cables, Power Cord, Clear PVC Tubing (10'), CPC Fitting (0.25"), Straight Fitting (1), Elbow Fitting (1), Tubing Clamps (6), F/AIR Filter Canister (1), Fan Filter (1) , Imaging Blanket, Alcohol Wipes Pearl Users Guide Pearl Imager Installation/Training Computer (min: Intel Core i5-8500, 4.1 GHz, 8 GB DDR4 RAM, 256 GB HDD, Windows 10, network card ) Monitor 24" TFT Widescreen	please inquire
<b>9430-03</b> contains:	<b>Pearl Trilogy Premium</b> Pearl Trilogy Imager Standard (9430-01) and Image Studio Software 25 User License*	please inquire
<b>9430-05</b> contains:	<b>Pearl Trilogy Premium Suite</b> Pearl Trilogy Imager Standard (9430-01) and Image Studio Software 25 User License* 9000-100 SmartFlow Anesthesia Suite 9000-140 Pearl Imaging Cart * All additional User Licenses refer to one Instrument Serial Number	please inquire
<b>9000-100</b>	<b>SmartFlow Anesthesia Suite</b>	15.143 €

## Pearl Accessories

Part No.	Description	€
9300-20	Pearl Docking Station	13.992 €
9300-21	Pearl Imaging Bed	6.533 €
9300-22	Pearl Clean Box	13.992 €
9000-140	Pearl Imaging Cart	12.619 €
9000-822	SmartFlow Rebreather, large	6.205 €
9000-802	Mouse Nose Cone, large (3 pieces)	361 €
9000-803	Mouse Nose Cone, small (3 pieces)	361 €
9000-804	Nose Cone Plugs, large (6 pieces)	239 €
9300-800	Nose Cone Plugs, small (6 pieces)	260 €
300-08568	Rat Nose Cone	102 €
9300-801	Rubber Cord for Imaging Bed (10')	211 €
9300-810	Imaging Blankets (25 Blankets)	260 €
9300-820	Charcoal Filter (8 Filters)	335 €
9300-821	HEPA-Filter Assembly	414 €
9300-830	Clean Box Cleaning Kit	414 €
	contains: # 9300-831 (Heavy Duty Optical Wipes), # 9300-832 (Optical Tissue), # 9300-833 (Optical Cleaner).	
9300-831	Heavy Duty Optical Wipes (300 wipes) Durable Technicloth wipers for use with Optical Cleaner for heavy duty cleaning of the window or entire Clean Box.	209 €
9300-832	Optical Tissue (150 tissues) Premium Grade Optical Tissue for use with Optical Cleaner for cleaning of the window or entire Clean Box.	209 €
9300-833	Optical Cleaner (3 Pack) Formula MC Glass Cleaner for use with Optical Tissue for cleaning of the window or entire Clean Box.	209 €
9300-834	Clean Box Cover Slip	532 €
9300-835	Glass Replacement Kit	1.547 €
9957-046	Desiccant Replacement Kit	410 €
9300-840	Dye Decontamination Kit contains: 200 Povidone-Iodine Pads, 400 Alcohol Prep Pads	217 €
9300-841	Povidone Iodine Refills, 400 Pads	145 €
9300-842	Alcohol Preparation Pad Refills, 400 Pads	145 €
9300-23	Organ Tray Base	3.125 €
9300-836	Disposable Organ Trays (5 pcs)	169 €
9300-837	Disposable Organ Trays (25 pcs)	644 €
9300-850	Compound Injection Clips (25 pcs)	183 €
9300-851	Compound Injection Clips (100 pcs)	525 €



## MPX Blotting System

Part No.	Description	€
	<b>MPX<sup>M</sup> (Multiplexer) Blotting System</b>	
	Helps to use your antibodies at optimised conditions. No cutting, stripping and re-probing.	
921-00000	MPX Blotting System	671 €
921-00110	MPX Clamp Nut (2 Ea.)	103 €
921-00120	MPX Membrane Cushion (Pack of 10)	71 €
921-00200	MPX Comb, One-Lane plus One Marker Lane	108 €
921-00202	MPX Comb, Two-Lane plus One Marker Lane	108 €
921-00204	MPX Comb, Three-Lane plus One Marker Lane	108 €
921-00206	MPX Comb, Four-Lane plus One Marker Lane	108 €

## Computer and Monitors

<b>616-06141</b>	<b>Computer</b>	<b>please inquire</b>
	The PC has the following min. requirements:	
	Intel Core i5-8500 Processor, 4.1 GHz, 8 GB DDR4 RAM, 256 GB HDD, NVIDIA GeForce Graphics, Gigabit Networking Card, DVD Writer, Windows 10 Pro, 3 Year Basic Warranty - Next Business Day	
616-06144T	Monitor 24" TFT Widescreen	please inquire

## Imager Training

Part No.	Description	€
9201-50	Odyssey Training/Retraining (per day)	2.500 €
9300-50	Pearl Imager Training/Retraining (per day)	2.500 €

## Empiria Studio - Data Analysis Software & Licenses

Part No.	Description	€
	<b>Empiria Studio for Odyssey DLx and XF</b>	
9141-500	Empiria Studio Analysis Software - 1 User License	299 €
9141-505	Empiria Studio Analysis Software - 5 User Licenses	1.125 €
9141-505IS	Empiria Studio Analysis Software - 5 User Licenses and Image Studio Analysis Software - 1 user license	1.375 €
9141-507	Empiria Studio Analysis Software - 10 user licenses	1.875 €
9141-510	Empiria Studio Analysis Software - 25 User Licenses	3.750 €
9141-510IS	Empiria Studio Analysis Software - 25 User Licenses Image Studio Analysis Software - 1 user license	3.850 €
9141-500U	Empiria Studio Analysis Software - Upgrade Package	99 €



**Free 120 Day Trial: [licor.com/TryEmpiria](http://licor.com/TryEmpiria)**

## IRDye conjugated Secondary Antibodies

Part No.	Description	€
<b>Secondary Antibodies for the 800 nm Channel</b>		
925-32210	IRDye 800CW Goat Anti-Mouse IgG (H+L), 0.1 mg	127 €
925-32211	IRDye 800CW Goat Anti-Rabbit IgG (H+L), 0.1 mg	127 €
925-32212	IRDye 800CW Donkey Anti-Mouse IgG (H+L), 0.1 mg	127 €
925-32213	IRDye 800CW Donkey Anti-Rabbit IgG (H+L), 0.1 mg	127 €
925-32214	IRDye 800CW Donkey Anti-Goat IgG (H+L), 0.1 mg	127 €
925-32218	IRDye 800CW Donkey Anti-Chicken IgG (H+L), 0.1 mg	127 €
925-32219	IRDye 800CW Goat Anti-Rat IgG (H+L), 0.1 mg	127 €
925-32232	IRDye 800CW Goat Anti-Human IgG (H+L), 0.1 mg	127 €
925-32411	IRDye 800CW Donkey Anti-Guinea Pig IgG (H+L), 0.1 mg	127 €
925-32280	IRDye 800CW Goat Anti-Mouse IgM ( $\mu$ Chain Specific), 0.1 mg	127 €
<b>Secondary Antibodies for the 700 nm Channel</b>		
925-68070	IRDye 680RD Goat Anti-Mouse IgG (H+L), 0.1 mg	127 €
925-68071	IRDye 680RD Goat Anti-Rabbit IgG (H+L), 0.1 mg	127 €
925-68072	IRDye 680RD Donkey Anti-Mouse IgG (H+L), 0.1 mg	127 €
925-68073	IRDye 680RD Donkey Anti-Rabbit IgG (H+L), 0.1 mg	127 €
925-68074	IRDye 680RD Donkey Anti-Goat IgG (H+L), 0.1 mg	127 €
925-68075	IRDye 680RD Donkey Anti-Chicken IgG (H+L), 0.1 mg	127 €
925-68076	IRDye 680RD Goat Anti-Rat IgG (H+L), 0.1 mg	127 €
925-68077	IRDye 680RD Donkey Anti-Guinea Pig IgG (H+L), 0.1 mg	127 €
925-68078	IRDye 680RD Goat Anti-Human IgG (H+L), 0.1 mg	127 €
925-68180	IRDye 680RD Goat Anti-Mouse IgM ( $\mu$ Chain Specific), 0.1 mg	127 €
925-68020	IRDye 680LT Goat Anti-Mouse IgG (H+L), 0.1 mg	127 €
925-68021	IRDye 680LT Goat Anti-Rabbit IgG (H+L), 0.1 mg	127 €
925-68022	IRDye 680LT Donkey Anti-Mouse IgG (H+L), 0.1 mg	127 €
925-68023	IRDye 680LT Donkey Anti-Rabbit IgG (H+L), 0.1 mg	127 €
925-68024	IRDye 680LT Donkey Anti-Goat IgG (H+L), 0.1 mg	127 €
925-68028	IRDye 680LT Donkey Anti-Chicken IgG (H+L), 0.1 mg	127 €
925-68029	IRDye 680LT Goat Anti-Rat IgG (H+L), 0.1 mg	127 €
925-68030	IRDye 680LT Donkey Anti-Guinea Pig IgG (H+L), 0.1 mg	127 €
925-68032	IRDye 680LT Goat Anti-Human IgG (H+L), 0.1 mg	127 €
925-68080	IRDye 680LT Goat Anti-Mouse IgM ( $\mu$ Chain Specific), 0.1 mg	127 €

## IRDye conjugated Secondary Antibodies

Part No.	Description	€
<b>Secondary Antibodies for the 800 nm Channel</b>		
926-32210	IRDye 800CW Goat Anti-Mouse IgG (H+L), 0.5 mg	214 €
926-32211	IRDye 800CW Goat Anti-Rabbit IgG (H+L), 0.5 mg	214 €
926-32212	IRDye 800CW Donkey Anti-Mouse IgG (H+L), 0.5 mg	214 €
926-32213	IRDye 800CW Donkey Anti-Rabbit IgG (H+L), 0.5 mg	214 €
926-32214	IRDye 800CW Donkey Anti-Goat IgG (H+L), 0.5 mg	214 €
926-32218	IRDye 800CW Donkey Anti-Chicken IgG (H+L), 0.5 mg	214 €
926-32219	IRDye 800CW Goat Anti-Rat IgG (H+L), 0.5 mg	214 €
926-32232	IRDye 800CW Goat Anti-Human IgG (H+L), 0.5 mg	214 €
926-32411	IRDye 800CW Donkey Anti-Guinea Pig IgG (H+L), 0.5 mg	214 €
926-32350	IRDye 800CW Goat anti-Mouse IgG <sub>1</sub> Specific, 0.5 mg	330 €
926-32351	IRDye 800CW Goat anti-Mouse IgG <sub>2a</sub> Specific, 0.5 mg	330 €
926-32352	IRDye 800CW Goat anti-Mouse IgG <sub>2b</sub> Specific, 0.5 mg	330 €
926-32280	IRDye 800CW Goat Anti-Mouse IgM ( $\mu$ Chain Specific), 0.5 mg	214 €
926-32202	IRDye 800CW Rabbit Anti-HRP IgG (H+L), 0.5 mg	214 €
<b>Secondary Antibodies for the 700 nm Channel</b>		
<i>IRDye 680RD conjugates are highly versatile and our recommended choice</i>		
926-68070	IRDye 680RD Goat Anti-Mouse IgG (H+L), 0.5 mg	214 €
926-68071	IRDye 680RD Goat Anti-Rabbit IgG (H+L), 0.5 mg	214 €
926-68072	IRDye 680RD Donkey Anti-Mouse IgG (H+L), 0.5 mg	214 €
926-68073	IRDye 680RD Donkey Anti-Rabbit IgG (H+L), 0.5 mg	214 €
926-68074	IRDye 680RD Donkey Anti-Goat IgG (H+L), 0.5 mg	214 €
926-68075	IRDye 680RD Donkey Anti-Chicken IgG (H+L), 0.5 mg	214 €
926-68076	IRDye 680RD Goat Anti-Rat IgG (H+L), 0.5 mg	214 €
926-68077	IRDye 680RD Donkey Anti-Guinea Pig IgG (H+L), 0.5 mg	214 €
926-68078	IRDye 680RD Goat Anti-Human IgG (H+L), 0.5 mg	214 €
926-68180	IRDye 680RD Goat Anti-Mouse IgM ( $\mu$ Chain Specific), 0.5 mg	214 €
<i>IRDye 680LT conjugates are highly sensitive; protocol optimisation might be required</i>		
926-68020	IRDye 680LT Goat Anti-Mouse IgG (H+L), 0.5 mg	214 €
926-68021	IRDye 680LT Goat Anti-Rabbit IgG (H+L), 0.5 mg	214 €
926-68022	IRDye 680LT Donkey Anti-Mouse IgG (H+L), 0.5 mg	214 €
926-68023	IRDye 680LT Donkey Anti-Rabbit IgG (H+L), 0.5 mg	214 €
926-68024	IRDye 680LT Donkey Anti-Goat IgG (H+L), 0.5 mg	214 €
926-68028	IRDye 680LT Donkey Anti-Chicken IgG (H+L), 0.5 mg	214 €
926-68029	IRDye 680LT Goat Anti-Rat IgG (H+L), 0.5 mg	214 €
926-68032	IRDye 680LT Goat Anti-Human IgG (H+L), 0.5 mg	214 €
926-68030	IRDye 680LT Donkey Anti-Guinea Pig IgG (H+L), 0.5 mg	214 €
926-68050	IRDye 680LT Goat anti-Mouse IgG <sub>1</sub> Specific, 0.5 mg	330 €
926-68051	IRDye 680LT Goat anti-Mouse IgG <sub>2a</sub> Specific, 0.5 mg	330 €
926-68052	IRDye 680LT Goat anti-Mouse IgG <sub>2b</sub> Specific, 0.5 mg	330 €
926-68080	IRDye 680LT Goat Anti-Mouse IgM ( $\mu$ Chain Specific), 0.5 mg	214 €



## IRDye Labeling Reagents & VRDye Conjugates

Part No.	Description	€
<b><i>Infrared Dyes for the 800 nm Channel</i></b>		
929-70020	IRDye 800CW Infrared Dye NHS ester (0.5 mg)	167 €
929-70021	IRDye 800CW Infrared Dye NHS ester (5.0 mg)	945 €
929-70022	IRDye 800CW Infrared Dye NHS ester (50.0 mg)	5.345 €
929-70023	IRDye 800CW Infrared Dye NHS ester (100.0 mg)	7.407 €
929-70024	IRDye 800CW Infrared Dye NHS ester (250.0 mg)	9.616 €
929-70025	IRDye 800CW Infrared Dye NHS ester (500.0 mg)	12.497 €
929-70026	IRDye 800CW Infrared Dye NHS ester (1.0 g)	16.263 €
929-80020	IRDye 800CW Maleimide (0.5 mg)	167 €
929-80021	IRDye 800CW Maleimide (5.0 mg)	945 €
929-80050	IRDye 800CW Maleimide (50.0 mg)	5.345 €
929-72020	IRDye 800RS Infrared Dye NHS ester (0.5 mg)	167 €
929-72021	IRDye 800RS Infrared Dye NHS ester (5.0 mg)	945 €
<b><i>Infrared Dyes for the 700 nm Channel</i></b>		
929-70050	IRDye 680RD Infrared Dye NHS ester (0.5 mg)	167 €
929-70051	IRDye 680RD Infrared Dye NHS ester (5.0 mg)	945 €
929-70052	IRDye 680RD Infrared Dye NHS ester (50.0 mg)	6.275 €
929-71050	IRDye 680RD Maleimide (0.5 mg)	167 €
929-71051	IRDye 680RD Maleimide (5.0 mg)	945 €
929-71010	IRDye 680LT Infrared Dye NHS ester (0.5 mg)	167 €
929-71011	IRDye 680LT Infrared Dye NHS ester (5.0 mg)	1.164 €
929-71500	IRDye 680LT Infrared Dye NHS ester (50.0 mg)	6.275 €
929-71008	IRDye 680LT Maleimide (0.5 mg)	167 €
929-71009	IRDye 680LT Maleimide (5.0 mg)	945 €
<b><i>Fluorescently labeled secondary antibodies for Odyssey M</i></b>		
926-49010	VRDye 490 Goat anti-Mouse IgG, (H + L) 0.5 mg	240 €
926-49020	VRDye 490 Goat anti-Rabbit IgG, (H + L) 0.5 mg	240 €
926-54010	VRDye 549 Goat anti-Mouse IgG, (H + L) 0.5 mg	240 €
926-54020	VRDye 549 Goat anti-Rabbit IgG, (H + L) 0.5 mg	240 €
926-54012	VRDye™ 549 Donkey anti-Mouse IgG Secondary Antibody	240 €
926-54013	VRDye™ 549 Donkey anti-Rabbit IgG Secondary Antibody	240 €
926-54014	VRDye™ 549 Donkey anti-Goat IgG Secondary Antibody	240 €
926-49012	VRDye™ 490 Donkey anti-Mouse IgG Secondary Antibody	240 €
926-49013	VRDye™ 490 Donkey anti-Rabbit IgG Secondary Antibody	240 €
926-49014	VRDye™ 490 Donkey anti-Goat IgG Secondary Antibody	240 €
926-54030	VRDye™ 549 Streptavidin	202 €
926-49030	VRDye™ 490 Streptavidin	202 €

## IRDye & VRDye Labeling Kits

Part No.	Description	€
<b><i>Infrared Labelling Kits for the 800 nm Channel</i></b>		
928-38040	IRDye 800CW High MW Protein Labeling Kit	579 €
928-38042	IRDye 800CW Low MW Protein Labeling Kit	636 €
928-38044	IRDye 800CW Microscale Protein Labeling Kit	475 €
<b><i>Infrared Labelling Kits for the 700 nm Channel</i></b>		
928-38072	IRDye 680RD High MW Protein Labeling Kit	579 €
928-38074	IRDye 680RD Low MW Protein Labeling Kit	636 €
928-38076	IRDye 680RD Microscale Protein Labeling Kit	475 €
928-38066	IRDye 680LT High MW Protein Labeling Kit	579 €
928-38068	IRDye 680LT Low MW Protein Labeling Kit	636 €
928-38070	IRDye 680LT Microscale Protein Labeling Kit	475 €
<b><i>Fluorescence Labeling Kits for other detection systems</i></b>		
928-49040	VRDye 490 High MW Protein Labeling Kit	579 €
928-49044	VRDye 490 Microscale Protein Labeling Kit	475 €
928-54040	VRDye 549 High MW Protein Labeling Kit	579 €
928-54044	VRDye 549 Microscale Protein Labeling Kit	475 €
<b><i>Click Chemistry</i></b>		
929-50000	IRDye 800CW DBCO, 0.5 mg	494 €
929-55000	IRDye 800CW DBCO, 5.0 mg	945 €
929-55050	IRDye 800CW DBCO, 50.0 mg	5.345 €
929-50005	IRDye 680RD DBCO, 0.5 mg	494 €
929-55005	IRDye 680RD DBCO, 5.0 mg	945 €
929-60000	IRDye 800CW Azide, 0.5 mg	494 €
929-65000	IRDye 800CW Azide, 5.0 mg	945 €
929-75050	IRDye 800CW Azide, 50.0 mg	5.345 €
929-60002	IRDye 800CW Alkyne, 0.5 mg	494 €
929-65002	IRDye 800CW Alkyne, 5.0 mg	945 €
929-65050	IRDye 800CW Alkyne, 50.0 mg	5.345 €
929-60005	IRDye 680RD Azide, 0.5 mg	494 €
929-65005	IRDye 680RD Azide, 5.0 mg	945 €
929-60006	IRDye 680RD Alkyne, 0.5 mg	494 €
929-65006	IRDye 680RD Alkyne, 5.0 mg	945 €

## Protein Marker, Stains

Part No.	Description	€
<b>Protein Marker</b>		
928-40000	Protein Molecular Weight Marker, 500µl (10 - 250 kDa)	240 €
928-50000	Chameleon™ Vue Pre-stained Protein Ladder, 250µl	159 €
928-60000	Chameleon Duo Pre-stained Protein Ladder, 500µl	239 €
928-90000	Chameleon Kit Pre-stained Protein Ladder, 2x 250µl	239 €
928-70000	Chameleon 700 Pre-stained Protein Ladder, 500µl	188 €
928-80000	Chameleon 800 Pre-stained Protein Ladder, 500µl	188 €
<b>Streptavidin</b>		
925-32230	IRDye 800CW Streptavidin (0.1 mg)	127 €
926-32230	IRDye 800CW Streptavidin (0.5 mg)	202 €
925-68079	IRDye 680RD Streptavidin (0.1 mg)	127 €
926-68079	IRDye 680RD Streptavidin (0.5 mg)	202 €
925-68031	IRDye 680LT Streptavidin (0.1 mg)	127 €
926-68031	IRDye 680LT Streptavidin (0.5 mg)	202 €
<b>IRDye Quencher</b>		
929-70030	IRDye QC-1 NHS Ester (0.5 mg)	313 €
929-70031	IRDye QC-1 NHS Ester (5.0 mg)	1.376 €
<b>Cell &amp; Protein Stains</b>		
929-90010	CellVue Burgundy (100 nmol, 10 ml of Diluent C)	391 €
929-90020	CellVue NIR815 (100 nmol, 10 ml of Diluent C)	391 €
929-90001	CellVue Dilluent C (6x 10 ml)	360 €
929-90030	PSVue 794 Reagent Kit	765 €
<b>EMSA Reagents</b>		
829-07921	p53 IRDye 700 (1.25 nmol)	85 €
829-07922	STAT3 IRDye 700 (1.25 nmol)	85 €
829-07923	CREB IRDye 700 (1.25 nmol)	85 €
829-07924	NFkB IRDye 700 (1.25 nmol)	85 €
829-07925	AP-1 IRDye 700 (1.25 nmol)	85 €
829-07926	Sp-1 IRDye 700 (1.25 nmol)	85 €
829-07929	HIF-1 IRDye 700 (1.25 nmol)	85 €
829-07910	Odyssey EMSA Buffer Kit (100 EMSAs) <small>Components: 10x Binding Buffer, 25 mM DTT, 2.5% Tween-20, Poly(dI-dC) Tris EDTA, Salmon Sperm DNA, 50% Glycerol, 1% NP-40, 1 M KCl, 100 mM MgCl, 200 mM EDTA (pH 8.0), 10x Orange Loading Dye</small>	316 €
927-10100	10x Orange Loading Dye (3 tubes à 1.0 ml)	90 €

## Normalisation Reagents, Primary Antibodies, Chemi IR kits

Part No.	Description	€
<b>Total Protein Stains for Normalisation</b>		
926-11010	Revert 700 Total Protein Stain Kit 100ml, incl. Stain, Wash and Reversal Solution*	155 €
926-11016	Revert 700 Total Protein Stain Kit 250ml, incl. Stain, Wash and Reversal Solution*	290 €
926-11015	Revert 700 Stain and Wash Solution Kit 100 ml, incl. Stain and Wash Solution	133 €
926-11011	Revert 700 Total Protein Stain, 100ml	115 €
926-11021	Revert 700 Total Protein Stain, 250ml	211 €
926-10010	Revert 520 Total Protein Stain Kit 200ml, incl. Stain, Wash and Reversal Solution*	178 €
926-10016	Revert 520 Total Protein Stain Kit 500ml, incl. Stain, Wash and Reversal Solution*	333 €
926-10015	Revert 520 Stain and Wash Solution Kit 200 ml, incl. Stain and Wash Solution	153 €
926-10011	Revert 520 Total Protein Stain, 200ml	132 €
926-10021	Revert 520 Total Protein Stain, 500ml	242 €
<b>Loading Indicators</b>		
926-20002	Odyssey 28kD Loading Indicator - 800nm	144 €
926-20004	Odyssey 28kD Loading Indicator - 700nm	144 €
<b>Primary Antibodies</b>		
926-42210	$\beta$ -Actin Rabbit monoclonal antibody, 0.1 ml	455 €
926-42211	$\beta$ -Tubulin Rabbit polyclonal antibody, 0.1 ml	455 €
926-42212	$\beta$ -Actin Mouse monoclonal antibody, 0.1 ml	455 €
926-42213	$\alpha$ -Tubulin Mouse monoclonal antibody, 0.1 ml	455 €
926-42214	COX IV Rabbit monoclonal antibody, 0.1 ml	455 €
926-42215	Vinculin Rabbit Monoclonal Antibody, 0.1 ml	455 €
926-42216	GAPDH Rabbit Monoclonal Antibody, 0.1 ml	455 €
926-42218	Histone H3 Mouse Monoclonal Antibody, 0.1 ml	455 €
926-42219	Histone H3 Rabbit Monoclonal Antibody, 0.1 ml	455 €

## BrightSite Small Animal Imaging Agents

Part No.	Description	€
<b>Probes</b>		
926-08446	IRDye 800CW EGF Optical Probe (20 nmol, lyophilized)	528 €
926-08946	IRDye 800CW 2-DG Optical Probe (100 nmol, lyophilized)	528 €
926-09889	IRDye 800CW RGD Optical Probe (15 nmol, lyophilized)	528 €
926-50401	IRDye 800CW PEG Contrast Agent (15 nmol, lyophilized)	528 €
926-09375	IRDye 800CW BoneTag Optical Probe (40 nmol, lyophilized)	528 €
926-09374	IRDye 680RD BoneTag Optical Probe (40 nmol, lyophilized)	528 €
<b>Control Probes</b>		
929-08972	IRDye 800CW Carboxylate (100 nmol, lyophilized)	203 €
929-09406	IRDye 800CW Carboxylate (5 mg)	829 €
929-70122	IRDye 800CW Carboxylate (50 mg)	2.504 €
929-70123	IRDye 800CW Carboxylate (100 mg)	3.254 €
929-70124	IRDye 800CW Carboxylate (250 mg)	5.239 €
929-70125	IRDye 800CW Carboxylate (500 mg)	6.837 €
929-70126	IRDye 800CW Carboxylate (1 g)	8.891 €
929-08980	IRDye 680RD Carboxylate (100 nmol, lyophilized)	203 €

## Kits, Buffers & more

Part No.	Description	€
<b><i>Protein Sample Loading Buffer</i></b>		
928-40004	4x Protein Sample Loading Buffer (15 ml)	49 €
<b><i>Western Blotting Membranes &amp; Kits</i></b>		
926-31090	Odyssey Nitrocellulose Membranes (10 sheets, 0.22 µm, 7 x 8.5 cm)	86 €
926-31092	Roll of Nitrocellulose Membranes (1 roll, 0.22 µm, 30 cm x 3 m) *	281 €
926-31097	Protein Sample Loading Buffer and PVDF Membrane Kit 4x Protein Sample Loading Buffer (15ml), 1 roll of Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 26.5 cm x 3.75 m)	559 €
829-32080	PBS Blocking Buffer and PVDF Membrane Kit, large 10x 500 ml bottles of Intercept Blocking Buffer (PBS), 1 roll of Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 26.5 cm x 3.75 m)	1.483 €
926-32100	TBS Blocking Buffer and PVDF Membrane Kit, large 10x 500 ml bottles of Intercept Blocking Buffer (TBS), 1 roll of Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 26.5 cm x 3.75 m)	1.483 €
926-32098	PBS Blocking Buffer and PVDF Membrane Kit, small 1x 500 ml bottle of Intercept Blocking Buffer (PBS), 1 roll of Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 26.5 cm x 3.75 m)	650 €
926-32099	TBS Blocking Buffer and PVDF Membrane Kit, small 1x 500 ml bottle of Intercept Blocking Buffer (TBS), 1 roll of Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 26.5 cm x 3.75 m)	650 €
<b><i>Stripping Buffer</i></b>		
926-92000	WesternSure ECL Stripping Buffer (100 ml, 5x concentrated)	159 €
928-40028	NewBlot IR Stripping Buffer (100 ml, 5x concentrated)	159 €
928-40030	New Blot™ Nitro Stripping Buffer (100 ml, 5x concentrated) *	245 €
928-40032	New Blot™ PVDF Stripping Buffer (100 ml, 5x concentrated) *	247 €
<b><i>Intercept Blocking Buffer</i></b>		
927-60001	Intercept (TBS) Blocking Buffer, 500 ml	136 €
927-60003	Intercept (TBS) Blocking Buffer, 3x 500 ml	365 €
927-60010	Intercept (TBS) Blocking Buffer, 10x 500 ml	1.073 €
927-60025	Intercept (TBS) Blocking Buffer, 25x 500 ml	2.351 €
927-60050	Intercept (TBS) Blocking Buffer, 50x 500 ml	4.108 €
927-70001	Intercept (PBS) Blocking Buffer, 500 ml	136 €
927-70003	Intercept (PBS) Blocking Buffer, 3x 500 ml	365 €
927-70010	Intercept (PBS) Blocking Buffer, 10x 500 ml	1.073 €
927-70025	Intercept (PBS) Blocking Buffer, 25x 500 ml	2.351 €
927-70050	Intercept (PBS) Blocking Buffer, 50x 500 ml	4.108 €
927-80001	Intercept (TBS) Protein-Free Blocking Buffer, 500 ml	136 €
927-80003	Intercept (TBS) Protein-Free Blocking Buffer, 3x 500 ml	365 €
927-80010	Intercept (TBS) Protein-Free Blocking Buffer, 10x 500 ml	1.073 €
927-90001	Intercept (PBS) Protein-Free Blocking Buffer, 500 ml	136 €
927-90003	Intercept (PBS) Protein-Free Blocking Buffer, 3x 500 ml	365 €
927-90010	Intercept (PBS) Protein-Free Blocking Buffer, 10x 500 ml	1.073 €



## Antibody Diluents & Incubation Boxes

Part No.	Description	€
<b>Intercept Antibody Diluents</b>		
927-65001	Intercept T20 (TBS) Antibody Diluent, 500ml	144 €
927-65003	Intercept T20 (TBS) Antibody Diluent, 3x 500ml	384 €
927-65010	Intercept T20 (TBS) Antibody Diluent, 10x 500ml	1.130 €
927-75001	Intercept T20 (PBS) Antibody Diluent, 500ml	144 €
927-75003	Intercept T20 (PBS) Antibody Diluent, 3x 500ml	384 €
927-75010	Intercept T20 (PBS) Antibody Diluent, 10x 500ml	1.130 €
927-66003	Intercept T20 (TBS) Blocking Buffer & Antibody Diluent Kit (contains: Intercept (TBS) Blocking Buffer, 500ml and Intercept (TBS) Antibody Diluent, 2x 500ml)	365 €
927-76003	Intercept T20 (PBS) Blocking Buffer & Antibody Diluent Kit (contains: Intercept (PBS) Blocking Buffer, 500ml and Intercept (PBS) Antibody Diluent, 2x 500ml)	365 €
927-85001	Intercept T20 (TBS) Protein-Free Antibody Diluent, 500ml	144 €
927-85003	Intercept T20 (TBS) Protein-Free Antibody Diluent, 3x 500ml	384 €
927-85010	Intercept T20 (TBS) Protein-Free Antibody Diluent, 10x 500ml	1.130 €
927-95001	Intercept T20 (PBS) Protein-Free Antibody Diluent, 500ml	144 €
927-95003	Intercept T20 (PBS) Protein-Free Antibody Diluent, 3x 500ml	384 €
927-95010	Intercept T20 (PBS) Protein-Free Antibody Diluent, 10x 500ml	1.130 €
<b>Blot Storage</b>		
929-95100	Western Blot Storage Bags, 40 pc	38 €
<b>Black Western Incubation Boxes</b>		
<i>also available in pink and transparent, please check our website</i>		
Small - Size: 7.2 x 4.9 x 3.0 cm		
929-97101	Western Incubation Box Small One Pack	18 €
929-97105	Western Incubation Box Small Five Pack	55 €
929-97110	Western Incubation Box Small Ten Pack	86 €
Medium - Size: 8.9 x 6.5 x 2.8 cm		
929-97201	Western Incubation Box Medium One Pack	21 €
929-97205	Western Incubation Box Medium Five Pack	61 €
929-97210	Western Incubation Box Medium Ten Pack	100 €
Large - Size: 11.5 x 8.8 x 2.8 cm		
929-97301	Western Incubation Box Large One Pack	26 €
929-97305	Western Incubation Box Large Five Pack	77 €
929-97310	Western Incubation Box Large Ten Pack	133 €
Extra Large - Size: 15.2 x 10.1 x 3.1 cm		
929-97401	Western Incubation Box Extra Large One Pack	27 €
929-97405	Western Incubation Box Extra Large Five Pack	91 €
929-97410	Western Incubation Box Extra Large Ten Pack	154 €

## Western & In-Cell Western Kits

Part No.	Description	€
<b>Western Blotting Kits</b>		
926-32081	Western Blotting Kit I RD (PBS/PVDF) IRDye 800CW Goat Anti-Mouse Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Rabbit Secondary Antibody (0.1 mg), Intercept Blocking Buffer (PBS), 500 ml, 10 Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 10 x 10 cm)	456 €
926-32082	Western Blotting Kit II RD (PBS/PVDF) IRDye 800CW Goat Anti-Rabbit Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Mouse Secondary Antibody (0.1 mg), Intercept Blocking Buffer (PBS), 500 ml, 10 Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 10 x 10 cm)	456 €
926-32083	Western Blotting Kit III RD (PBS/Nitro) IRDye 800CW Goat Anti-Mouse Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Rabbit Secondary Antibody (0.1 mg), Intercept Blocking Buffer (PBS), 500 ml, 10 Odyssey Nitrocellulose Membranes (0.22 µm, 7 x 8.5 cm)	456 €
926-32084	Western Blotting Kit IV RD (PBS/Nitro) IRDye 800CW Goat Anti-Rabbit Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Mouse Secondary Antibody (0.1 mg), Intercept Blocking Buffer (PBS), 500 ml, 10 Odyssey Nitrocellulose Membranes (0.22 µm, 7 x 8.5 cm)	456 €
926-35010	Western Blotting Kit V RD (TBS/PVDF) IRDye 800CW Goat Anti-Mouse Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Rabbit Secondary Antibody (0.1 mg), Intercept Blocking Buffer (TBS), 500 ml, 10 Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 10 x 10 cm)	456 €
926-35011	Western Blotting Kit VI RD (TBS/PVDF) IRDye 800CW Goat Anti-Rabbit Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Mouse Secondary Antibody (0.1 mg), Intercept Blocking Buffer (TBS), 500 ml, 10 Millipore ImmobilonTM-FL PVDF Membranes (0.45 µm, 10 x 10 cm)	456 €
926-35014	Western Blotting Kit VII RD (TBS/Nitro) IRDye 800CW Goat Anti-Mouse Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Rabbit Secondary Antibody (0.1 mg), Intercept Blocking Buffer (TBS), 500 ml, 10 Odyssey Nitrocellulose Membranes (0.22 µm, 7 x 8.5 cm)	456 €
926-35015	Western Blotting Kit VIII RD (TBS/Nitro) IRDye 800CW Goat Anti-Rabbit Secondary Antibody (0.1 mg), IRDye 680RD Goat Anti-Mouse Secondary Antibody (0.1 mg), Intercept Blocking Buffer (TBS), 500 ml, 10 Odyssey Nitrocellulose Membranes (0.22 µm, 7 x 8.5 cm)	456 €
<b>In-Cell Western Kits</b>		
926-42091	CellTag™ 700 Stain ICW Kit I IRDye 800CW Goat Anti Mouse Antibody (0.5 mg, lyophilized), Intercept Blocking Buffer (500 ml), CellTag 700 Stain (2x 10 nmol). Optimized for: 20 96-well plates or 10 384-well plates	411 €
926-42092	CellTag™ 700 Stain ICW Kit II IRDye 800CW Goat Anti Rabbit Antibody (0.5 mg, lyophilized), Intercept Blocking Buffer (500 ml), CellTag 700 Stain (2x 10 nmol). Optimized for: 20 96-well plates or 10 384-well plates	411 €
926-42093	GAM/GAR CellTag™ 520 Stain ICW Kit III IRDye 800CW Goat Anti Mouse and IRDye 680RD Goat anti Rabbit Antibodies (0.5 mg, lyophilized), Intercept Blocking Buffer (500 ml), CellTag 520 Stain (2x 50 nmol). Optimized for: 20 96-well plates or 10 384-well plates	492 €
926-42094	GAR/GAM CellTag™ 520 Stain ICW Kit IV IRDye 800CW Goat Anti Rabbit and IRDye 680RD Goat anti Mouse Antibodies (0.5 mg, lyophilized), Intercept Blocking Buffer (500 ml), CellTag 520 Stain (2x 50 nmol). Optimized for: 20 96-well plates or 10 384-well plates	492 €
926-41090	CellTag™ 700 Stain, 2 x 10 nmol	211 €
926-41094	CellTag™ 520 Stain, 2 x 50 nmol	296 €
928-40022	Sapphire700 Stain (100 µl)	143 €
<b>Quick Western Kits</b>		
926-69100	Quick Western Kit - IRDye 680RD IRDye 680RD Detection Reagent, Intercept Blocking Buffer	478 €
<b>Microplates</b>		
926-19156	96-well ICW Microplate, 8-pack (back with clear bottom)	126 €
926-19157	96-well ICW Microplate, 32-pack (back with clear bottom)	377 €

## Service Contracts / Extended Warranties

Part No.	Description	€
<b>Basic Extended Warranties</b>		
3340-80E	Odyssey M Basic Extended Warranty, 1 year	6.313 €
9142-80E	Odyssey DLx Basic Extended Warranty, 1 year	4.387 €
2802-80E	Odyssey XF Basic Extended Warranty, 1 year	3.210 €
9400-80E	Pearl Basic Extended Warranty, 1 year	4.387 €

The "**Basic Extended Warranty**" includes web and phone support, live software webinars, LI-COR certified spare parts, labour and travel costs.

<b>Standard Extended Warranties</b>		
3340-82E	Odyssey M Standard Extended Warranty, 1 year	6.741 €
9142-82E	Odyssey DLx Standard Extended Warranty, 1 year	4.815 €
2802-82E	Odyssey XF Standard Extended Warranty, 1 year	3.638 €
9400-82E	Pearl Standard Extended Warranty, 1 year	4.815 €

The "**Standard Extended Warranty**" includes web and phone support, live software webinars, a 10% discount code\* for reagent and accessory orders placed with LI-COR directly, LI-COR certified spare parts, labour and travel costs. (\* Multipackages Blocking buffer and incubation boxes excluded)

<b>Premium Support Contracts</b>		
3340-84	Odyssey M Premium Support Contract, 1 year	8.346 €
9142-84	Odyssey DLx Premium Support Contract, 1 year	6.420 €
2802-84	Odyssey XF Premium Support Contract, 1 year	5.029 €
9400-84	Pearl Premium Support Contract, 1 year	6.420 €

The "**Premium Support Contract**" includes web and phone support, live software webinars, a 10% discount code\* for reagent and accessory orders placed with LI-COR directly, one license of Empiria Studio Western blot analysis software, one annual on-site preventative maintenance check, LI-COR certified spare parts, labour and travel costs. (\* Multipackages Blocking buffer and incubation boxes excluded)

### Services

Preventative Maintenance	please inquire
Verify Normal Operation Test / Re-qualification for warranty or support contract coverage	please inquire

## Shipping Charges

Description	€
Standard delivery (DPD)	12,90 €
* Courier surcharge for dangerous goods delivery	40 €
<small>Parts categorised as dangerous goods (IATA) are indicated</small>	

**Free Shipping for orders above 750€ when ordered online**  
**[www.licor.com/bio/store](http://www.licor.com/bio/store)**

## Contact Information

LI-COR Biosciences GmbH  
Siemensstraße 25A

D-61352 Bad Homburg

Phone: +49 (0) 6172-1717771

Fax: +49 (0) 6172-1717799

Email: [bio-eu@licor.com](mailto:bio-eu@licor.com)

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

V14.12.2022 (Valid as of 01.01.2023)

©2023 LI-COR inc. LI-COR, Odyssey, C-DiGit, Pearl, IRDye, VRDye, BrightSite, CellTag, WesternSure, Image Studio, Empiria Studio, Intercept

are trademarks or registered trademarks of LI-COR, inc.

All other trademarks belong to their respective owners and are used by permission.

SmartFlow Anesthesia System: LI-COR in cooperation with Euthanex Corporation.

### All information without engagement.

## Standard Terms & Conditions of Sale

### § 1. General remarks, scope

(1) Unless agreed otherwise in writing in the individual case, the following terms of business shall apply to all current and future business relations between ourselves and the purchaser.

(2) We do not effect legal transactions with private consumers. Consequently, these General Terms of Business have no validity in respect of transactions with private end-users.

(3) Even if known, any departing, conflicting or supplementary General Terms of Business shall not form part of the contract unless their validity is expressly agreed in writing.

### § 2. Conclusion of contract

(1) Our quotations are always non-binding.

(2) By ordering goods the customer makes a binding declaration that he wishes to acquire the goods ordered. We shall be entitled to accept the contract offer inherent in the order within two weeks of receipt by us by issuing a written order confirmation. The supply contract shall not exist until we have issued a written order confirmation or the order has been executed.

(3) If the customer orders the goods electronically we will confirm receipt of the order without delay. The confirmation of receipt shall not form an order confirmation. The confirmation of receipt may be combined with the order confirmation.

(4) The contract is concluded subject to the reservation that we are supplied correctly and punctually by our own suppliers. This shall only apply in the event that we are not responsible for the non-delivery.

(5) The customer will be informed without delay if it is not possible to execute the order. Any payment will be reimbursed without delay.

(6) If the customer orders the goods electronically the text of the contract will be stored by us and sent to the customer electronically (by e-mail) on request along with the present General Terms of Business.

### § 3. Delivery and payment

(1) We will always endeavour to comply with the delivery dates and deadlines stated in the order confirmation. Nonetheless, details of delivery and performance dates are provided for guidance only and are non-binding on us, unless we have given an express assurance in writing that a delivery date is binding. If a binding assured delivery date is not met for reasons for which the customer is responsible, e.g. incomplete or non-punctual performance of his contractual obligations, in particular incomplete and non-punctual forwarding of all necessary approvals, documentation, samples and drawings, then the assertion against us of any claim for damages shall be excluded.

(2) If the failure to meet the deadline is due to unforeseeable events for which we are not responsible, e.g. strike, measures by authorities, operating or transport disruption, fire or water damage, war, energy or raw materials shortage or force majeure, we shall be released from compliance with the delivery dates and prices.

## Standard Terms & Conditions of Sale

(3) Should we be in default of delivery for reasons for which we are responsible, the customer may only demand payment of damages instead of performance if he has set us a reasonable subsequent period in which to perform the service and this subsequent period has expired without result.

(4) If performance or delivery is delayed or cancelled at the customer's request or for reasons for which the customer is responsible we shall be entitled to claim any expenses that we have thereby incurred from the customer in damages and to charge the customer for costs incurred, e.g. storage charges, as appropriate.

(5) The scope of supply shall be determined by our written order confirmation.

(6) We reserve the right to make technical changes – in particular construction and design changes – based on improvements in technology or changed legislative requirements within the confines of what is reasonable, even after conclusion of the contract, provided that the fitness of the object supplied for the intended purpose is not thereby impaired. The same shall apply to changes in the scope of supply.

(7) Partial deliveries shall be made only with the express consent of the customer; the higher carriage costs thereby incurred will be invoiced in addition.

### § 4. Packaging and shipping, transfer of risk

(1) Packing and the choice of packaging materials shall be our responsibility. The choice of shipping method shall be at our discretion. If the customer wishes express carriage, however, he shall pay the additional costs thereby incurred.

(2) Performance and delivery will be effected with the usual level of due care unless the customer has given very specific instructions when placing or confirming the order. We shall not be liable for damage arising therefrom.

(3) Should supply include erection and assembly, the risk shall transfer to the customer at the time of acceptance.

### § 5. Prices and payment terms

(1) The selling price quoted is for delivery ex works excluding statutory value-added tax in each case. Supply and invoicing shall be carried out at the list prices (plus value-added tax and carriage costs) and conditions effective on the shipping date.

(2) Our prices are always calculated in euros, the price in euros being binding even if payment of the purchase price in a currency other than euros is agreed. Changes in the exchange rate with the euro which occur after the contract has been concluded relative to the exchange rate on the date the contract is concluded shall be for the purchaser's account.

(3) The purchase price shall be payable immediately following issue of the invoice, within 30 days without deduction.

(4) On expiry of 30 days from issue of the invoice the customer shall be in arrears. Whilst he is in arrears the customer shall pay interest on the debt at 8% above the interest base rate. We reserve the right to substantiate and claim greater damages for the arrears in individual cases.

(5) The customer shall be entitled to offset amounts only if his counter-claims have been legally established or recognised by us.

(6) Should we become aware following conclusion of the contract of circumstances which suggest that the customer may be unable to pay, we shall be entitled to refuse to supply the goods and to set the customer a reasonable period in which to pay the purchase price or provide adequate security. In this event supply of the goods purchased shall take place step by step with payment of the purchase price or provision of the security.

(7) If the customer is in arrears with a payment or there are grounds to doubt his ability to pay we shall be entitled to declare any deferred invoice amounts due immediately.

(8) If payment of the purchase price in instalments has been agreed we shall be entitled to declare the entire outstanding purchase price due immediately should the customer be in arrears with the payment of an instalment.

### § 6. Reservation of ownership

(1) We reserve right of ownership of the goods sold until full settlement of all claims from a current business relationship.

(2) The customer shall bear the risk for the goods supplied by us and shall be obliged to insure them adequately against fire, water and theft risks at replacement value.

(3) The customer shall be obliged to inform us without delay of any access to the goods by third parties, for example in the event of attachment, and of any damage to or destruction of the goods. The customer shall notify us without delay of any change in ownership of the goods or any change in his own address. The customer shall be liable for the costs arising from any access by third parties to the goods.

(4) In the event of breach of contract by the customer, in particular in the event of default on payment or infringement of one of the above obligations, we shall be entitled to withdraw from the contract and to demand return of the goods.

### § 7. Warranty

(1) Our warranty for defects in the goods supplied by us shall initially provide for repair or replacement at our discretion.

(2) If the remedial measures are unsuccessful the customer may at his discretion demand a reduction in payment (reduction) or annulment of the contract (rescission). However, the customer shall have no right of rescission in the event of minor infringement of contractual liabilities on our part, in particular in the case of defects which reduce the value of the object only insignificantly.

(3) Obvious defects must be indicated to us in writing within a period of two weeks of receipt of the goods; otherwise the assertion of any warranty claim shall be excluded. Dispatch within this period shall be sufficient to comply with the deadline. The customer shall bear the full burden of proof for all claim conditions, in particular the defect itself, the date on which the defect was found and the punctuality of the complaint.

(4) If the customer chooses rescission from the contract following failure to remedy a defect he shall have no further claim to damages on account of the defect.

(5) If the customer chooses compensation following failure of remedial work the goods shall remain with the customer if this is reasonable for him. The compensation shall be limited to the difference between the purchase price and the value of the defective object. This shall not apply if the breach of contract was caused maliciously by us.

## Standard Terms & Conditions of Sale

- (6) The warranty period shall be one year from delivery of the goods. This shall not apply if the customer has not indicated the defect to us punctually (item 3 of this provision).
- (7) Only the product description shall be deemed agreed in respect of the nature of the goods, unless it has been expressly agreed between ourselves and the customer in writing that a certain nature of the goods is binding.
- (8) Our liability shall be limited to the legal warranty with the scope described above. We shall not provide the customer with warranties in the legal sense. Manufacturers' warranties in respect of the wearing parts used in our products shall be unaffected by this.
- (9) No warranty is given in respect of the deterioration of wearing parts which occurs as a result of the usual and proper use of our equipment.
- (10) Our equipment and products are intended solely for laboratory and research purposes. We shall not be liable for any damage arising from failure to observe the relevant accompanying assembly, operating, maintenance and inspection instructions. Neither shall we be liable should the customer carry out repairs, install other parts or make other technical changes to the equipment supplied by us without our consent. Neither shall we be liable for damage arising from the use of our equipment domestically or on humans or animals.
- (11) In the event of the sale of used equipment or used spare parts only the guarantees agreed in writing with authorised members of staff shall apply. There shall be no other warranty in respect of this equipment and these parts.
- (12) Assignment of compensation and warranty claims by the customer to third parties is impermissible.
- (13) We shall be liable for consequential damage of defects only in cases of grossly negligent or wilful breach of our contractual liabilities.
- (14) The advice given by us in respect of applied technology is given to the best of our knowledge. However, we shall be liable for any resulting damage only in cases of grossly negligent or deliberately incorrect advice.

### **§ 8. Assembly**

- (1) Assembly of our equipment shall always be carried out on the customer's site by our authorised technical staff.
- (2) In the event that assembly is to be carried out by the customer himself the customer shall ensure that the technical conditions specified by us (e.g. PC system requirements) are met and shall confirm this to us in writing on request. We accept no liability for damage arising from the fact that the necessary technical requirements are not provided by the customer. Any additional costs and expenses arising therefrom shall be for the customer's account.

### **§ 9. Miscellaneous, concluding provisions**

- (1) The law of the Federal Republic of Germany shall apply to all legal relations with our customers. The provisions of the UN Convention on the Sale of Goods shall not apply.
- (2) Special agreements shall require our written consent. This shall also apply to any departure from the present terms of business.
- (3) The place of performance for the customer's payment and other obligations and the sole legal venue for all disputes arising from this contract shall be our place of business in each case. This shall also apply if the customer has no general legal venue in Germany or his domicile or usual place of residence is not known at the time the claim is brought.
- (4) Should individual provisions of the contract with the customer – including these General Terms of Business – be or become ineffective in full or in part, the effectiveness of the other provisions shall be unaffected. The fully or partly ineffective ruling shall be replaced by a ruling which comes as close as possible to the financial result of the ineffective provision.

ORDERING - Fax: +49 (0)6172 1717799 - eMail: gmbh@licor.com  
ONLINE ORDERING at [www.licor.com/bio/store](http://www.licor.com/bio/store)

***New Empiria Studio Data Analysis Software  
Free 120 Day Trial at [licor.com/TryEmpiria](http://licor.com/TryEmpiria)***



*Expert-Level Analysis Made Simple. You'll Never Look Back!*

*On-Demand Western Blot Education Portal for Your Entire Lab  
Register now for free at*

**[www.lambdau.net](http://www.lambdau.net)**



**Lambda U**



Ordering  
Visit our online catalog  
Fax: +49 (0) 6172 - 1717799  
Email: [gmbh@licor.com](mailto:gmbh@licor.com)

***LI-COR***<sup>®</sup>