

Biocytin hydrazide *CAS 102743-85-1*

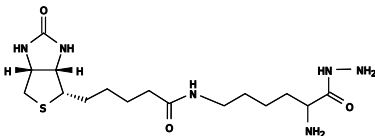
Catalog Number: 3086

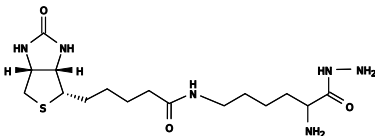
Unit Size: 25 mg

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid white
Molecular Weight	386.51
Soluble In	Water
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

This water-soluble biocytin hydrazide is used for the selective non-radioactive detection of glycoconjugates. The method involves either chemical (periodate-induced) or enzymatic (via galactose oxidase) oxidation of glycoconjugates, the resultant aldehyde groups are then labeled with biocytin hydrazide, followed by interaction with an avidin-based enzyme probe.