

Cy7 biotin conjugateCatalog number: 3106
Unit size: 5 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Soluble In	Water

Spectral Properties

Excitation Wavelength	756 nm
Emission Wavelength	779 nm

Applications

Cy7 biotin conjugate demonstrates excellent avidin binding and strong NIR 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. Cy7 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 short wavelength emission. Cy7 Biotin is a water-soluble reagent and its near infrared fluorescence is pH independent from pH 4 to pH 10. A flexible PEG spacer between biotin moiety and fluorescent tag minimizes the steric hindrance involved in binding to avidin, streptavidin or neutravidin.