

Cyanine 5B monoacid

Catalog Number: 924

Unit Size: 5 mg

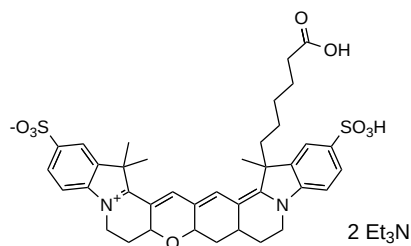
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	937.27
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	671 nm
Emission Wavelength	682 nm

Applications

Cyanine 5B monoacid is a cyanine dye used to label biological molecules for fluorescence imaging and other fluorescence-based biochemical analysis. It is commonly used for labeling peptides, proteins, and oligonucleotides. Cyanine 5B acid is a modified Cyanine 5 dye that offers improved photostability compared to conventional Cy5 fluorophores. It contains a carboxylic acid functional group and serves as a precursor for preparing reactive derivatives for bioconjugation. Cyanine 5B acid is suitable for applications including fluorescence imaging and other fluorescence-based detection methods.