

LI-COR Biotechnology Pricelist



 United Kingdom

LI-COR[®]

Content

	Page
INSTRUMENTS & ACCESSORIES	
Odyssey® M	3
Odyssey® DLx	4
Odyssey® XF, D-DiGit™ Gel Scanner	5
C-DiGit® Blot Scanner & WesternSure® Reagents	6
Pearl® Imager Family- Small Animal <i>in vivo</i> Imaging Systems	7
Pearl Accessories	8
MPX™ (Multiplexer) Blotting System	9
COMPUTER SYSTEMS & IMAGER TRAINING	
Training for Odyssey Family Instruments & Pearl	9
ANALYSIS SOFTWARE & LICENSES	
Application and Data Analysis Software for Odyssey Family Instruments & Pearl	9
IRDye® CONJUGATES	
IRDye Secondary Antibodies	10
IRDye® AND VRDye™ LABELING REAGENTS	
IRDye NHS Esters and Maleimide	12
VRDye Secondary Antibodies	12
IRDye and VRDye Labeling Kits	13
Click Chemistry	13
NORMALISATION REAGENTS, PROTEIN MARKER, PRIMARY ANTIBODIES etc.	
Protein Marker	14
Streptavidin	14
IRDye Quencher	14
EMSA Reagents	14
Cell & Protein Stains	14
Normalisation Reagents, Primary Antibodies	15
BrightSite™ SMALL ANIMAL IMAGING AGENTS	
IRDye Cancer Probes	15
IRDye Structural & Control Probes	15
KITS, BUFFERS & MORE	
Western Kits, Stripping Buffer, Sample Buffer	16
Intercept Blocking Buffer	16
Intercept Antibody Diluents, Blot Storage	17
INCUBATION BOXES	17
WESTERN & IN-CELL WESTERN KITS	18
SERVICE CONTRACTS / EXTENDED WARRANTIES	19
CONTACT INFORMATION	20
STANDARD TERMS & CONDITIONS OF SALE	20
SHIPPING CHARGES	22

Odyssey Family Instruments & Accessories

Part No.	Description	£
3340-BA 3350-BA	Odyssey M (Fluorescence)- Data Integrity Bundle, Basic Odyssey M (Chemiluminescence/Fluorescence)- Data Integrity Bundle, Basic	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	Odyssey M (Fluorescence)- Data Integrity Bundle, Standard Odyssey M (Chemiluminescence/Fluorescence)- Data Integrity Bundle, Standard	please inquire
contains:	Odyssey M Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training Computer 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 Monitor 24" TFT Widescreen Monitor	
3340-PRE 3350-PRE	Odyssey M (Fluorescence)- Data Integrity Bundle, Premium Odyssey M (Chemiluminescence/Fluorescence)- Data Integrity Bundle, Premium	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 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 Monitor 24" TFT Widescreen Monitor	

Odyssey M Accessories

926-18971	Microscope Slide Holder	£1,483
926-18972	Microplate Insert	£1,483
926-70003	Silicone Mat (19 x 26 cm)	£95

Odyssey Family Instruments & Accessories

Part No.	Description	£
9142-BA	Odyssey DLx - Data Integrity Bundle, Basic	please inquire
contains:	Odyssey DLx Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training	
9142-STA	Odyssey DLx - Data Integrity Bundle, Standard	please inquire
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	
9142-PRE	Odyssey DLx - Data Integrity Bundle, Premium	please inquire
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)	£55
926-72200	6 MP Alignment Guide (for multiple plates)	£63
926-70000	Silicone Mat (13 x 13 cm)	£54
926-70001	Silicone Mat (22 x 22cm)	£127
926-71000	4" Soft Roller	£66
926-71804	Odyssey Pen 4-Pack, black fine-liner for membranes	£44
929-97901	Mouse Nose Cone	£98
929-97902	Rat Nose Cone	£98

Odyssey Family Instruments & Accessories, D-DiGit

Part No.	Description	£
2802-BA	Odyssey XF - Data Integrity Bundle, Basic	please inquire
contains:	Odyssey XF Infrared Imaging System and Accessory Pack Empiria Studio Analysis Software (10 licenses), LI-COR Acquisition Software Installation & Training	
2802-STA	Odyssey XF - Data Integrity Bundle, Standard	please inquire
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	
2802-PRE	Odyssey XF - Data Integrity Bundle, Premium	please inquire
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	
Odyssey XF/Fc Accessories		
926-40020	Odyssey XF/Fc Imaging Tray, 5 pack	£74
9957-046	Desiccant Replacement Kit	£373
3500-00	D-DiGit Gel Scanner	£4,070
3500-80	D-DiGit One Year Extended Warranty	£289
3500-81	D-DiGit Two Years Extended Warranty <i>Only available at the initial D-DiGit purchase</i>	£519
3600-10	DiGit Duo	£10,310
	contains: C-DiGit Blot Scanner, D-DiGit Gel Scanner, Laptop PC	



C-DiGit Blot Scanner

Part No.	Description	£
3600-00	C-DiGit Blot Scanner	£6,808
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	£1,819
3600-500U	C-DiGit Image Studio Analysis Software - Upgrade Package	£978
3600-501	C-DiGit Image Studio Analysis Software - additional 1 user license	£143
3600-505	C-DiGit Image Studio Analysis Software - additional 5 user license	£389
3600-510	C-DiGit Image Studio Analysis Software - additional 25 user license	£719
3600-80	C-DiGit One Year Extended Warranty	£450
3600-81	C-DiGit Two Years Extended Warranty <i>Only available at the initial C-DiGit purchase</i>	£814
616-06141LCDG	Laptop PC for the C-DiGit	£1,467

WesternSure Chemiluminescent Reagents

Part No.	Description	
HRP Conjugates		
926-80010	WesternSure Goat anti-Mouse HRP, 1.0 ml	£139
926-80011	WesternSure Goat anti-Rabbit HRP, 1.0 ml	£139
Chemiluminescence Substrates		
926-95000	WesternSure PREMIUM Chemiluminescent Substrate, 100 ml kit	£422
926-95010	WesternSure PREMIUM Chemiluminescent Substrate, 20 ml kit	£121
Protein Marker		
926-91000	WesternSure Pen	£105
926-98000	WesternSure Pre-Stained Chemiluminescent Protein Ladder	£231

Pearl Instruments

Part No.	Description	£
<p>9430-01 contains:</p>	<p>Pearl Trilogy Standard Pearl Trilogy Small Animal Imaging System 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</p>	<p>please inquire</p>
<p>9430-03 contains:</p>	<p>Pearl Trilogy Premium Pearl Trilogy Imager Standard (9430-01) and Image Studio Software 25 User License*</p>	<p>please inquire</p>
<p>9430-05 contains:</p>	<p>Pearl Trilogy Premium Suite 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</p>	<p>please inquire</p>
<p>9000-100</p>	<p>SmartFlow Anesthesia Suite</p>	<p>£13,767</p>

Pearl Accessories

Part No.	Description	£
9300-20	Pearl Docking Station	£12,720
9300-21	Pearl Imaging Bed	£5,940
9300-22	Pearl Clean Box	£12,720
9000-140	Pearl Imaging Cart	£11,472
9000-822	SmartFlow Rebreather, large	£5,641
9000-802	Mouse Nose Cone, large (3 pieces)	£329
9000-803	Mouse Nose Cone, small (3 pieces)	£329
9000-804	Nose Cone Plugs, large (6 pieces)	£218
9300-800	Nose Cone Plugs, small (6 pieces)	£237
300-08568	Rat Nose Cone	£93
9300-801	Rubber Cord for Imaging Bed (10')	£192
9300-810	Imaging Blankets (25 Blankets)	£237
9300-820	Charcoal Filter (8 Filters)	£305
9300-821	HEPA-Filter Assembly	£377
9300-830	Clean Box Cleaning Kit	£377
	contains: # 9300-831 (Heavy Duty Optical Wipes), # 9300-832 (Optical Tissue), # 9300-833 (Optical Cleaner).	
9300-831	Heavy Duty Optical Wipes (300 wipes)	£190
	Durable Technicloth wipers for use with Optical Cleaner for heavy duty cleaning of the window or entire Clean Box.	
9300-832	Optical Tissue (150 tissues)	£190
	Premium Grade Optical Tissue for use with Optical Cleaner for cleaning of the window or entire Clean Box.	
9300-833	Optical Cleaner (3 Pack)	£190
	Formula MC Glass Cleaner for use with Optical Tissue for cleaning of the window or entire Clean Box.	
9300-834	Clean Box Cover Slip	£484
9300-835	Glass Replacement Kit	£1,407
9957-046	Desiccant Replacement Kit	£373
9300-840	Dye Decontamination Kit	£198
	contains: 200 Povidone-Iodine Pads, 400 Alcohol Prep Pads	
9300-841	Povidone Iodine Refills, 400 Pads	£132
9300-842	Alcohol Preparation Pad Refills, 400 Pads	£132
9300-23	Organ Tray Base	£2,841
9300-836	Disposable Organ Trays (5 pcs)	£154
9300-837	Disposable Organ Trays (25 pcs)	£586
9300-850	Compound Injection Clips (25 pcs)	£167
9300-851	Compound Injection Clips (100 pcs)	£478

MPX Blotting System

Part No.	Description	£
MPX (Multiplexer) Blotting System		
Helps to use your antibodies at optimised conditions. No cutting, stripping and re-probing.		
921-00000	MPX Blotting System	£610
921-00110	MPX Clamp Nut (2 Ea.)	£94
921-00120	MPX Membrane Cushion (Pack of 10)	£65
921-00200	MPX Comb, One-Lane plus One Marker Lane	£99
921-00202	MPX Comb, Two-Lane plus One Marker Lane	£99
921-00204	MPX Comb, Three-Lane plus One Marker Lane	£99
921-00206	MPX Comb, Four-Lane plus One Marker Lane	£99

Computer and Monitors

616-06141	Computer	please inquire
The PC has the following min. configuration: 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)	£2,060
9300-50	Pearl Imager Training/Retraining (per day)	£2,060

Empiria Studio - Data Analysis Software & Licenses

Part No.	Description	£
Empiria Studio for Odyssey CLx and Fc		
9141-500	Empiria Studio Analysis Software - 1 User License	£272
9141-505	Empiria Studio Analysis Software - 5 User Licenses	£1,023
9141-505IS	Empiria Studio Analysis Software - 5 User Licenses and Image Studio Analysis Software - 1 user license	£1,250
9141-507	Empiria Studio Analysis Software - 10 user licenses	£1,705
9141-510	Empiria Studio Analysis Software - 25 User Licenses	£3,410
9141-510IS	Empiria Studio Analysis Software - 25 User Licenses Image Studio Analysis Software - 1 user license	£3,500
9141-500U	Empiria Studio Analysis Software - Upgrade Package	£90




Free 120 Day Trial: licor.com/TryEmpiria

IRDye conjugated Secondary Antibodies

Part No.	Description	£
Secondary Antibodies for the 800 nm Channel		
925-32210	IRDye 800CW Goat Anti-Mouse IgG (H+L), 0.1 mg	£116
925-32211	IRDye 800CW Goat Anti-Rabbit IgG (H+L), 0.1 mg	£116
925-32212	IRDye 800CW Donkey Anti-Mouse IgG (H+L), 0.1 mg	£116
925-32213	IRDye 800CW Donkey Anti-Rabbit IgG (H+L), 0.1 mg	£116
925-32214	IRDye 800CW Donkey Anti-Goat IgG (H+L), 0.1 mg	£116
925-32218	IRDye 800CW Donkey Anti-Chicken IgG (H+L), 0.1 mg	£116
925-32219	IRDye 800CW Goat Anti-Rat IgG (H+L), 0.1 mg	£116
925-32232	IRDye 800CW Goat Anti-Human IgG (H+L), 0.1 mg	£116
925-32411	IRDye 800CW Donkey Anti-Guinea Pig IgG (H+L), 0.1 mg	£116
925-32280	IRDye 800CW Goat Anti-Mouse IgM (μ Chain Specific), 0.1 mg	£116
Secondary Antibodies for the 700 nm Channel		
925-68070	IRDye 680RD Goat Anti-Mouse IgG (H+L), 0.1 mg	£116
925-68071	IRDye 680RD Goat Anti-Rabbit IgG (H+L), 0.1 mg	£116
925-68072	IRDye 680RD Donkey Anti-Mouse IgG (H+L), 0.1 mg	£116
925-68073	IRDye 680RD Donkey Anti-Rabbit IgG (H+L), 0.1 mg	£116
925-68074	IRDye 680RD Donkey Anti-Goat IgG (H+L), 0.1 mg	£116
925-68075	IRDye 680RD Donkey Anti-Chicken IgG (H+L), 0.1 mg	£116
925-68076	IRDye 680RD Goat Anti-Rat IgG (H+L), 0.1 mg	£116
925-68077	IRDye 680RD Donkey Anti-Guinea Pig IgG (H+L), 0.1 mg	£116
925-68078	IRDye 680RD Goat Anti-Human IgG (H+L), 0.1 mg	£116
925-68180	IRDye 680RD Goat Anti-Mouse IgM (μ Chain Specific), 0.1 mg	£116
925-68020	IRDye 680LT Goat Anti-Mouse IgG (H+L), 0.1 mg	£116
925-68021	IRDye 680LT Goat Anti-Rabbit IgG (H+L), 0.1 mg	£116
925-68022	IRDye 680LT Donkey Anti-Mouse IgG (H+L), 0.1 mg	£116
925-68023	IRDye 680LT Donkey Anti-Rabbit IgG (H+L), 0.1 mg	£116
925-68024	IRDye 680LT Donkey Anti-Goat IgG (H+L), 0.1 mg	£116
925-68028	IRDye 680LT Donkey Anti-Chicken IgG (H+L), 0.1 mg	£116
925-68029	IRDye 680LT Goat Anti-Rat IgG (H+L), 0.1 mg	£116
925-68030	IRDye 680LT Donkey Anti-Guinea Pig IgG (H+L), 0.1 mg	£116
925-68032	IRDye 680LT Goat Anti-Human IgG (H+L), 0.1 mg	£116
925-68080	IRDye 680LT Goat Anti-Mouse IgM (μ Chain Specific), 0.1 mg	£116

IRDye conjugated Secondary Antibodies

Part No.	Description		£
Secondary Antibodies for the 800 nm Channel			
			
926-32210	IRDye 800CW Goat Anti-Mouse IgG (H+L), 0.5 mg		£195
926-32211	IRDye 800CW Goat Anti-Rabbit IgG (H+L), 0.5 mg		£195
926-32212	IRDye 800CW Donkey Anti-Mouse IgG (H+L), 0.5 mg		£195
926-32213	IRDye 800CW Donkey Anti-Rabbit IgG (H+L), 0.5 mg		£195
926-32214	IRDye 800CW Donkey Anti-Goat IgG (H+L), 0.5 mg		£195
926-32218	IRDye 800CW Donkey Anti-Chicken IgG (H+L), 0.5 mg		£195
926-32219	IRDye 800CW Goat Anti-Rat IgG (H+L), 0.5 mg		£195
926-32232	IRDye 800CW Goat Anti-Human IgG (H+L), 0.5 mg		£195
926-32411	IRDye 800CW Donkey Anti-Guinea Pig IgG (H+L), 0.5 mg		£195
926-32350	IRDye 800CW Goat anti-Mouse IgG ₁ Specific, 0.5 mg		£300
926-32351	IRDye 800CW Goat anti-Mouse IgG _{2a} Specific, 0.5 mg		£300
926-32352	IRDye 800CW Goat anti-Mouse IgG _{2b} Specific, 0.5 mg		£300
926-32280	IRDye 800CW Goat Anti-Mouse IgM (μ Chain Specific), 0.5 mg		£195
926-32202	IRDye 800CW Rabbit Anti-HRP IgG (H+L), 0.5 mg		£195
Secondary Antibodies for the 700 nm Channel			
<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		£195
926-68071	IRDye 680RD Goat Anti-Rabbit IgG (H+L), 0.5 mg		£195
926-68072	IRDye 680RD Donkey Anti-Mouse IgG (H+L), 0.5 mg		£195
926-68073	IRDye 680RD Donkey Anti-Rabbit IgG (H+L), 0.5 mg		£195
926-68074	IRDye 680RD Donkey Anti-Goat IgG (H+L), 0.5 mg		£195
926-68075	IRDye 680RD Donkey Anti-Chicken IgG (H+L), 0.5 mg		£195
926-68076	IRDye 680RD Goat Anti-Rat IgG (H+L), 0.5 mg		£195
926-68077	IRDye 680RD Donkey Anti-Guinea Pig IgG (H+L), 0.5 mg		£195
926-68078	IRDye 680RD Goat Anti-Human IgG (H+L), 0.5 mg		£195
926-68180	IRDye 680RD Goat Anti-Mouse IgM (μ Chain Specific), 0.5 mg		£195
<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		£195
926-68021	IRDye 680LT Goat Anti-Rabbit IgG (H+L), 0.5 mg		£195
926-68022	IRDye 680LT Donkey Anti-Mouse IgG (H+L), 0.5 mg		£195
926-68023	IRDye 680LT Donkey Anti-Rabbit IgG (H+L), 0.5 mg		£195
926-68024	IRDye 680LT Donkey Anti-Goat IgG (H+L), 0.5 mg		£195
926-68028	IRDye 680LT Donkey Anti-Chicken IgG (H+L), 0.5 mg		£195
926-68029	IRDye 680LT Goat Anti-Rat IgG (H+L), 0.5 mg		£195
926-68032	IRDye 680LT Goat Anti-Human IgG (H+L), 0.5 mg		£195
926-68030	IRDye 680LT Donkey Anti-Guinea Pig IgG (H+L), 0.5 mg		£195
926-68050	IRDye 680LT Goat anti-Mouse IgG ₁ Specific, 0.5 mg		£300
926-68051	IRDye 680LT Goat anti-Mouse IgG _{2a} Specific, 0.5 mg		£300
926-68052	IRDye 680LT Goat anti-Mouse IgG _{2b} Specific, 0.5 mg		£300
926-68080	IRDye 680LT Goat Anti-Mouse IgM (μ Chain Specific), 0.5 mg		£195

IRDye Labeling Reagents & VRDye™ Conjugates

Part No.	Description	£
<i>Infrared Dyes for the 800 nm Channel</i>		
929-70020	IRDye 800CW Infrared Dye NHS ester (0.5 mg)	£152
929-70021	IRDye 800CW Infrared Dye NHS ester (5.0 mg)	£860
929-70022	IRDye 800CW Infrared Dye NHS ester (50.0 mg)	£4,860
929-70023	IRDye 800CW Infrared Dye NHS ester (100.0 mg)	£6,734
929-70024	IRDye 800CW Infrared Dye NHS ester (250.0 mg)	£8,742
929-70025	IRDye 800CW Infrared Dye NHS ester (500.0 mg)	£11,361
929-70026	IRDye 800CW Infrared Dye NHS ester (1.0 g)	£14,785
929-80020	IRDye 800CW Maleimide (0.5 mg)	£152
929-80021	IRDye 800CW Maleimide (5.0 mg)	£860
929-80050	IRDye 800CW Maleimide (50.0 mg)	£4,860
929-72020	IRDye 800RS Infrared Dye NHS ester (0.5 mg)	£152
929-72021	IRDye 800RS Infrared Dye NHS ester (5.0 mg)	£860
<i>Infrared Dyes for the 700 nm Channel</i>		
929-70050	IRDye 680RD Infrared Dye NHS ester (0.5 mg)	£152
929-70051	IRDye 680RD Infrared Dye NHS ester (5.0 mg)	£860
929-70052	IRDye 680RD Infrared Dye NHS ester (50.0 mg)	£5,705
929-71050	IRDye 680RD Maleimide (0.5 mg)	£152
929-71051	IRDye 680RD Maleimide (5.0 mg)	£860
929-71010	IRDye 680LT Infrared Dye NHS ester (0.5 mg)	£152
929-71011	IRDye 680LT Infrared Dye NHS ester (5.0 mg)	£1,059
929-71500	IRDye 680LT Infrared Dye NHS ester (50.0 mg)	£5,705
929-71008	IRDye 680LT Maleimide (0.5 mg)	£152
929-71009	IRDye 680LT Maleimide (5.0 mg)	£860
<i>Fluorescently labeled secondary antibodies for Odyssey M</i>		
926-49010	VRDye 490 Goat anti-Mouse IgG, (H + L) 0.5 mg	£219
926-49020	VRDye 490 Goat anti-Rabbit IgG, (H + L) 0.5 mg	£219
926-54010	VRDye 549 Goat anti-Mouse IgG, (H + L) 0.5 mg	£219
926-54020	VRDye 549 Goat anti-Rabbit IgG, (H + L) 0.5 mg	£219
926-54012	VRDye™ 549 Donkey anti-Mouse IgG Secondary Antibody	£219
926-54013	VRDye™ 549 Donkey anti-Rabbit IgG Secondary Antibody	£219
926-54014	VRDye™ 549 Donkey anti-Goat IgG Secondary Antibody	£219
926-49012	VRDye™ 490 Donkey anti-Mouse IgG Secondary Antibody	£219
926-49013	VRDye™ 490 Donkey anti-Rabbit IgG Secondary Antibody	£219
926-49014	VRDye™ 490 Donkey anti-Goat IgG Secondary Antibody	£219
926-54030	VRDye™ 549 Streptavidin	£184
926-49030	VRDye™ 490 Streptavidin	£184

IRDye & VRDye Labelling Kits

Part No.	Description	£
<i>Infrared Labelling Kits for the 800 nm Channel</i>		
928-38040	IRDye 800CW High MW Protein Labeling Kit	£527
928-38042	IRDye 800CW Low MW Protein Labeling Kit	£579
928-38044	IRDye 800CW Microscale Protein Labeling Kit	£432
<i>Infrared Labelling Kits for the 700 nm Channel</i>		
928-38072	IRDye 680RD High MW Protein Labeling Kit	£527
928-38074	IRDye 680RD Low MW Protein Labeling Kit	£579
928-38076	IRDye 680RD Microscale Protein Labeling Kit	£432
928-38066	IRDye 680LT High MW Protein Labeling Kit	£527
928-38068	IRDye 680LT Low MW Protein Labeling Kit	£579
928-38070	IRDye 680LT Microscale Protein Labeling Kit	£432
<i>Fluorescence Labelling Kits for other detection systems</i>		
928-49040	VRDye 490 High MW Protein Labeling Kit	£527
928-49044	VRDye 490 Microscale Protein Labeling Kit	£432
928-54040	VRDye 549 High MW Protein Labeling Kit	£527
928-54044	VRDye 549 Microscale Protein Labeling Kit	£432
<i>Click Chemistry</i>		
929-50000	IRDye 800CW DBCO, 0.5 mg	£450
929-55000	IRDye 800CW DBCO, 5.0 mg	£860
929-55050	IRDye 800CW DBCO, 50.0 mg	£4,860
929-50005	IRDye 680RD DBCO, 0.5 mg	£450
929-55005	IRDye 680RD DBCO, 5.0 mg	£860
929-60000	IRDye 800CW Azide, 0.5 mg	£450
929-65000	IRDye 800CW Azide, 5.0 mg	£860
929-75050	IRDye 800CW Azide, 50.0 mg	£4,860
929-60002	IRDye 800CW Alkyne, 0.5 mg	£450
929-65002	IRDye 800CW Alkyne, 5.0 mg	£860
929-65050	IRDye 800CW Alkyne, 50.0 mg	£4,860
929-60005	IRDye 680RD Azide, 0.5 mg	£450
929-65005	IRDye 680RD Azide, 5.0 mg	£860
929-60006	IRDye 680RD Alkyne, 0.5 mg	£450
929-65006	IRDye 680RD Alkyne, 5.0 mg	£860

Protein Marker, Stains

Part No.	Description	£
Protein Marker		
928-40000	Protein Molecular Weight Marker, 500µl (10 - 250 kDa)	£219
928-50000	Chameleon™ Vue Pre-stained Protein Ladder, 250µl	£145
928-60000	Chameleon Duo Pre-stained Protein Ladder, 500µl	£218
928-90000	Chameleon Kit Pre-stained Protein Ladder, 2x 250µl	£218
928-70000	Chameleon 700 Pre-stained Protein Ladder, 500µl	£171
928-80000	Chameleon 800 Pre-stained Protein Ladder, 500µl	£171
Streptavidin		
925-32230	IRDye 800CW Streptavidin (0.1 mg)	£116
926-32230	IRDye 800CW Streptavidin (0.5 mg)	£184
925-68079	IRDye 680RD Streptavidin (0.1 mg)	£116
926-68079	IRDye 680RD Streptavidin (0.5 mg)	£184
925-68031	IRDye 680LT Streptavidin (0.1 mg)	£116
926-68031	IRDye 680LT Streptavidin (0.5 mg)	£184
IRDye Quencher		
929-70030	IRDye QC-1 NHS Ester (0.5 mg)	£285
929-70031	IRDye QC-1 NHS Ester (5.0 mg)	£1,251
Cell & Protein Stains		
929-90010	CellVue Burgundy (100 nmol, 10 ml of Diluent C)	£356
929-90020	CellVue NIR815 (100 nmol, 10 ml of Diluent C)	£356
929-90001	CellVue Diluent C (6x 10 ml)	£328
929-90030	PSVue 794 Reagent Kit	£696
EMSA Reagents		
829-07921	p53 IRDye 700 (1.25 nmol)	£78
829-07922	STAT3 IRDye 700 (1.25 nmol)	£78
829-07923	CREB IRDye 700 (1.25 nmol)	£78
829-07924	NFkB IRDye 700 (1.25 nmol)	£78
829-07925	AP-1 IRDye 700 (1.25 nmol)	£78
829-07926	Sp-1 IRDye 700 (1.25 nmol)	£78
829-07929	HIF-1 IRDye 700 (1.25 nmol)	£78
829-07910	Odyssey EMSA Buffer Kit (100 EMSAs) 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	£288
927-10100	10x Orange Loading Dye (3 tubes à 1.0 ml)	£82

Normalisation Reagents, Primary Antibodies

Part No.	Description	£
Total Protein Stains for Normalisation		
926-11010	Revert 700 Total Protein Stain Kit 100ml, incl. Stain, Wash and Reversal Solution*	£141
926-11016	Revert 700 Total Protein Stain Kit 250ml, incl. Stain, Wash and Reversal Solution*	£264
926-11015	Revert 700 Stain and Wash Solution Kit 100 ml, incl. Stain and Wash Solution	£121
926-11011	Revert 700 Total Protein Stain, 100ml	£105
926-11021	Revert 700 Total Protein Stain, 250ml	£192
926-10010	Revert 520 Total Protein Stain Kit 200ml, incl. Stain, Wash and Reversal Solution*	£162
926-10016	Revert 520 Total Protein Stain Kit 500ml, incl. Stain, Wash and Reversal Solution*	£303
926-10015	Revert 520 Stain and Wash Solution Kit 200 ml, incl. Stain and Wash Solution	£140
926-10011	Revert 520 Total Protein Stain, 200ml	£120
926-10021	Revert 520 Total Protein Stain, 500ml	£220
Loading Indicators		
926-20002	Odyssey 28kD Loading Indicator - 800nm	£131
926-20004	Odyssey 28kD Loading Indicator - 700nm	£131
Primary Antibodies		
926-42210	β -Actin Rabbit monoclonal antibody, 0.1 ml	£414
926-42211	β -Tubulin Rabbit polyclonal antibody, 0.1 ml	£414
926-42212	β -Actin Mouse monoclonal antibody, 0.1 ml	£414
926-42213	α -Tubulin Mouse monoclonal antibody, 0.1 ml	£414
926-42214	COX IV Rabbit monoclonal antibody, 0.1 ml	£414
926-42215	Vinculin Rabbit Monoclonal Antibody, 0.1 ml	£414
926-42216	GAPDH Rabbit Monoclonal Antibody, 0.1 ml	£414
926-42218	Histone H3 Mouse Monoclonal Antibody, 0.1 ml	£414
926-42219	Histone H3 Rabbit Monoclonal Antibody, 0.1 ml	£414

BrightSite Small Animal Imaging Agents

Part No.	Description	£
Probes		
926-08446	IRDye 800CW EGF Optical Probe (20 nmol, lyophilized)	£480
926-08946	IRDye 800CW 2-DG Optical Probe (100 nmol, lyophilized)	£480
926-09889	IRDye 800CW RGD Optical Probe (15 nmol, lyophilized)	£480
926-50401	IRDye 800CW PEG Contrast Agent (15 nmol, lyophilized)	£480
926-09375	IRDye 800CW BoneTag Optical Probe (40 nmol, lyophilized)	£480
926-09374	IRDye 680RD BoneTag Optical Probe (40 nmol, lyophilized)	£480
Control Probes		
929-08972	IRDye 800CW Carboxylate (100 nmol, lyophilized)	£185
929-09406	IRDye 800CW Carboxylate (5 mg)	£754
929-70122	IRDye 800CW Carboxylate (50 mg)	£2,277
929-70123	IRDye 800CW Carboxylate (100 mg)	£2,959
929-70124	IRDye 800CW Carboxylate (250 mg)	£4,763
929-70125	IRDye 800CW Carboxylate (500 mg)	£6,216
929-70126	IRDye 800CW Carboxylate (1 g)	£8,083
929-08980	IRDye 680RD Carboxylate (100 nmol, lyophilized)	£185

Kits, Buffers & more

Part No.	Description	£
<i>Protein Sample Loading Buffer</i>		
928-40004	4x Protein Sample Loading Buffer (15 ml)	£45
<i>Western Blotting Membranes & Kits</i>		
926-31090	Odyssey Nitrocellulose Membranes (10 sheets, 0.22 µm, 7 x 8.5 cm)	£79
926-31092	Roll of Nitrocellulose Membranes (1 roll, 0.22 µm, 30 cm x 3 m) *	£256
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)	£509
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)	£1,349
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)	£1,349
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)	£591
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)	£591
<i>Stripping Buffer</i>		
926-92000	WesternSure ECL Stripping Buffer (100 ml, 5x concentrated)	£145
928-40028	NewBlot IR Stripping Buffer (100 ml, 5x concentrated)	£145
928-40030	New Blot™ Nitro Stripping Buffer (100 ml, 5x concentrated) *	£223
928-40032	New Blot™ PVDF Stripping Buffer (100 ml, 5x concentrated) *	£225
<i>Intercept Blocking Buffer</i>		
927-60001	Intercept (TBS) Blocking Buffer, 500 ml	£124
927-60003	Intercept (TBS) Blocking Buffer, 3x 500 ml	£332
927-60010	Intercept (TBS) Blocking Buffer, 10x 500 ml	£976
927-60025	Intercept (TBS) Blocking Buffer, 25x 500 ml	£2,138
927-60050	Intercept (TBS) Blocking Buffer, 50x 500 ml	£3,735
927-70001	Intercept (PBS) Blocking Buffer, 500 ml	£124
927-70003	Intercept (PBS) Blocking Buffer, 3x 500 ml	£332
927-70010	Intercept (PBS) Blocking Buffer, 10x 500 ml	£976
927-70025	Intercept (PBS) Blocking Buffer, 25x 500 ml	£2,138
927-70050	Intercept (PBS) Blocking Buffer, 50x 500 ml	£3,735
927-80001	Intercept (TBS) Protein-Free Blocking Buffer, 500 ml	£124
927-80003	Intercept (TBS) Protein-Free Blocking Buffer, 3x 500 ml	£332
927-80010	Intercept (TBS) Protein-Free Blocking Buffer, 10x 500 ml	£976
927-90001	Intercept (PBS) Protein-Free Blocking Buffer, 500 ml	£124
927-90003	Intercept (PBS) Protein-Free Blocking Buffer, 3x 500 ml	£332
927-90010	Intercept (PBS) Protein-Free Blocking Buffer, 10x 500 ml	£976

Antibody Diluents & Incubation Boxes

Part No.	Description	£
Intercept Antibody Diluents		
927-65001	Intercept T20 (TBS) Antibody Diluent, 500ml	£131
927-65003	Intercept T20 (TBS) Antibody Diluent, 3x 500ml	£350
927-65010	Intercept T20 (TBS) Antibody Diluent, 10x 500ml	£1,028
927-75001	Intercept T20 (PBS) Antibody Diluent, 500ml	£131
927-75003	Intercept T20 (PBS) Antibody Diluent, 3x 500ml	£350
927-75010	Intercept T20 (PBS) Antibody Diluent, 10x 500ml	£1,028
927-66003	Intercept T20 (TBS) Blocking Buffer & Antibody Diluent Kit (contains: Intercept (TBS) Blocking Buffer, 500ml and Intercept (TBS) Antibody Diluent, 2x 500ml)	£332
927-76003	Intercept T20 (PBS) Blocking Buffer & Antibody Diluent Kit (contains: Intercept (PBS) Blocking Buffer, 500ml and Intercept (PBS) Antibody Diluent, 2x 500ml)	£332
927-85001	Intercept T20 (TBS) Protein-Free Antibody Diluent, 500ml	£131
927-85003	Intercept T20 (TBS) Protein-Free Antibody Diluent, 3x 500ml	£350
927-85010	Intercept T20 (TBS) Protein-Free Antibody Diluent, 10x 500ml	£1,028
927-95001	Intercept T20 (PBS) Protein-Free Antibody Diluent, 500ml	£131
927-95003	Intercept T20 (PBS) Protein-Free Antibody Diluent, 3x 500ml	£350
927-95010	Intercept T20 (PBS) Protein-Free Antibody Diluent, 10x 500ml	£1,028
Blot Storage		
929-95100	Western Blot Storage Bags, 40 pc	£35
Black Western Incubation Boxes		
Small - Size: 7.2 x 4.9 x 3.0 cm		
929-97101	Western Incubation Box Small One Pack	£17
929-97105	Western Incubation Box Small Five Pack	£50
929-97110	Western Incubation Box Small Ten Pack	£79
Medium - Size: 8.9 x 6.5 x 2.8 cm		
929-97201	Western Incubation Box Medium One Pack	£20
929-97205	Western Incubation Box Medium Five Pack	£56
929-97210	Western Incubation Box Medium Ten Pack	£91
Large - Size: 11.5 x 8.8 x 2.8 cm		
929-97301	Western Incubation Box Large One Pack	£24
929-97305	Western Incubation Box Large Five Pack	£70
929-97310	Western Incubation Box Large Ten Pack	£121
Extra Large - Size: 15.2 x 10.1 x 3.1 cm		
929-97401	Western Incubation Box Extra Large One Pack	£25
929-97405	Western Incubation Box Extra Large Five Pack	£83
929-97410	Western Incubation Box Extra Large Ten Pack	£140

Western & In-Cell Western Kits

Part No.	Description	£
Western Blotting Kits		
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)	£415
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)	£415
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)	£415
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)	£415
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)	£415
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)	£415
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)	£415
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)	£415
CellTag™ In-Cell Western Kits		
926-42091	GAM 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	£374
926-42092	GAR 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	£374
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	£448
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	£448
926-41090	CellTag™ 700 Stain, 2 x 10 nmol	£192
926-41094	CellTag™ 520 Stain, 2 x 50 nmol	£270
928-40022	Sapphire700 Stain (100 µl)	£130
Quick Western Kits		
926-69100	Quick Western Kit - IRDye 680RD IRDye 680RD Detection Reagent, Intercept Blocking Buffer	£435
Microplates		
926-19156	96-well ICW Microplate, 8-pack (back with clear bottom)	£115
926-19157	96-well ICW Microplate, 32-pack (back with clear bottom)	£343

Service Contracts / Extended Warranties

Part No.	Description	£
Basic Extended Warranties		
3340-80E	Odyssey M Basic Extended Warranty, 1 year	£5,885
9142-80E	Odyssey DLx Basic Extended Warranty, 1 year	£3,959
2802-80E	Odyssey XF Basic Extended Warranty, 1 year	£2,889
9400-80E	Pearl Basic Extended Warranty, 1 year	£3,959

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

Standard Extended Warranties		
3340-82E	Odyssey M Standard Extended Warranty, 1 year	£6,206
9142-82E	Odyssey DLx Standard Extended Warranty, 1 year	£4,280
2802-82E	Odyssey XF Standard Extended Warranty, 1 year	£3,210
9400-82E	Pearl Standard Extended Warranty, 1 year	£4,280

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)

Premium Support Contracts		
3340-84	Odyssey M Premium Support Contract, 1 year	£7,597
9142-84	Odyssey DLx Premium Support Contract, 1 year	£5,671
2802-84	Odyssey XF Premium Support Contract, 1 year	£4,601
9400-84	Pearl Premium Support Contract, 1 year	£5,671

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

Contact Information

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road, Cambridge CB4 0WS
Phone: +44 (0) 1223-422104
Email: bio-eu@licor.com

www.licor.com/bio

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. Definitions

For the purposes of the Contract the terms listed below shall bear the following meanings:-

"Buyer" the person, firm or company purchasing the Products under the Contract.

"Contract" the contract or contracts for the purchase of the Products by the Buyer comprising these Terms and the Order.

"LI-COR" LI-COR Biosciences UK Limited, its assignees and successors in title.

"Order" the Buyer's order for the Products as accepted by LI-COR and attached to these Terms.

"Products" the Products to be purchased by the Buyer under the Contract according to the Specifications.

"Specifications" the requirements as to category, quality and quantity of the Products (if any) detailed in the Order.

"Terms" these terms and conditions of sale.

§ 2. General provisions, scope

(1) The Products are sold upon these Terms and no agent or representative of LI-COR has any authority to vary or omit these Terms or any part thereof. Any terms and conditions printed on the Buyer's order forms or in the correspondence of either party or elsewhere or implied (whether in writing or verbally) by trade, custom or practice or course of dealing, are binding only insofar as they do not vary from these Terms and have been specifically agreed in writing by an authorised representative of LI-COR and any purported provisions to the contrary are hereby expressly excluded.

(2) These Terms together with any particulars incorporated in the Order constitute the entire agreement between the parties concerning the supply of the Products and supersede any prior promises, undertakings, contracts, agreements or implications (whether written or oral). The Buyer hereby acknowledges that it has not entered into the Contract in reliance upon any representation other than those which have been reduced to writing and expressly included in the Contract.

(3) LI-COR does not effect legal transactions with private consumers. Consequently, these Terms have no validity in respect of and are not intended for transactions with private end-users.

§ 3. Formation of the Contract

(1) Any quotations issued by LI-COR are always non-binding and do not constitute offers capable of acceptance, and LI-COR reserves the right to withdraw or revise such quotations at any time before accepting the Order.

(2) By submitting the Order the Buyer makes a binding declaration that it wishes to acquire the Products on these Terms. LI-COR shall be entitled to accept the contract offer inherent in the Order within two weeks of receipt by LI-COR by issuing a written Order confirmation. This Contract shall not exist until LI-COR has issued such written Order confirmation or the Order has been executed in full.

(3) If the Buyer submits the Order electronically LI-COR will confirm receipt of the Order without delay unless precluded from doing so by malfunctions on its website. The confirmation of receipt shall not form an acceptance of the Order which shall follow thereafter.

(4) LI-COR accepts the Order subject to being supplied correctly and punctually by its own suppliers. Accordingly, LI-COR shall be in no manner liable for non-delivery caused by the non-fulfilment of LI-COR's orders by LI-COR's suppliers.

(5) The Buyer will be informed without delay if it is not possible to execute the Order. Any payment will be reimbursed as soon as reasonably possible.

(6) If the Buyer submits the Order electronically the text of the Contract will be stored by LI-COR and sent to the Buyer electronically (by e-mail) on request along with a copy of these Terms.

§ 4. Delivery and payment

(1) LI-COR will endeavour to comply with any delivery dates and deadlines stated in the Order confirmation provided that all necessary information has been supplied in a timely fashion by the Buyer with the Order. However, time shall not be of the essence in respect of the performance of LI-COR's obligations under the Contract, and details of delivery and performance dates are provided for guidance only and are non-binding on LI-COR.

(2) The Buyer's sole remedy in the event of non-delivery which is not the fault of the Buyer shall be the reimbursement of sums paid for the Order.

(3) Delivery shall be at the premises of the Buyer or at such other premises as are designated by the Buyer in writing and shall be at the cost and risk of the Buyer.

(4) The costs of carriage and packaging will be charged to the Buyer on the invoice for the Products at the time of despatch. All packaging is non-returnable unless otherwise stated in writing.

(5) Each delivery and the Order shall constitute a separate Contract under these Terms and LI-COR's failure to make any delivery in full, on time or at all shall not vitiate the Contract as a whole.

(6) If performance or delivery is delayed or cancelled at the Buyer's request or for reasons for which the Buyer is responsible LI-COR shall be entitled to claim any expenses which it has already incurred to date fulfilling the Order and the Buyer shall promptly reimburse LI-COR for such costs upon receiving an invoice for the same.

Standard Terms & Conditions of Sale

(7) The scope of supply shall be determined by LI-COR's written Order confirmation.

(8) LI-COR reserves the right to make technical changes to the Products – in particular construction and design changes – based on improvements in technology or changed legislative requirements within the confines of what is reasonable, even after formation of the Contract, provided that the fitness of the Products for their intended purpose is not thereby materially impaired. The same shall apply to changes in the scope of supply.

(9) If partial deliveries are necessary, the higher carriage costs thereby incurred will be invoiced in addition and paid for by the Buyer.

§ 5. Packaging and shipping, transfer of risk

(1) Packing and the choice of packaging materials shall be LI-COR's responsibility. The choice of shipping method shall be at LI-COR's discretion. If the Buyer wishes express carriage, however, it shall pay the additional costs thereby incurred.

(2) Save as provided in sub-paragraph (6) below, risk in the Products shall pass to the Buyer at the point they leave LI-COR's premises. Property in the Products remains LI-COR's and shall only pass to the Buyer upon full payment by the Buyer of all sums due (including any interest thereon). In the event of the Products being sold by the Buyer in such manner as to pass to a third party a valid title to the Products whilst any such sums are due as aforesaid, the Buyer shall be the trustee for LI-COR of the proceeds of such sale or to the claim for such proceeds and the Buyer shall place such proceeds in a separate bank account. LI-COR's rights under this paragraph shall attach to the proceeds of such sale. Nothing herein shall constitute the Buyer the agent of LI-COR for the purposes of any such sub-sale.

(3) Before full payment being made as aforesaid, LI-COR may at any time enter the Buyer's premises and repossess and remove the Products therefrom (and dispose of the same in any manner it may decide) and before such payment the Buyer shall keep such Products as fiduciary agent and bailee of LI-COR separate and identifiable for this purpose.

(4) Notwithstanding the provisions of this paragraph the Buyer shall be permitted to sell or use the Products with the prior written consent of LI-COR. Provided that such consent may be immediately terminated by LI-COR in writing at any time and shall automatically terminate if the Buyer adopts a resolution for it to be wound up or if a petition is presented for the appointment of an administrator or a receiver or an administrative receiver is appointed in respect of any part of the Buyer's undertaking or assets or if the Buyer is unable to pay its debts within the meaning of Section 121 of the Insolvency Act 1986 (or any re-enactment of further enactment thereof) unless LI-COR otherwise expressly agrees in writing to the continuation of such right of sale/use.

(5) The Buyer shall insure and keep insured the Products to their full value against all normal commercial risks until the date that property in the Products passes from LI-COR and shall whenever requested by LI-COR produce a copy of the policy of insurance.

(6) Should supply include erection and assembly by LI-COR, the risk shall transfer to the Buyer at the time of acceptance.

§ 6. 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) LI-COR's prices are always calculated in £ Sterling, the price in £ Sterling being binding even if payment of the purchase price in a currency other than £ Sterling is agreed. Changes in the exchange rate with the £ Sterling which occur after the Contract has been formed relative to the exchange rate on the date the Contract is formed 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 Buyer shall be in arrears. Whilst the Buyer is in arrears the Buyer shall pay interest on the outstanding amount before and after judgment at an annual rate of 8% above the base rate from time to time of the Bank of England (such interest to accrue on a day to day basis from the due date of payment until receipt by LI-COR of the full amount). The Buyer shall furthermore indemnify LI-COR against all costs and expenses (including reasonable legal costs and expenses on a full indemnity basis) incurred or sustained by LI-COR in recovering sums due in each case without prejudice to any other rights or remedies available to LI-COR. LI-COR also reserves the right to substantiate and claim greater damages for the arrears in individual cases.

(5) If the Buyer is in arrears as aforesaid LI-COR may furthermore suspend the delivery of Products under the Order (and any other) until payment in full of the outstanding amount (including interest accrued).

(6) Should LI-COR become aware following formation of the Contract of circumstances which suggest that the Buyer may be unable to pay, LI-COR shall be entitled to refuse to supply the Products and to set the Buyer a reasonable period in which to pay the purchase price or provide adequate security. In this event supply of the Products purchased shall take place step by step with payment of the purchase price or provision of the security in such manner as LI-COR shall require.

(7) If the Buyer is in arrears with a payment under the Contract or there are grounds to doubt its ability to pay LI-COR shall be entitled to declare any deferred invoice amounts (under the Order or any other) due immediately.

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

§ 7. Warranty

(1) LI-COR warrants that the Products shall be reasonably fit for the purpose for which they are supplied and shall conform with the Specification. All other warranties are hereby excluded to the fullest extent permitted by law.

(2) Defects must be indicated to LI-COR in writing within a period of two weeks of receipt of the Products or (in the case of latent defects) within two weeks of discovery, otherwise the right to assert of any warranty claim shall be excluded. The Buyer 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. The Buyer's sole remedy in such event shall be replacement of the defective Products or reimbursement, as the parties shall agree in writing.

(3) The warranty period shall be one year from delivery of the Products. This period and the right to claim under the warranty shall not apply if the Buyer has not notified the defect to LI-COR within the period required under sub-paragraph (2) above.

(4) LI-COR's liability shall be limited to the legal warranty with the scope described in this paragraph 7. Manufacturers' warranties in respect of the components and software used in the Products shall be unaffected by this.

(5) No warranty is given in respect of the deterioration of components of Products which occurs as a result of the usual and proper use of our equipment or normal wear and tear.

(6) The Products are intended solely for laboratory and research purposes. LI-COR shall not be liable for any damage arising from failure to observe the relevant accompanying assembly, storage, operating, maintenance and inspection instructions. Neither shall LI-COR be liable should the Buyer carry out repairs, install other parts or make other technical changes to the Products without LI-COR's written consent or use the Products contrary to LI-COR's instructions. Neither shall LI-COR be liable for damage arising from the use of the Products domestically or on humans or animals.

(7) In the event of the sale of used Products or used spare parts only the guarantees agreed in writing with authorised representatives of LI-COR shall apply. No other warranty shall be given in respect of such Products and these parts.

(8) Assignment of compensation and warranty claims by the customer to third parties is impermissible. Nothing in the Contract should be construed as conferring rights on any third party, whether under the Contracts (Rights of Third Parties) Act 1999 or otherwise and no third party shall be entitled to enforce the performance of any obligations under the Contract.

(9) LI-COR has no liability in contract, tort (including negligence) or otherwise for any loss of use, profits, revenue, goodwill, or savings, loss under current or future contracts, administrative costs or disaster recovery costs or for any consequential or indirect loss or damage of any nature under the Contract.

(10) The advice given by LI-COR in respect of the Products is given to the best of LI-COR's knowledge. However, LI-COR shall be liable for any resulting damage only in cases of grossly negligent or deliberately incorrect advice.

(11) Nothing in these Terms should be construed as excluding or limiting LI-COR's liability in negligence for death, personal injury or fraud or otherwise insofar as any exclusion or limitation of liability is void, prohibited or unenforceable by law. All of paragraph 7 should be construed as subject to this sub-paragraph (11).

(12) The Buyer shall indemnify and hold harmless LI-COR against all liabilities, losses, claims and expenses incurred by LI-COR by reason of any breach by the Buyer of the terms of the Contract or by reason of any use of the Products (whether directly or indirectly and whether in isolation or compounded with other products) by the Buyer or any person to whom the Products were supplied by the Buyer.

§ 8. Assembly

(1) Assembly of the Products shall always be carried out on the Buyer's site by LI-COR's authorised technical staff only unless LI-COR permits otherwise in writing.

(2) In the event that assembly is to be carried out by the Buyer itself the Buyer shall ensure that the technical conditions specified by LI-COR (e.g. PC system requirements) are met and shall confirm this to LI-COR in writing on request. LI-COR accepts no liability for damage arising from the fact that the necessary technical requirements are not met by the Buyer. Any additional costs and expenses arising therefrom shall be for the Buyer's account.

§ 9. Force Majeure

LI-COR shall not be liable or responsible for any loss or damage caused by delay in the performance or non-performance of its obligations hereunder due to the occurrence of circumstances beyond its control.

Standard Terms & Conditions of Sale

§10. General

(1) The Buyer shall not without the prior written consent of LI-COR assign, transfer or sub-contract the benefit or the burden of the Contract or any part thereof. LI-COR may assign or transfer its rights and obligations under the Contract.

(2) Notices under the Contract are to be served in writing and may be served by fax or first class pre-paid post to the recipient's party's address as it appears in the Contract. Notices served by fax shall be deemed received immediately following transmission (subject to issue of a valid transmission slip) and notices served by post shall be deemed to have been received 2 business days after despatch (in the case of UK post) and 6 business days after despatch (in the case of air mail).

(3) The Buyer shall have no right of set off under the Contract.

§11. Law and Jurisdiction

(1) Any controversy or claim of whatsoever nature arising out of or relating in any manner whatsoever to the Contract or any breach of any terms of the Contract shall be governed by and construed in all respects in accordance with the laws of England. The place of the Contract's performance shall be England.

(2) Each party hereby irrevocably acknowledges and agrees that the Courts of England shall have exclusive jurisdiction to resolve any controversy or claim of whatsoever nature arising out of or relating in any manner to the Contract any terms of the Contract or any breach of the Contract or any such terms.

Shipping Charges

Description	£
Standard delivery within 3 working days Without engagement. Shipping charges may vary dependent on parcel weight and size.	£25
* Courier surcharge for dangerous goods delivery Parts categorised as dangerous goods (IATA) are indicated	£85

Free Shipping for orders above £750 when ordered online

www.licor.com/bio/store

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

Notes

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

***New Empiria Studio Data Analysis Software
Free 120 Day Trial at 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



Lambda U

LI-COR[®]