

**LI-COR®***Experience Excellence*

BrightSite™

Small Animal Imaging Agents

Exceptional near-infrared fluorescent probes for targeting of tumors, bone, lymphatics, and more

Every agent is carefully validated with cultured cell assays, microscopy, *in vivo* imaging of animal models, and histology to ensure high affinity and specificity.

- Ready-to-use probes allow you to begin animal studies immediately
- Enjoy the sensitivity and signal-to-noise advantages of near-infrared fluorescence
- Simply administer the agent, and image with any small animal imaging equipment with appropriate 680 nm or 800 nm filter sets
- Compatible with most small animal imaging systems, including instruments from LI-COR Biosciences (Odyssey® and Pearl® Impulse), Caliper (CRi), Perkin Elmer (VisEn), and Carestream

Available Imaging Agents:

- **IRDye® 2-DG Optical Probe**
 - This 2-Deoxyglucose optical agent targets increased glucose metabolism and GLUT transporter expression in cancer cells. www.licor.com/2dg
- **IRDye EGF Optical Probes**
 - This probe uses recombinant human EGF to target the EGF receptor, which is overexpressed in approximately 80% of solid tumors. www.licor.com/egf
- **IRDye RGD Optical Probes**
 - The Arg-Gly-Asp tripeptide motif in this probe targets overexpression of integrins by tumor cells or during tumor angiogenesis. www.licor.com/rgd
- **IRDye PEG Contrast Agent**
 - This versatile compound is used as a general tumor contrast agent by taking advantage of the enhanced permeability and retention phenomenon. Additionally, the agent is excellent for imaging surface vasculature and lymphatics. www.licor.com/peg
- **IRDye HA (Hyaluronan) Optical Probes**
 - Hyaluronan (HA) has an important structural role in the extracellular matrix. The HA probe targets CD44 and several other receptors in the lymphatic system and lymph nodes. www.licor.com/ha
- **IRDye BoneTag™ Optical Probes**
 - These calcium-chelating probes incorporate into bone and are used to image bone structure and remodeling. Combine with a second, spectrally distinct probe for simultaneous dual-probe imaging. www.licor.com/bonetag

Mechanism of Probe Retention

P/N	Product	Description	Amt (nmol)	MW (kDa)	Dose (IV)nmol	Ex _{max}	Em _{max}	
Cell Surface Protein Binding								
926-08446	IRDye® 800CW EGF	Recombinant human EGF probe; targets overexpression of EGF receptor	20	7.3	1	779	795	
926-08946	IRDye 800CW 2-DG	Glucose analog; targets GLUT receptor overexpression, increased glucose metabolism	100	1.3	10-20	774	791	
926-09889	IRDye 800CW RGD	RGD (Arg-Gly-Asp); targets overexpression of integrins by tumor cells or during angiogenesis	15	1.7	1	776	792	
926-50001	IRDYE 800CW HA	Hyaluronan probe; targets CD44 and other receptors in lymphatic system and lymph nodes	20	8-40	2	775	794	
926-08445	IRDye 680RD EGF	Recombinant human EGF probe; targets overexpression of EGF receptor	25	7.3	2	674	698	
926-09888	IRDye 680RD RGD	RGD (Arg-Gly-Asp); targets overexpression of integrins by tumor cells or during angiogenesis	25	1.7	2	673	695	
926-50000	IRDye 680RD HA	Hyaluronan probe; targets CD44 and other receptors in lymphatic system and lymph nodes	50	8-40	4	671	695	
NEW!	929-90030	PSVue® 794	A near-infrared fluorescent probe for detection of apoptotic cells, necrotic cells, bacteria, and other anionic membranes			-	794	810

Calcium Chelating

926-09374	IRDye 680RD BoneTag™	Calcium-binding probe; images bone structure and remodeling	40	-	2	675	697
926-09375	IRDye 800CW BoneTag	Calcium-binding probe; images bone structure and remodeling	40	-	2	780	795

Vascular/Lymphatic Contrast

926-50401	IRDye 800CW PEG	Polyethylene glycol probe; general tumor contrast agent, also used to image surface vasculature and lymphatics	15	30-60	1	778	806
-----------	-----------------	--	----	-------	---	-----	-----

Cell Membrane

929-90010	CellVue® Burgundy	Lipophilic fluorescent dye; stably integrates into cell membrane	-	-	-	683	707
929-90020	CellVue NIR815	Lipophilic fluorescent dye; stably integrates into cell membrane	-	-	-	786	814

Near-Infrared Fluorophores and Labeling Kits

929-70021	IRDye 800CW NHS Ester	Labels primary and secondary amines, such as lysine residues in proteins	-	-	-	778	794
929-80021	IRDye 800CW Maleimide	Labels free sulfhydryl (-SH) groups, such as cysteine residues in proteins	-	-	-	778	794
929-09406	IRDye 800CW Carboxylate	Non-reactive form of dye; used as control for dye retention	-	-	-	778	794
929-70051	IRDye 680RD NHS Ester	Labels primary and secondary amines, such as lysine residues in proteins	-	-	-	680	694
929-71051	IRDye 680RD Maleimide	Labels free sulfhydryl (-SH) groups, such as cysteine residues in proteins	-	-	-	680	694
929-08980	IRDye 680RD Carboxylate	Non-reactive form of dye; used as control for dye retention	-	-	-	680	694
928-38040	IRDye 800CW						
	Labeling Kit, High MW	NHS ester reactive dye; for labeling of 3 X 1 mg protein (45-200 kDa)	-	-	-	778	794
928-38042	IRDye 800CW						
	Labeling Kit, Low MW	NHS ester reactive dye; for labeling of 3 X 1 mg protein (14-45 kDa)	-	-	-	778	794
928-38044	IRDye 800CW						
	Labeling Kit, Microscale	NHS ester reactive dye; for labeling of 3 X 100 µg protein (14-200 kDa)	-	-	-	778	794
928-38072	IRDye 680RD						
	Labeling Kit, High MW	NHS ester reactive dye; for labeling of 3 X 1 mg protein (45-200 kDa)	-	-	-	680	694
928-38074	IRDye 680RD						
	Labeling Kit, Low MW	NHS ester reactive dye; for labeling of 3 X 1 mg protein (14-45 kDa)	-	-	-	680	694
928-38076	IRDye 680RD						
	Labeling Kit, Microscale	NHS ester reactive dye; for labeling of 3 X 100 µg protein (14-200 kDa)	-	-	-	680	694

Application	IRDye EGF	IRDye RGD	IRDye 2-DG	IRDye PEG	IRDye HA	IRDye BoneTag	CellVue	PSVue
Tumor Imaging	•	•	•	•	•			
Metabolic Imaging			•					
Inflammation/Arthritis		•	•					
Vasculature (Contrast)				•				
Lymphatic Imaging				•	•			
Lymph Node Imaging				•	•			
Structural Imaging						•		
Cell Trafficking							•	
Apoptosis								•

LI-COR is an ISO 9001 registered company. ©2012 LI-COR, Inc. Specifications subject to change. LI-COR, IRDye, Odyssey, Pearl, BrightSite, and BoneTag are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR 4647 Superior St. • P.O. Box 4000 • Lincoln, Nebraska 68504
 LI-COR Biosciences North America: 800-645-4267 / 402-467-0700 • FAX: 402-467-0819 • Technical Support: 800-645-4260
 LI-COR GmbH, Germany: Serving Europe, Africa, and the Middle East: +49 (0) 6172 17 17 771
 LI-COR Ltd, UK: Serving UK, Ireland and Scandinavia: +44 (0) 1223 422104
 In other countries, contact LI-COR Biosciences or a local LI-COR distributor: <http://www.licor.com/distributors>
www.licor.com/bio