

D-Biotin ethylenediamine

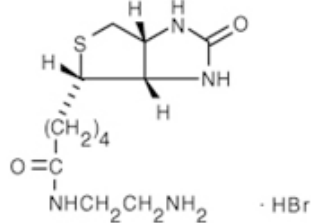
Catalog Number	Packaging Size
B007	25 mg

Storage upon receipt: -20°C

Introduction

The primary aliphatic amine of biotin ethylenediamine can be reversibly coupled to aldehydes and ketones to form a Schiff base – which can be reduced to a stable amine derivative by sodium borohydride (NaBH₄) or sodium cyanoborohydride (NaCNBH₃) to form new biotinylated probes. Carboxylic acids of proteins and other water-soluble biopolymers can be coupled to this molecule in aqueous solution using water-soluble carbodiimides such as EDAC.

Specifications

Molecular Formula:	C ₁₂ H ₂₃ BrN ₄ O ₂ S	
Molecular Weight:	367.30	
CAS Number:	216299-38-6	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Biotinylation reagent

References:

1. [Biotinylation reagents for the study of cell surface proteins.](#)
Elia G,
Proteomics (2008) 8:4012-4024
2. [Using laser scanning confocal microscopy as a guide for electron microscopic study: a simple method for correlation of light and electron microscopy.](#)
Sun XJ, Tolbert LP, Hildebrand JG
J Histochem Cytochem (1995) 43:329-335
3. Intracellular Dye-Filling *in Vivo* Combined with Light and Electron Microscopic Immunocytochemistry.
Pilowsky PM, Llewellyn-Smith IJ
Neuroscience Protocols, Wouterlood FG, Ed. 1993; (93-050-20):na pp. 01-16