

D-Biotin-XX, SE

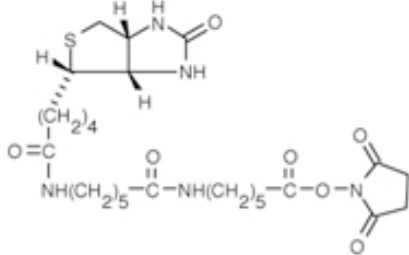
Catalog Number	Packaging Size
B005	50 mg

Storage upon receipt: -20°C

Introduction

The amine-reactive **biotin-XX, SE** can be used to attach this important hapten to biomolecules of interest for subsequent detection with streptavidin, avidin or NeutrAvidin® biotin-binding protein. This molecule has a 14 atom spacer designated by "XX" to facilitate reactivity with the molecule and eventual accessibility of the biotin to avidin or streptavidin.

Specifications

Molecular Formula:	C ₂₆ H ₄₁ N ₅ O ₇ S	
Molecular Weight:	567.70	
CAS Number:	89889-52-1	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Biotinylation reagent

References:

1. [Plasticity of B cell receptor internalization upon conditional depletion of clathrin.](#)
Stoddart A, Jackson AP, Brodsky FM
Mol Biol Cell (2005) 16:2339-2348
2. [Acetylcholinesterase dynamics at the neuromuscular junction of live animals.](#)
Krejci E, Martinez-Pena y Valenzuela I, Ameziane R, Akaaboune M
J Biol Chem (2006) 281:10347-10354
3. [Custom fluorescent-nucleotide synthesis as an alternative method for nucleic acid labeling.](#)
Henegariu O, Bray-Ward P, Ward DC
Nat Biotechnol (2000) 18:345-348