

Limitation of Liability and Limited Use Label License

LI-COR IRDye® 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 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. Buyer shall not determine the structure or otherwise reverse engineer this product. 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 DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE INFRARED DYES AND REAGENT PRODUCTS. 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. 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.

IRDye 800CW, IRDye 700DX, IRDye 800PA, IRDye 700PA and IRDye infrared dyes are covered by U.S. and foreign patents and patents pending. LI-COR, IRDye, Odyssey, Aeries, and In-Cell Western are trademarks or registered trademarks of LI-COR, Inc.

LI-COR 4647 Superior Street • P.O. Box 4000 • Lincoln, Nebraska 68504 USA
Biosciences Technical Support: 800-645-4260 • North America: 800-645-4267 •
International: 402-467-0700 • LI-COR GmbH*: +49 (0) 6172 17 17 771
LI-COR UK Ltd.: +44 (0)1223 422104 • www.licor.com

* Serving Austria, Czech Republic, Germany, Hungary, Slovakia, and Switzerland

Doc #988-08897

LI-COR® IRDye®

Infrared Imaging Reagents

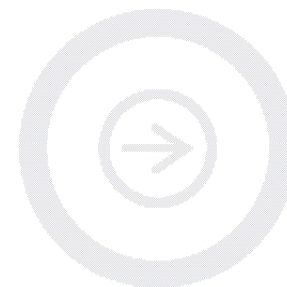
2X Protein Loading Buffer

Product Number: 928-40004

Quantity: 30 mL

Shipping: RT

Storage: RT



Published October 2006. The most recent version of this pack insert is posted at <http://biosupport.licor.com/support>.

LI-COR
Biosciences

FORMULATION

62.5 mM	Tris-HCl, pH 6.8
25%	Glycerol
2%	SDS
0.1% (w/v)	Orange G

USE

2X Protein Loading Buffer is used as a loading buffer for protein gel electrophoresis. This loading buffer is not visualized when imaged on the Odyssey® Infrared Imaging System. The buffer is optimized for use with SDS-PAGE and Tris-glycine-SDS running buffer.

INSTRUCTIONS FOR USE

1. Add reducing agent
 - a. Add 50 μ l of β -mercaptoethanol to 950 μ l of Protein Loading Buffer.
 - b. Do not store Protein Loading Buffer with β -mercaptoethanol for more than two weeks.
2. Dilute Sample
 - a. The recommended concentration for electrophoresis is 1X.
 - b. For every 1 μ l of sample, add 1 μ l of Protein Loading Buffer.

STORAGE

Keep in a tightly sealed tube.

Store at room temperature.

Expiration is one year from date received..