

EZ-Link[®] Psoralen-PEG₃-Biotin

29986

0744.4

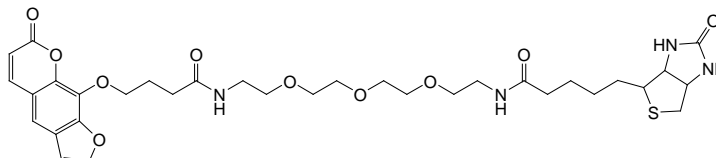
Number

29986

Description**EZ-Link Psoralen-PEG₃-Biotin, 5mg**

Molecular Weight: 688.79

Spacer Arm: 36.86Å



Storage: This product is shipped at ambient temperature. Upon receipt store product at 4°C protected from light and moisture.

Introduction

Thermo Scientific EZ-Link Psoralen-PEG₃-Biotin contains a water-soluble polyethylene glycol (PEG) spacer arm that produces better solubility of biotinylated molecules than aliphatic long-chain spacers. The extended spacer provides maximum accessibility for detection of molecules labeled with biotin binding proteins. Unlike other aliphatic, long-chain biotinylation reagents, the PEG spacer will not become buried or hidden in hydrophobic pockets. Also, modifications performed with this unique biotin derivative will not decrease the solubility of biotinylated molecules, but actually promotes conjugate solubility.

The photoreactive psoralen group provides much greater covalent insertion yields than phenyl azide-containing compounds, making this reagent a superior alternative to typical photoactivatable biotin derivatives. The psoralen tricyclic planar ring system intercalates into double-stranded DNA, RNA, PCR products, cDNA, oligonucleotides and to a lesser extent single-stranded DNA. Psoralen-PEG₃-Biotin can be used with gel-purified oligonucleotides because agarose does not interfere with conjugation. High-yield conjugation occurs after a brief exposure to UV light (>350nm, 10-30 minutes). The nucleic acid labeling site of the psoralen group (the 5,6 double bond in thymine residues) does not interfere with subsequent hybridization reactions, creating a highly sensitive direct labeling alternative to radioactive probes.

Example Procedure for Biotinylating DNA

The following protocol is an adaptation from Henriksen, *et al.* For best results, empirically optimize each specific application and experimental system.

1. Adjust DNA or RNA to the desired concentration (e.g., 20-100µg/mL) in TE buffer (10mM Tris, 1mM EDTA, pH 7.4).

Note: The psoralen group can crosslink the two strands of double-stranded DNA. Before labeling, heat-denature double-stranded DNA to produce single-stranded DNA. To denature, boil DNA for 5 minutes and then place tube in a dry ice/ethanol bath to quickly cool the denatured DNA.

2. In subdued indirect light, dissolve Psoralen-PEG₃-Biotin in DMF or DMSO to 20mM (~1.38mg/100µL).
3. Add the biotin solution to the DNA or RNA and mix well. In a standard reaction, the final concentration of Psoralen-PEG₃-Biotin is ~200µM (1:100 dilution of stock solution prepared in Step 2). A serial dilution of the Psoralen-PEG₃-Biotin stock solution may be required when using small reaction volumes.
4. Irradiate the open reaction tube from above using a long wavelength UV light source for at least 10-30 minutes. If desired, keep the sample cooled on ice during irradiation.
5. To remove non-reacted biotin, precipitate sample with 0.2M potassium acetate and two volumes of ethanol. After centrifugation, wash pellet with 70% ethanol and allow it to dry. Dissolve the biotinylated sample in water or buffer.

Determination of Biotin Incorporation

Biotin incorporation can be estimated using the HABA (4'-hydroxyazobenzene-2-carboxylic acid) method. In solution, the HABA dye binds avidin, forming a complex with maximal absorption at 500 nm. When biotin is added to the solution, its higher affinity for avidin displaces the HABA and the absorption at 500 nm decreases proportionately. The absorbance of the HABA-avidin solution is measured before and after adding the biotin-containing sample. The change in absorbance relates to the amount of biotin in the sample. The Thermo Scientific Pierce Biotin Quantitation Kit (Product No. 28005) contains a biotinylated protein control and a premix of HABA and avidin supplied in convenient No-Weigh™ Microtube packaging.

Related Thermo Scientific Products

95034, 95035	3UV Lamps (254nm, 302nm, 365nm settings)
69576	Slide-A-Lyzer® MINI Dialysis Unit Kit, 10K MWCO, 0.1mL, 10/pkg
66382, 66807	Slide-A-Lyzer Dialysis Cassette Kits, 10K MWCO, for 0.5-3mL and 3-12mL sample volumes, respectively
28005	Pierce® Biotin Quantitation Kit
20347	Streptavidin Agarose Resin, 2mL
20228	Pierce Monomeric Avidin Kit
21115	Pierce Biotinylated Protein Interaction Pull-Down Kit
21126	Streptavidin, Horseradish Peroxidase Conjugated, 1mg
21324	Streptavidin, Alkaline Phosphatase Conjugated, 1mg
15120	Streptavidin Coated Strip Plates, 5 plates
21330	EZ-Link NHS-PEG ₄ -Biotin, 25mg
21331	EZ-Link Sulfo-NHS-SS-Biotin, 100mg
21325	EZ-Link NHS-Chromogenic Biotin, 10mg

General References

Henriksen, U., *et al.* (1991). Azidobenzoyl-, azidoacridinyl-diazocyclopentadienyl-carbonyl-, and 8-propyloxypsoralen photobiotinylation reagents. Syntheses and photoreactions with DNA and protein. *Photochem Photobiol A Chem* **57**:331-42.

Oser, A., *et al.* (1988). Sensitive non-radioactive dot-blot hybridization using DNA probes labeled with chelate group substituted psoralen and quantitative detection by europium ion fluorescence. *Nucl Acid Res* **16**:1181-96.

Wassarman, D.A. (1993). Psoralen crosslinking of small RNAs *in vitro*. *Molec Biol Reports* **17**:143-51.

Product Reference

Zhang, Y., *et al.* (2001). Reproducible and inexpensive probe preparation for oligonucleotides arrays. *Nucl Acid Res* **29**(13):E66-6.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.