

APC-Biotin Conjugate

Catalog Number: 3091

Unit Size: 1 mg

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	651 nm
Emission Wavelength	660 nm

Applications

The bifunctional APC-Biotin conjugate demonstrates excellent avidin binding and strong fluorescence upon binding to a protein or nucleic acid target. It has avidin-binding properties similar to native biotin in terms of high affinity, fast association, and non-cooperative binding to avidin and streptavidin tetramers. APC-Biotin conjugate can be used for detecting and quantifying biotin binding sites of avidin, streptavidin or neutravidin. APC (allophycocyanin), isolated from *Spirulina* sp. blue-green algae, might be one of the brightest NIR fluorophores. The APC-Biotin conjugate is prepared from crosslinked APC, thus has higher stability.