

D-Biocytn

Catalog Number	Packaging Size
B010	100 mg

Storage upon receipt: -20°C

Introduction

Biocytn can be used as microinjectable anterograde and transneuronal tracers that is fixable with aldehyde-based fixatives. Additionally the primary amine on the lysine can be reversibly coupled to aldehydes and ketones to form a Schiff base – which can be reduced to a generate stable amine derivative by sodium borohydride (NaBH₄) or sodium cyanoborohydride (NaCNBH₃). 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 ₂₈ N ₄ O ₄ S	
Molecular Weight:	372.48	
CAS Number:	576-19-2	
Storage Conditions:	-20°C	
Shipping Condition:	Room Temperature	

Applications

Anterograde and transneuronal tracers

References:

1. [Spike resonance properties in hippocampal O-LM cells are dependent on refractory dynamics.](#)
Kispersky TJ, Fernandez FR, Economo MN, White JA,
J Neurosci (2012) 32:3637-3651
2. [Light exposure induces short- and long-term changes in the excitability of retinorecipient neurons in suprachiasmatic nucleus.](#)
LeSauter J, Silver R, Cloues R, Witkovsky P,
J Neurophysiol (2011) 106:576-588
3. [Differential GABAB-receptor-mediated effects in perisomatic- and dendrite-targeting parvalbumin interneurons.](#)
Booker SA, Gross A, Althof D, Shigemoto R, Bettler B, Frotscher M, Hearing M, Wickman K,
Watanabe M, Kulik Á, Vida I,
J Neurosci (2013) 33:7961-7974