

D-Biotin, SE

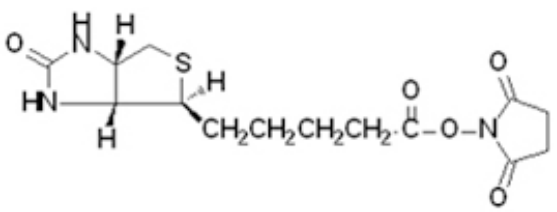
Catalog Number	Packaging Size
B003	100 mg

Storage upon receipt: -20°C

Introduction

The amine-reactive **D-biotin succinimidyl ester** can be used to attach this important hapten to biomolecules of interest for subsequent detection with streptavidin, avidin or NeutrAvidin® biotin-binding protein.

Specifications

Molecular Formula:	C ₁₄ H ₁₉ N ₃ O ₅ S	
Molecular Weight:	341.38	
CAS Number:	35013-72-0	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Biotinylation reagent

References:

1. [Zymogen/enzyme discrimination using peptide chloromethyl ketones.](#)
Williams EB, Krishnaswamy S, Mann KG
J Biol Chem (1989) 264:7536-7545
2. [Clathrin-mediated endocytosis and recycling of autocrine motility factor receptor to fibronectin fibrils is a limiting factor for NIH-3T3 cell motility.](#)
Le PU, Benlimame N, Lagana A, Raz A, Nabi IR
J Cell Sci (2000) 113 (Pt 18):3227-3240
3. Bioluminescent Homogeneous Enzyme Binding Assay for Biotin Using Luciferase as a Label.
Arakawa H, Maeda M, Tsuji A
Anal Lett (1992) 25:1055-1055