

Product Information Sheet

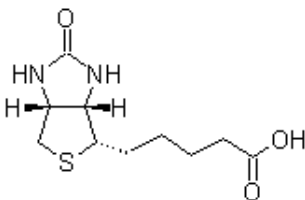
Ordering Information

Product Number: 3001
Product Name: Biotin *CAS# 58-85-5*
Unit Size: 1 g
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 244.31

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: N/A

Emission Wavelength: N/A

Application Notes

Biotin binds with high affinity to both avidins and streptavidins. Biotin and its derivatives can be conjugated to many biomolecules without significantly altering the biological activity of the target molecules since biotin is relatively a small molecule. A biopolymer (such as proteins) can react with several molecules of biotin that, in turn, can each bind one avidin. This characteristic greatly increases the sensitivity of many biological assays. Biotin derivatives are widely used for biological detections and purification. Our biotin is highly purified and tested for avidin-binding affinity.