

Product Number

4200-60 (IRDye® 700)

4000-45 (IRDye 800)

Quantity: 100 µL

Storage: -20°C

Revised: March, 2006

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.

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

Doc #988-06994

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

50-700 BP Sizing Standard

Product Description

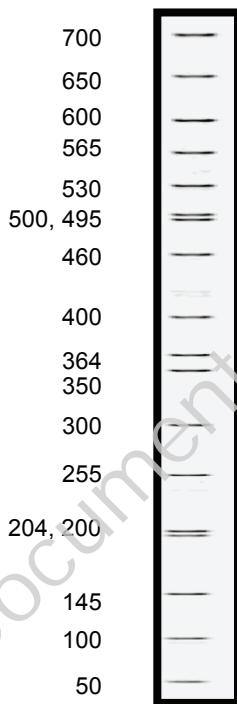
The 50-700 bp size standard is composed of 18 IRDye®-labeled DNA fragments with equal banding intensities in 90% formamide solution with bromophenol blue. The fragments cover a size range of 50 to 700 bp (see figure below). The fragment sizes were selected to allow rapid identification of bands and accurate fragment sizing over the entire range with image analysis programs like Gene ImagIR.

Protocol

- Immediately prior to use, the product should be denatured at 95°C for 1-2 minutes, then placed on ice for 10 minutes.
- Load 1 µL for 48 well combs and 0.8 µL for 64-well combs.

Precautions

- The IRDye labeled DNA should be stored in darkness. For best performance, minimize exposure to light.
- This sizing standard is not recommended for 41 cm or 66 cm gels; in these cases, higher molecular weight bands may exhibit multiple bands.



Licensing

The size standard is produced under license using PCR technology described in U.S. Patent 4,683,202 owned by Roche Molecular Systems, Inc. No license to use the PCR process is conveyed expressly or by implication to the purchaser by the purchase of this product.

These products are offered for research and development purposes only. They cannot be used to manufacture derivatives or products for re-sale.