

D-Biocytn hydrazide

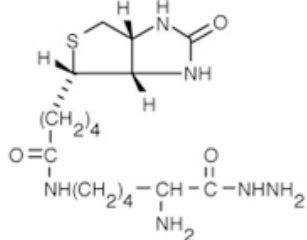
Catalog Number	Packaging Size
B012	25 mg

Storage upon receipt: -20°C

Introduction

Biocytn hydrazide can be used as microinjectable anterograde and transneuronal tracers that is fixable with aldehyde-based fixatives. Additionally, the primary aliphatic amine of this molecule 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 ₁₆ H ₃₀ N ₆ O ₃ S	
Molecular Weight:	386.51	
CAS Number:	102743-85-1	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Anterograde and transneuronal tracers

References:

1. 5-HT_{4(a)} receptors avert opioid-induced breathing depression without loss of analgesia.
Manzke T, Guenther U, Ponimaskin EG, Haller M, Dutschmann M, Schwarzacher S, Richter DW
Science (2003) 301:226-229
2. [Enzyme-modulated cleavage of dsDNA for supramolecular design of biosensors.](#)
Yan F, Sadik OA
Anal Chem (2001) 73:5272-5280
3. [Capture of peptides with N-terminal serine and threonine: a sequence-specific chemical method for Peptide mixture simplification.](#)
Chelius D, Shaler TA
Bioconjug Chem (2003) 14:205-211