

Product Information Sheet

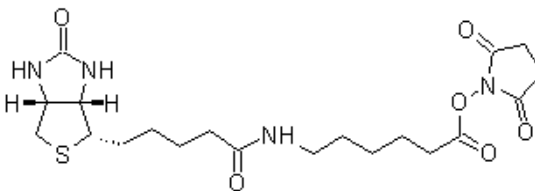
Ordering Information

Product Number: 3010
Product Name: Biotin-X, succinimidyl ester *CAS# 72040-63-2*
Unit Size: 25 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 454.54

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: N/A
Emission Wavelength: N/A

Application Notes

This biotin has a spacer arm of ~22 Å that reduces steric hindrance when binding several biotinylated molecules to one avidin complex; This reagent penetrates cell membranes because there is no charged group.