

Product Number
829-05570

M13 Forward (-29) IRDye® 700 Primer

For LI-COR DNA Analysis Systems

Storage: -20°C

Revised: February 2010

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.

© 2007 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR and IRDye are registered trademarks of LI-COR, Inc., in the United States and other countries. All other trademarks belong to their respective owners.

Doc #988-06984

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

Dilution Procedure

Add 1000 µl sterile ice-cold 1XTE Buffer to the 1.0 nmol of lyophilized primer. A tube of 1XTE Buffer is included with your primer order. Mix well until all of the primer is dissolved (un-dissolved primer should be easily visible as a dark blue pellet). Aliquot into 20 sterile microcentrifuge tubes at 50 µl/tube.

This dilution will produce a primer solution with a 1.0 µM (1.0 pmol/µl) concentration. This primer solution may be used directly in LI-COR sequencing protocols.

1XTE Buffer Solution

The 1XTE included with the primer is 10 mM Tris, 1 mM EDTA, pH = 7.4 (± .2). Additional 1XTE can be made by diluting 10 µl of 100XTE Buffer (Sigma Chemical Company, Catalog #T9285) to a 1 mL total volume with sterile distilled water. Store at 4°C.

Measuring Concentration

The concentration of the primer solution may be verified by measurement of a dilution of the final solution on a UV/Vis absorption spectrophotometer equipped with a microcell. Since EDTA interferes with measurement of absorbance at 260 nm, the concentration should be determined based on the IRDye 700 peak at 685 nm. The molar absorptivity of IRDye 700 at 685 nm in aqueous solution is 170,000. The concentration may be calculated using the following formula:

$$\frac{(\text{Absorbance at 685 nm}) \times (\text{Dilution Factor})}{(170,000 \times \text{Cell path length (cm)})} = \text{Molar concentration of the original solution}$$

Example: If, using a 1 cm path length cuvette, the absorbance at 685 nm equals 0.10 for a 1:2 dilution of the final primer solution, the concentration of the original solution can be determined as follows:

$$(0.10 \times 2) / (170,000 \times 1) = 1.2 \times 10^{-6} \text{ M} = 1.2 \mu\text{M}$$

Storage Conditions

Minimize exposure to light. Store tubes at -20°C in a light-tight container or foil shipping bag. Minimizing freeze/thaw cycles may help preserve the integrity of the fluorescent primer.

Sequence: 5'-CAC GAC GTT GTA AAA CGA C-3'

Tm: 58.1°C

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