

## D-Biotin-PEG3, SE

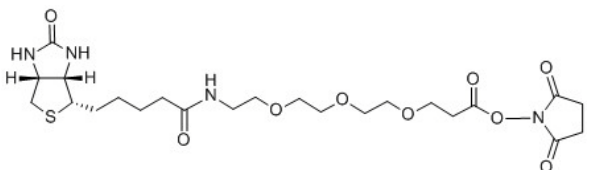
Catalog Number	Packaging Size
<b>B006</b>	<b>25 mg</b>

**Storage upon receipt:** -20°C

### Introduction

The amine-reactive **biotin-PEG3, SE** is a pegylated, water-soluble reagent for simple and efficient biotin labeling of antibodies, proteins and other primary amine-containing macromolecules for subsequent detection with streptavidin, avidin or NeutrAvidin® biotin-binding protein. This molecule has a PEG3 spacer to improve water solubility and facilitate reactivity with the molecule and eventual accessibility of the biotin to avidin or streptavidin.

### Specifications

<b>Molecular Formula:</b>	C <sub>23</sub> H <sub>36</sub> N <sub>4</sub> O <sub>9</sub> S	
<b>Molecular Weight:</b>	544.62	
<b>CAS Number:</b>	-	
<b>Storage Conditions:</b>	-20°C	
<b>Shipping Condition:</b>	Room Temperature	

### Applications

Biotinylation reagent

### References:

1. A method for gene enrichment based on the avidin-biotin interaction. Application to the Drosophila ribosomal RNA genes.  
Manning, J., et al.  
Biochemistry (1977) 16:1364-70
2. A one-step procedure for biotinylation and chemical crosslinking of lymphocyte surface and intracellular membrane-associated molecules  
Altin, J.G., et al.  
Anal Biochem (1995) 224:382-9
3. Immunoaffinity isolation of membrane antigens with biotinylated monoclonal antibodies and immobilized streptavidin matrices  
Updyke, T.V. and Nicolson, G.L.  
J Immunol Meth (1984) 73:83-95