

Cyanine 5B maleimide

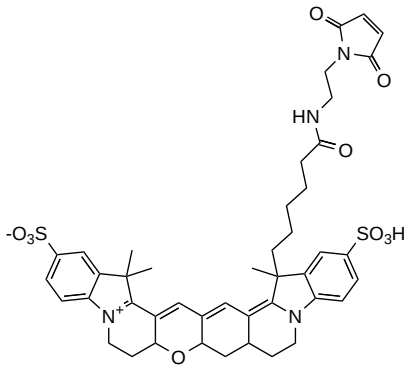
Catalog Number: 926

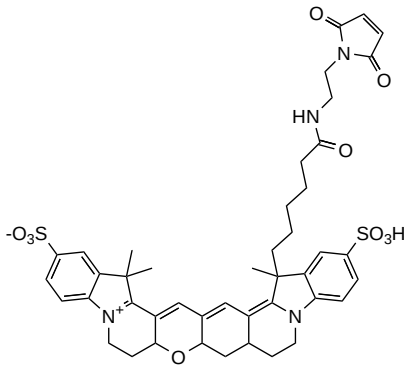
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	857
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	671 nm
Emission Wavelength	682 nm

Applications

Cyanine 5B maleimide 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 proteins and peptides containing accessible thiol groups.

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