

Product Number

926-50000

Storage: -20°C
prior to reconstitution;
4°C after reconstitution

Revised: July 2011

Updates available at:

<http://biosupport.licor.com>

**Limitation of Liability and
Limited Use Label License**

LI-COR IRDye optical probes are offered for research purposes only and are not intended for human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the amount of product purchased and the components of the product in research conducted by the buyer (whether the buyer is a not-for-profit, academic or for-profit entity). The buyer shall not sell or otherwise transfer this product, its components, or materials made there from to any third party. Buyer shall not use this product or its components for commercial purposes. The term "commercial purposes" shall mean any activity by a party for consideration and may include, but is not limited to, use of the product or its components (i) in manufacturing, (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label license for LI-COR IRDye infrared dyes. Inquiries regarding the licensing of one or more IRDye reagents should be submitted by e-mail to busdev@licor.com.

LI-COR DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE INFRARED DYES AND REAGENTS. LI-COR MAKES NO WARRANTIES OF ANY KIND, EITHER EXPRESS OR IMPLIED, AS TO ANY MATTER INCLUDING, BUT NOT LIMITED TO, WARRANTY OF FITNESS FOR PURPOSE, OR MERCHANTABILITY OR RESULTS OBTAINED FROM USE OF IRDYE INFRARED DYES AND REAGENTS. IN NO EVENT SHALL LI-COR BE LIABLE FOR LOST PROFITS, CONSEQUENTIAL, EXEMPLARY, SPECIAL, DIRECT, INCIDENTAL, OR PUNITIVE DAMAGES, OR ATTORNEY FEES, EVEN IF LI-COR HAD BEEN ADVISED OF, KNEW OR SHOULD HAVE KNOWN, OF THE POSSIBILITIES THEREOF. NO EMPLOYEE, AGENT OR REPRESENTATIVE OF LI-COR HAS THE AUTHORITY TO BIND LI-COR TO ANY ORAL REPRESENTATION OR WARRANTY EXCEPT AS SPECIFICALLY SET FORTH HEREIN.

© 2011 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR, Pearl, BrightSite, and IRDye are trademarks and registered trademarks of LI-COR, Inc., in the United States and other countries. Pearl Imager and IRDye reagents are covered by U.S. patents pending.

Doc # 988-11826

LI-COR

Biosciences

4647 Superior Street • P.O. Box 4000
Lincoln, Nebraska 68504 USA
North America: 800-645-4267
International: 402-467-0700
FAX: 402-467-0819

LI-COR GmbH Germany, Serving Europe,
Middle East and Africa: +49 (0) 6172 17 17 771
LI-COR UK Ltd. UK, Serving UK, Ireland, and
Scandinavia: +44 (0) 1223 422104
All other countries, contact LI-COR
Biosciences or a local LI-COR distributor:
<http://www.licor.com/distributors>

www.licor.com

BrightSite™
Small Animal Imaging Agents

IRDye® 680RD HA Optical Probe

IRDye 680RD HA (hyaluronan) Optical Probe from LI-COR® Biosciences is a near-infrared labeled optical imaging agent for small animal imaging.

Description

Hyaluronan (hyaluronic acid; HA) is an extracellular matrix glycosaminoglycan formed from disaccharide units containing N-acetylglucosamine and glucuronic acid. HA interacts with several cell surface receptors, namely CD44, RHAMM (receptor for hyaluronan mediated motility; CD168), LYVE-1 (lymphatic vessel endothelial HA receptor-1), HARE (hyaluronan receptor for endocytosis), layilin, and Toll-4. It binds to proteoglycans in cartilage and other tissues and fills an important structural role in the organization of the extracellular matrix. The agent is catabolized by receptor-mediated endocytosis and lysosomal degradation after transport via lymph to lymph nodes.

Material

The IRDye 680RD HA Optical Probe solution was passed through a 0.2 µm nylon membrane into a sterile polypropylene tube and lyophilized. The product is supplied as a powder lyophilized from water. The recommended individual dose per mouse (body weight ~25 g) is 4 nmol. Two tubes each containing 25 nmol of IRDye 680RD HA Optical Probe are provided.

Properties (in 1X PBS)

- Absorption maximum: 671 nm
- Emission maximum: 695 nm
- Appearance: Lyophilized solid

Storage and Handling

Upon receipt, immediately store at -20°C prior to reconstitution. When stored properly, this product is stable in the lyophilized state for 6 months. After reconstitution, store at 4°C. Reconstituted material should be used within 2 weeks. **Protect from light.**

Directions for Use

For systemic localization of lymph nodes: Inject 4 nmol intravenously via the tail vein. Reconstitute material in 0.625 mL of sterile 1X PBS for a final concentration of 0.04 nmol/µL. Allow the solution to sit at room temperature for 15 min to ensure the agent is completely dissolved. If desired, filter sterilize the solution through a 0.2 µm filter system.

For lymph vessel or nodal tracking in a specific region: Reconstitute the lyophilized solid in 125 µL 1X PBS to a final concentration of 0.2 nmol/µL. Inject ~ 0.6 nmol (3 µL) intradermally in region of interest.

- **In vivo Imaging:** It is important that unretained agent be allowed to clear from the animal's circulation prior to imaging. Suggested initial time frame for imaging is 24 hr post-injection; however, the timing required to achieve optimal signal-to-noise ratios will vary for each model system. IRDye 680RD HA does persist in lymph nodes for an extended period after IV administration (greater than 2 weeks). For best results, determine the optimal imaging time empirically. It is recommended the intradermal injection site be covered when imaging. This can be accomplished with the black drape provided with the Pearl® Impulse Imager.

Precautions

The probe is processed through the liver and excreted through the intestines, kidneys and bladder. This may cause increased background when imaging in the abdominal region.

Applications

Browse www.licor.com for posters, papers, and protocols related to the following applications.

- *In vitro* cell-based assays
- *In vivo* near-infrared optical animal imaging

References

1. Kuo, J.W. 2006. Practical Aspects of Hyaluronan Based Medical Products. CRC Press, Taylor & Francis Group; Boca Raton, FL.
2. Stern, R. (Ed) 2009. Hyaluronan in Cancer Biology. Academic Press, Elsevier, San Diego, CA.

Archived Document for Discontinued Product