

Product Number

**928-40014**

Quantity: 4 L

Storage: RT

Revised: September 2010

Updates available at:

<http://biosupport.licor.com>

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

LI-COR IRDye<sup>®</sup> Infrared Dyes and 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 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 and reagent products. Inquiries regarding the licensing of one or more IRDye infrared dyes should be submitted by e-mail to busdev@licor.com.

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.

© 2006 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR, IRDye and Odyssey are registered trademarks of LI-COR, Inc., in the United States and other countries. All other trademarks belong to their respective owners.

Doc # 988-08895

**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 and  
Africa: +49 (0) 6172 17 17 771  
LI-COR UK Ltd. UK, Serving UK, Ireland,  
Scandinavia: +44 (0) 1223 422104  
[www.licor.com](http://www.licor.com)

## Tris Glycine/Sodium Dodecyl Sulfate Buffer, 10X Liquid Concentrate

### Product Description

A 1X working solution (1:10 dilution of the 10X liquid concentrate) contains 0.025 M Tris base, 0.192 M Glycine, and 0.1% sodium dodecyl sulfate.

### Appearance

Clear, colorless liquid.

### Specifications

pH (1X, Water) at 25°C: 8.30 – 8.70.

### Storage

Store in provided container at room temperature; protect from moisture.

Expiration is one year from date received.

### Hazard

Irritant

Archived Document for Discontinued Product