

Cy5 biotin conjugate

Catalog number: 3100

Unit size: 5 mg

Product Details

Storage Conditions Freeze (<-15 °C)

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Blue solid

Molecular Weight 1353.82

Soluble In Water

Chemical Structure

Spectral Properties

Excitation Wavelength 651 nm

Emission Wavelength 670 nm

Applications

This bifunctional Cy5 biotin conjugate demonstrates excellent avidin binding and strong fluorescence. 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. Cy5 Biotin can be used for detecting and quantifying biotin binding sites of avidin, streptavidin or neutravidin. This reagent overcomes major shortcomings of commonly used Biotin-4-fluorescein, e.g., poor solubility in aqueous media and pH dependent absorbance and emission. Cy5 Biotin is a water-soluble reagent and its red fluorescence is pH independent from pH 4 to pH 10. A flexible PEG spacer between biotin moiety and fluorescent tag minimize steric hindrance involved in binding to avidin, streptavidin or neutravidin.