

Product Number

829-05352

(IRDye® 700 & IRDye 800)

Quantity: 100 - 125 Lanes

Storage: -20°C

Revised: February 2012

Updates available at:

<http://biosupport.licor.com>

**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 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.

© 2012 LI-COR, Inc. LI-COR is an ISO 9001 registered company. IRDye infrared dyes are covered by U.S. and foreign patents and patents pending. LI-COR, and IRDye are registered trademarks of LI-COR, Inc., in the United States and other countries.

Doc # 988-12812



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/bio

50-1500 bp Two-Color DNA Sizing Standard

Product Description

The two-color 50-1500 bp size standard is composed of 15 IRDye® DNA fragments, provided in a lyophilized form, that can be detected simultaneously in the 700 nm and 800 nm channels of a LI-COR DNA Analyzer. The DNA fragment sizes were selected to allow rapid identification of bands and accurate sizing over the entire range. Optional reconstitution volumes provide flexibility for the formulation that is best suited for your application.

Components

IRDye 700 DNA Size Standards

- 50-1500 bp size standard (2 tubes, lyophilized)
- Blue Stop Solution (1 tube, 0.25 mL) – contains 90% formamide in bromophenol blue

IRDye 800 DNA Size Standards

- 50-1500 bp size standard (2 tubes, lyophilized)
- Blue Stop Solution (1 tube, 0.25 mL) – contains 90% formamide in bromophenol blue

Protocol

Recommended Resuspension Volume

Desired Load Volume (µL / well)	Resuspension Volume (µL / tube)
1.0	50
0.8	40
0.6	30
0.5	25
0.4	20
0.2	10

- Resuspend one tube of IRDye 700 DNA size standard in desired **Resuspension Volume**, mix well and spin briefly.
- Transfer the entire volume into one tube of the IRDye 800 DNA size standard. Resuspend carefully by mixing and briefly spin again.
- Remove a small aliquot, denature the product at 95°C for 1-2 minutes.
- Place on ice for at least 10 minutes prior to loading.
- Discard any remaining denatured, small aliquots after loading.
- Store solution at -20°C until next use.

Load Volume Considerations

Comb type	Load volume (µL / well)
48 well	0.4 – 0.8
64 well	0.2 – 0.5

Stability

The lyophilized material is stable at -20°C for 12 months from the date of receipt. Once reconstituted as directed above, the material is stable at -20°C for 3 months.

Precautions

Protect from light.

