www.aatbio.com; E-Mail: info@aatbio.com

Product Information Sheet

Ordering Information

Product Number: 3004

Product Name: Biotin cadaverine

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: 442.50

Chemical Structure:

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

Application Notes

Biotin and biotin derivatives bind 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. This biotin derivative is carbonyl-reactive. It is used for colorimetric assays for Factor XII (when used with EDC) and cellular transglutaminase. It is also widely used for labeling peptides (carboxylic acid groups) and nucleotides (5´ phosphate groups) via the use of EDC.