

Cy3B PEG12 NHS ester

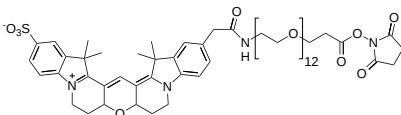
Catalog Number: 939

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	1216.48
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	560 nm
Emission Wavelength	571 nm

Applications

Cy3B PEG12 NHS ester is a bright, amine-reactive fluorescent labeling reagent that combines the superior photostability of the Cy3B fluorophore with the enhanced solubility and reduced steric hindrance of a PEG12 spacer. This compound features an N-hydroxysuccinimide (NHS) ester functional group, which reacts efficiently with primary amines on proteins, peptides, and other biomolecules under mild conditions. The inclusion of a PEG12 linker improves hydrophilicity and minimizes nonspecific interactions, making it especially suitable for applications that require high labeling specificity and minimal background signal. Cy3B is known for its high quantum yield and resistance to photobleaching, offering significant advantages over traditional Cy3 dyes in demanding applications such as fluorescence microscopy, flow cytometry, and single-molecule detection. With excitation and emission maxima around 560 nm and 571 nm, respectively, Cy3B PEG12 NHS ester is compatible with standard Cy3 filter sets and serves as a reliable choice for the development of robust, high-performance fluorescent probes.