

Product Number

**926-80000**

(2 x 250 nmol)

Storage: -20 °C

December, 2013

Updates available at:

<http://www.licor.com/packinserts>

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

LI-COR reagent products 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 a limited 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 therefrom 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 reagent products. LI-COR BIOSCIENCES DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE INFRARED DYES AND REAGENT PRODUCTS.

LI-COR BIOSCIENCES 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. 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.

© 2013 LI-COR, Inc. LI-COR is an ISO 9001 registered company. For patent information, visit [www.licor.com/patents](http://www.licor.com/patents).

LI-COR and ROSstar are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. ROSstar technology is licensed under US 2011/0070166.

Doc # 988-13229

**LI-COR**

4647 Superior Street • P.O. Box 4000  
Lincoln, Nebraska 68504 USA  
North America: 800-645-4267  
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](http://www.licor.com)

## ROSstar™ 800CW Probe

### Product Description

ROSstar 800CW is a hydrocyanine<sup>1-8</sup> based probe designed to detect extracellular reactive oxygen species (ROS) *in vivo*, specifically superoxide and hydroxyl radicals. The cell-impermeable probe can be used to detect oxidative stress using fluorescence imaging when using instrumentation with appropriate excitation and detection capabilities.

### Properties

**Concentration:** 1 nmol/μL (after reconstitution)

**Excitation Wavelength:** 780 nm

**Emissions Wavelength:** 810 nm

**Form:** Lyophilized

**Pack Size:** 2 x 250 nmol

**Appearance:** yellow to orange solid

### Storage and Handling

**Protect from light.** Upon receipt, immediately store at -20 °C. This product is stable in the lyophilized state for 3 months at -20 °C. Reconstituted product should be used immediately.

### Precautions

- ROSstar 800CW is light- and air-sensitive
- Open just before use – **Discard unused material**

### General Guidelines for Use

- Reconstitute material (1 tube containing 250 nmol) in 250 μL sterile water or 1x PBS for a final concentration of 1.0 nmol/μL. If desired, filter sterilize the solution through a 0.2 μm filter system prior to animal injection.
- The recommended individual dose per mouse (body weight ~20 g) will range from 30 - 50 nmol.  
*The dose may need further optimization, depending upon factors including animal model, mode of administration, instrument used for detection, and specialized applications.*
- Image 15 - 30 minutes after the injection of ROSstar 800CW.  
*Imaging optimization may be needed, depending upon factors including animal model, mode of administration, instrument used for detection, and specialized applications*
- For further information and applications of hydrocyanines in *in vivo* imaging of ROS, see references 1 - 8.

## Hydrocyanine References:

1. Kundu, K.; Knight, S. F.; Willett, N.; Lee, S.; Taylor, W. R.; Murthy, N. "Hydrocyanines: a class of fluorescent sensors that can image reactive oxygen species in cell culture, tissue, and *in vivo*", *Angewandte Chemie, International Edition* (2009), *48*(2), 299-303.
2. Withers, N. "Fluorescence imaging: Go with the glow", *Nature Chemistry*, (19 December 2008).
3. Kundu, K.; Knight, S. F.; Lee, S.; Taylor, W. R.; Murthy, N., "A Significant Improvement of the Efficacy of Radical Oxidant Probes by the Kinetic Isotope Effect", *Angewandte Chemie, International Edition* (2010), *49*(35), 6134-6138.
4. Selvam, S.; Kundu, K.; Templeton, K.; Murthy, N.; Garcia, A. "Minimally Invasive, Longitudinal Monitoring of Biomaterial-Associated Inflammation by Fluorescence Imaging", *Biomaterials*, (2011), *32*, 7785-7792.
5. Kim, J. Y.; Choi, W.; Kim, Y. H.; Tae, G. "Highly Selective *In Vivo* Imaging of Tumor as an Inflammation Site by ROS Detection Using Hydrocyanine-Conjugated, Functional Nano-Carriers", *Journal of Controlled Release* (2011), *156*, 398-405.
6. Xie, L.; Lin, A.; Kundu, K.; Murthy, N.; Lavenston, M.; Guldborg, M. "Quantitative Imaging of Cartilage and Bone Morphology, Reactive Oxygen Species, and Vascularization in A Rodent Model of Osteoarthritis", *Arthritis and Rheumatism*, (2012), *64*(6), 1899-1908.
7. Goodson, P.; Kumar, A.; Jain, Lucky; Kundu, Kousik; Murthy, Niren; Koval, Michael; Helms, My N. "Nadph oxidase regulates alveolar epithelial sodium channel activity and lung fluid balance in vivo via O<sup>2</sup> signaling", *American Journal of Physiology* (2012), *302*(2, Pt. 1), L410-L419.
8. Magalotti, S.; Gustafson, T. P.; Cao, Q.; Abendschein, D. F.; Pierce, R. A.; Berezin, M. Y.; Akers, W. J. Evaluation of Inflammatory Response to Acute Ischemia Using Near-Infrared Fluorescent Reactive Oxygen Sensors", *Molecular Imaging and Biology*, (2013), *15*(4), 423-430.

Archived Document for Research Product