

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, Aerius, and In-Cell Western are trademarks or registered trademarks of LI-COR, Inc.

LI-COR and IRDye are trademarks and registered trademarks of LI-COR, Inc. Expedite is a trademark of Applied Biosystems, Inc. All other trademarks belong to their respective owners.

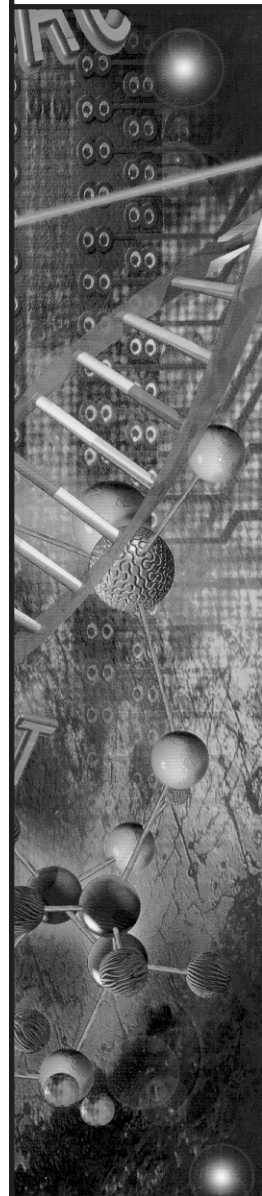
This product is sold for research purposes only. Any commercial use or resale of this product or derivative products (e.g., labeled primers) requires a separate license from LI-COR, inc. LI-COR and/or University of Nebraska patent numbers: U.S., 4,729,947; 5,207,880; 5,230,781; 5,571,388; 5,366,603; 5,360,523; 5,549,805; Canada, 1,230,161; Australia, 662,273; 670,622; Europe, 157,280; 533,302; Japan, 2007209. Other U.S. and foreign patent applications pending. All trademarks are the property of their respective owners.

LI-COR® 4647 Superior Street • P.O. Box 4000 • Lincoln, Nebraska 68504 USA
Biosciences
North America: 800-645-4267 • Technical Support: 800-645-4260
International: 402-467-0700 • FAX: 402-467-0819
E-mail: biohelp@licor.com • www.licor.com

Doc # 988-06938
Rev. 5 0609

DNA Synthesis With IRDye® 800 Phosphoramidite

Part #4000-33



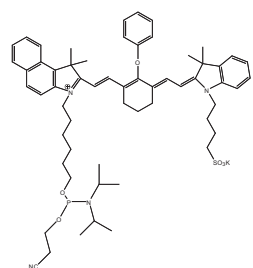
LI-COR®
Biosciences

Principle and Scope

The phosphoramidite of the Near IR fluorescent dye, IRDye® 800, can be used to label DNA molecules prepared in a DNA synthesizer. The dye is attached to the 5' end of the protected, support-bonded oligonucleotide via fast deprotection phosphoramidite chemistry. On syntheses at 200 nmol scale, typical crude yields of dye-labeled oligonucleotides are 65 to 100 nmol. Standard phosphoramidites and deprotection conditions cannot be used. This protocol was developed on an Applied Biosystems 8909 DNA synthesizer and works well using the Polygen DNA Synthesizer. For best results on other systems, protocol modification may be required.

Structure and Properties

Fig. 1. Structure of IRDye 800 Phosphoramidite



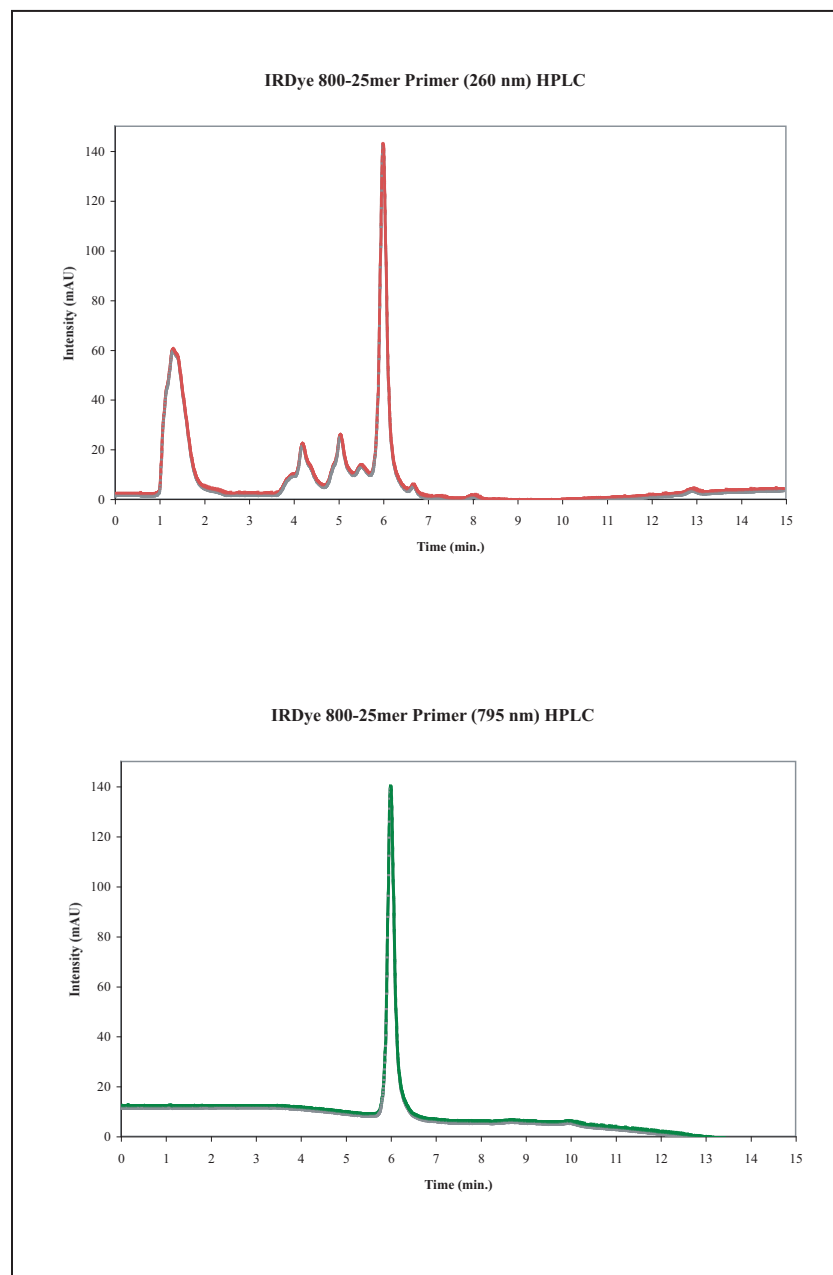
IRDye 800 Phosphoramidite

Absorption Max (λ_{max} , methanol)	787 nm
Absorption Max (aqueous oligonucleotide)	795 nm
Molar Absorptivity (ϵ , methanol)	270,000
Emission Max (methanol)	812 nm
Emission Max (aqueous oligonucleotide)	819 nm
Molecular Weight	999.30

Equipment and Reagents

- IRDye 800 phosphoramidite (LI-COR Biosciences, 4000-33)
- Polygen DNA-Synthesizer or Equivalent
- CPG column/beads, 500A, 200 nanomole (Proligo A322004-04, C322004-01, G322004-01, T321010-00, A302001-01, C302001-01, G302001-01, T301001-01 or equivalent)
- A, C, G, and T phosphoramidites (Proligo A112081-12, C112081-12, G112081-12, T111081-12 or equivalent)
- Synthesis Reagents: Oxidizing solution, Capping reagent solutions, Deblocking reagent solution, Activator solution, Acetonitrile (Glen Research 40-4132-52, Proligo L870020-06, L850020-4N, L820090-06, L880020-06, L810045-06 or equivalent)
- Ammonium Hydroxide
- Microcentrifuge, standard and filter tubes
- Gradient HPLC system, reverse phase C18, for product analysis and purification, if desired. Acetonitrile (HPLC grade), triethylammonium

Fig. 3 Typical Analytical HPLC of Crude Oligonucleotide



5. Primer Analysis and Purification by HPLC

The crude oligonucleotides are sufficiently pure for many purposes. Their purities can be checked with analytical HPLC, and preparative HPLC can be used when highest purity is needed. It is convenient to monitor at least two wavelengths with the HPLC detector. Detection at 795 nm shows just the dye-labeled components in the sample, while 260 or 280 nm will also reveal unlabeled products, such as the failure sequences of the synthesis. These HPLC conditions can be used for analysis and purification.

Solvents: A = 0.1 M triethylammonium acetate in
4% acetonitrile/96% water
B = 0.1 M triethylammonium acetate in
80% acetonitrile/20% water
Column: C18, 5 μm , 300 \AA , 10 \times 100 mm
Flow: 3.4 ml/min
Gradient: 30–40% B over 0.3 minutes
40–100% B from 0.3 – 8 minutes
100% B for 2 minutes
100–30% B for 5 minutes

Analytical HPLC. Dilute 1 μl of the filtered, deprotected oligonucleotide solution from step 3 with 99 μl of sterile, pure water. Inject 10 μl of the diluted solution for the analytical run. Typical chromatograms for both wavelengths are shown in Figure 3 for an oligonucleotide prepared with standard deprotection chemistry. The main peak (~6 min) is the desired product. Synthesis failure sequences are more hydrophilic, so these components elute earlier.

Preparative HPLC. If LC purification is desired, all or part of the filtered, deprotected oligonucleotide solution from step 3 is directly injected. Collect appropriate fractions and remove the solvents and the triethylammonium acetate buffer in a vacuum drying unit.

6. Solution Makeup

Dissolve the oligonucleotide in 1xTE (10 mM Tris, 1 mM EDTA, pH 7.4) to the desired concentration. For a 1.0 μM (1.0 pmol/ μL) solution, dilute the dye labeled oligonucleotide to obtain 0.27 absorbance units

Precautions

1. The oxidizing solution supplied by Glen Research (40-4132-52 or equivalent*) should be used. Harsher oxidizer solutions supplied by some manufacturers are not recommended.
2. All phosphoramidites are sensitive to moisture and oxygen, and the handling procedures described here are good practice in general. No special modifications of your standard procedures for similar products are likely to be necessary for IRDye 800 phosphoramidite. Protect the product from air and moisture. Use only dry, oxygen-free acetonitrile ("DNA synthesis grade") for preparation of the amidite solutions.
3. For best results, use the entire solution of IRDye 800 amidite within a day. There is some loss of coupling efficiency of the solutions on standing. If the solutions will not be used within a day, place the bottle in a jar with some Drierite and store at -20°C to prolong the working life of the amidite.
4. Flushing the lines of the synthesizer with acetonitrile (or methanol followed by acetonitrile) is generally sufficient to remove any traces of the dye.
5. The ammonia solution used in the deprotection of the oligonucleotides should be handled in a fumehood.

* Recipe: 4.3 g iodine, 90 ml water, 4 ml pyridine, and 900 ml THF

Procedure

1. Prepare a 0.1 M Solution of the IRDye 800 Phosphoramidite in Acetonitrile

- a. Add 1.00 ml of acetonitrile to the 100 μmol package to make the solution. It is good practice to provide a source of dry nitrogen or argon gas to displace the solvent that will be removed from the acetonitrile bottle. A needle connected to a stream of the gas and a bubbler is a convenient way to do this. Another approach is to use a second syringe barrel filled with desiccant for the vent, although this does not remove oxygen. Use a similar setup to allow for the escape of the gas displaced when the acetonitrile is added to the phosphoramidite vial.
- b. Choose a clean, dry syringe and needle of appropriate size. Pierce the seal of the acetonitrile reagent bottle with the needle and slowly withdraw slightly more than the desired amount of solvent.
- c. Do the next steps rapidly to minimize exposure of the acetonitrile to air. Remove the syringe and needle from the acetonitrile bottle. Invert the syringe and force any air bubbles out the needle. Push the excess acetonitrile from the syringe into an appropriate waste container. Inject the acetonitrile into the sealed IRDye 800 phosphoramidite vial. Remove the needles from the seals, and allow the IRDye 800 phosphoramidite to dissolve. Swirl the solution occasionally until no undissolved particles are observed on the sides of the vial.

2. DNA Synthesis Program

The basic synthesis protocol for your synthesizer can be used until the last step involving the dye amidite. A longer coupling time for the IRDye 800 phosphoramidite is recommended. A 5-10 minute increase in standard coupling time will improve final yields. Since the IRDye 800 phosphoramidite does not contain any trityl group, a "DMT On" mode should be used to avoid unnecessary exposure of the dye to the detritylation reagent.

3. Primer Deprotection

Note: All work with ammonium hydroxide (NH₄OH) should be performed in a fume hood. Methylamine deprotection methods do not work with IRDye phosphoramidites.

Place CPG beads from the synthesis column in a microcentrifuge tube (1.5–2.0 mL capacity). Add 400 µL of cold NH₄OH to each tube. Close tubes and mix the contents thoroughly on a vortex mixer. Allow the samples to stand in the dark for 1 hour at room temperature. Add an additional 400 µL of cold NH₄OH to each tube, close and mix thoroughly. Allow the samples to stand an additional 30 minutes in the dark at room temperature. Strip off the aqueous ammonia in a vacuum drying unit, then resuspend the dry pellet in 200 µL of sterile, ultrapure water. Decant solution from the beads or use a syringe filter to remove CPG beads from the solution and transfer to a new tube.

4. Measurement of Crude Synthesis Yield (Optional)

Dilute 1.0 µL of the filtered primer solution with 199.0 µL of sterile ultrapure water. Obtain the UV/vis absorption spectrum of the resulting solution (see Figure 2). The crude yield can then be calculated from the absorbance value at 795 nm (A_{795}), the absorptivity (270 µL / nmol / cm), the path length (b cm), the dilution factor (200) and the total primer solution volume (200 µL):

$$\text{Yield (nmol)} = A_{795} / (270 \times b) \times 200 \times 200$$

For example, a measured absorbance of 0.46 in a 1 cm cell would correspond to a crude yield of 68 nmol.

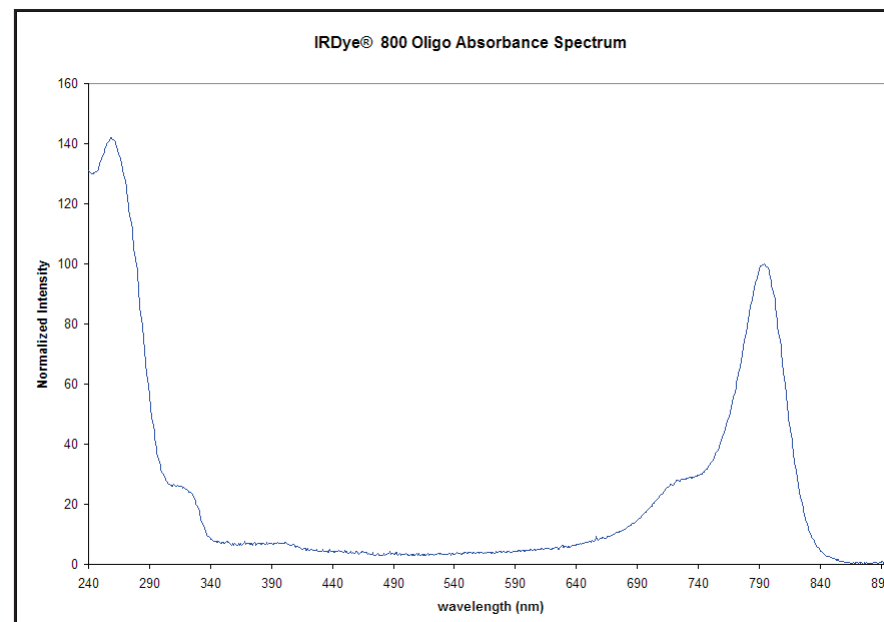


Fig. 2 Typical UV/Vis Spectrum