

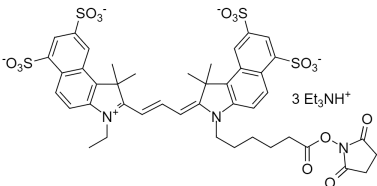
Cyanine 3.5 monosuccinimidyl ester [equivalent to Cy3.5® NHS ester]

Catalog number: 148
Unit size: 1 mg

Product Details

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

Chemical Properties

Appearance	Purple solid
Molecular Weight	1291.66
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	579 nm
Emission Wavelength	591 nm

Applications

A variety of cyanine 3.5 (Cy3.5®) dyes has been used to label biological molecules for fluorescence imaging and other fluorescence-based biochemical analysis. They are widely used for labeling peptides, proteins and oligos etc. Cy3.5® dyes have enhanced fluorescence upon binding to proteins. Cy3.5® NHS ester readily reacts with amino groups. Cy3® NHS ester readily reacts with amino groups. AAT Bioquest offers Cy dye NHS esters in the form of triethylammonium salts that are more soluble in DMSO and DMF than the corresponding potassium salts that are offered by some other vendors. The Cy dye triethylammonium salts have the same reactivity and give the conjugates identical to the the Cy dye potassium salts. Cy3.5® is the trademark of GE Healthcare.