



Optical Agent Application Guide

LI-COR® is a leading provider of NIR optical imaging agents for small animal research. These agents are labeled with the industry leading IRDye 800CW and IRDye 680 for sensitive, early and deep detection. Developed systematically for exceptional performance, these agents offer exceptional functionality with the simplicity of injecting and imaging.

	Part Number	IRDye 800CW EGF 926-08446	IRDye 800CW RGD 926-09889	IRDye 800CW 2-DG 926-08946	IRDye 800CW PEG 926-50401	IRDye 800CW HA 926-50001	IRDye 800CW BoneTag 926-09375	IRDye 680 BoneTag 926-09376	CellVue® Burgundy 929-90010	CellVue NIR815 929-90020
Mechanism of Binding	Receptor	✓	✓			✓				
	Transporter			✓						
	Calcium Chelating						✓	✓		
	Membrane binding								✓	✓
	Non binding Contrast Agent				✓					
Applications	Tumor imaging	✓	✓	✓	✓	✓				
	Metabolic imaging			✓						
	Inflammation/Arthritis		✓	✓						
	Vasculature (Contrast)				✓					
	Lymphatic Imaging				✓	✓				
	Lymph Node Imaging				✓	✓				
	Structural Imaging						✓	✓		
	Cell Trafficking								✓	✓
Fluorescent Dye Info	Excitation Maxima	779	776	774	778	775	780	680	683	786
	Emission Maxima	795	792	791	806	794	795	705	707	814
	IRDye 800CW	✓	✓	✓	✓	✓	✓			
	IRDye 680							✓		
	NIR 815									✓
	Burgundy								✓	
	Notes	For use with cells that overexpress EGFR	For use with cells that overexpress αvβ3 integrin receptors	For use with cells with increased glycolysis	Non-binding agent for vascular, lymphatic and general tumor	For targeted lymphatic/lymph node imaging and cells over-expressing CD44.	Bone binding agent for structural imaging	Bone binding agent for structural imaging	Cell Stain	Cell Stain
Weblink: www.licor.com	/egf	/rgd	/2dg	/peg	/ha	/bonetag	/bonetag	/cellvue	/cellvue	

982-11485 01/11

Information On Ordering:

Sales Phone: 402-467-0700 or 1-800-645-4267

Fax: 402-467-0819

E-Commerce URL: <http://www.licor.com/shoponline>

Technical Support Phone: 1-800-645-4260 or 402-467-0700