

Cyanine 5B monosuccinimidyl ester

Catalog Number: 925

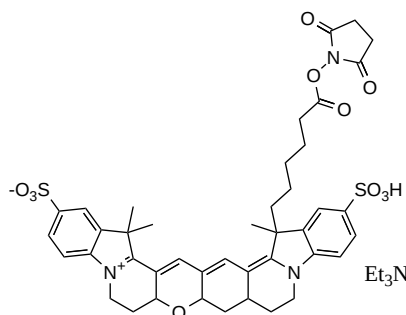
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	671 nm
Emission Wavelength	682 nm

Applications

Cyanine 5B monosuccinimidyl ester is a cyanine dye used to label biological molecules for fluorescence imaging and other fluorescence-based biochemical analysis. Cyanine 5B is a modified Cyanine 5 dye that offers improved photostability compared to conventional Cy5 fluorophores. It is commonly used for labeling peptides, proteins, and other amine-containing biomolecules through NHS ester chemistry.

Cyanine 5B monosuccinimidyl ester exhibits excitation and emission maxima of approximately 671 nm and 682 nm, respectively, making it compatible with standard red laser excitation systems. The NHS ester functional group reacts with primary amines to form stable covalent conjugates. This dye is suitable for applications including fluorescence imaging, flow cytometry, and other fluorescence-based detection methods.