

Biotin PEG4 succinimidyl ester

Catalog number: 3022
Unit size: 25 mg

Component	Storage	Amount (Cat No. 3022)
Biotin PEG4 succinimidyl ester	Freeze (< -15 °C)	1 vial (25 mg)

OVERVIEW

Biotin PEG4 succinimidyl ester is an amine-reactive biotin derivative that contains a long arm to increase its avidin-binding affinity. This PEG biotin reacts specifically and efficiently with primary amines (e.g., a side chain of lysine residues or aminosilane-coated surfaces) at pH 7-9 to form covalent bonds. The PEG spacer arm gives the reagent a long and flexible connection to minimize steric hindrance involved with binding to avidin molecules. The hydrophilic polyethylene glycol (PEG) spacer arm of this reagent imparts water solubility that is transferred to the labeled molecule, thus reducing aggregation of labeled proteins stored in solution. It is widely used to label a variety of biological molecules and samples.

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles

Protein stock solution (Solution A)

Mix 100 µL of a reaction buffer (e.g., 1 M sodium bicarbonate solution or 1 M phosphate buffer with pH ~9.0) with 900 µL of the target protein solution (e.g. antibody, protein concentration >2 mg/mL if possible) to give 1 mL protein labeling stock solution.

Note: The pH of the protein solution (Solution A) should be 8.5 ± 0.5 . If the pH of the protein solution is lower than 8.0, adjust the pH to the range of 8.0-9.0 using 1 M sodium bicarbonate solution or 1 M pH 9.0 phosphate buffer.

Note: The protein should be dissolved in 1X phosphate buffered saline (PBS), pH 7.2-7.4. If the protein is dissolved in Tris or glycine buffer, it must be dialyzed against 1X PBS, pH 7.2-7.4, to remove free amines or ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for protein precipitation.

Note: Impure antibodies or antibodies stabilized with bovine serum albumin (BSA) or gelatin will not be labeled well. The presence of sodium azide or thimerosal might also interfere with the conjugation reaction. Sodium azide or thimerosal can be removed by dialysis or spin column for optimal labeling results.

Note: The conjugation efficiency is significantly reduced if the protein concentration is less than 2 mg/mL. For optimal labeling efficiency the final protein concentration range of 2-10 mg/mL is recommended.

Biotin-PEG4 succinimidyl ester stock solution (Solution B)

Add anhydrous DMSO into the vial of Biotin-PEG4 succinimidyl ester to make a 10 mM stock solution. Mix well by pipetting or vortex.

Note: Prepare the Biotin-PEG4 succinimidyl ester (Solution B) before starting the conjugation. Use promptly. Extended storage of the Biotin-PEG4 succinimidyl ester stock solution may reduce its activity. Solution B can be stored in freezer for two weeks when kept from light and moisture. Avoid freeze-thaw cycles.

SAMPLE EXPERIMENTAL PROTOCOL

This labeling protocol was developed for the conjugate of Goat anti-mouse IgG with Biotin-PEG4 succinimidyl ester. You might need further optimization for your particular proteins.

Note: Each protein requires distinct Biotin-PEG4/protein ratio, which also depends on the properties of Biotin-PEG4. Over labeling of a protein could detrimentally affects its binding affinity while the protein conjugates of low Biotin-PEG4/protein ratio gives reduced sensitivity.

Run conjugation reaction

1. Use 10:1 molar ratio of Solution B (Biotin-PEG4)/Solution A (protein) as the starting point: Add 5 µL of the Biotin-PEG4 stock solution (Solution B, assuming the Biotin-PEG4 stock solution is 10 mM) into the vial of the protein solution (95 µL of Solution A) with effective shaking. The concentration of the protein is ~0.05 mM assuming the protein concentration is 10 mg/mL and the molecular weight of the protein is ~200KD.

Note: We recommend to use 10:1 molar ratio of Solution B (Biotin-PEG4)/Solution A (protein). If it is too less or too high, determine the optimal Biotin-PEG4/protein ratio at 5:1, 15:1 and 20:1 respectively.

2. Continue to rotate or shake the reaction mixture at room temperature for 30-60 minutes.

Purify the conjugation

The following protocol is an example of dye-protein conjugate purification by using a Sephadex G-25 column.

1. Prepare Sephadex G-25 column according to the manufacture instruction.
2. Load the reaction mixture (From "Run conjugation reaction") to the top of the Sephadex G-25 column.
3. Add PBS (pH 7.2-7.4) as soon as the sample runs just below the top resin surface.
4. Add more PBS (pH 7.2-7.4) to the desired sample to complete the column purification. Combine the fractions that contain the desired Biotin-PEG4-protein conjugate.

Note: For immediate use, the Biotin-PEG4-protein conjugate need be diluted with staining buffer, and aliquoted for multiple uses.

Note: For longer term storage, Biotin-PEG4-protein conjugate solution need be concentrated or freeze dried.

EXAMPLE DATA ANALYSIS AND FIGURES

Calculate DOS

You can calculate DOS using our Amplite® Colorimetric Biotin Quantification Kit (Cat# 5522).

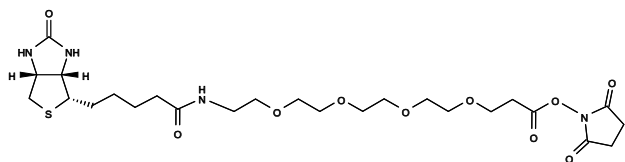


Figure 1. Chemical structure for Biotin PEG4 succinimidyl ester

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.