

Cy3B PEG8 methyltetrazine

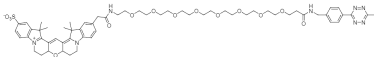
Catalog number: 947

Unit size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	1167.39
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	560 nm
Emission Wavelength	571 nm

Applications

Cy3® dye is one of the most common cyanine dyes used for labeling proteins, nucleic acids and other biological molecules. A variety of Cy3® 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. However, Cy3 suffers low fluorescence intensity. It is the least fluorescent dyes among all the Cy dyes. Cy3B is an improved version of Cy3 dyes with significantly increased fluorescence quantum yield and photostability. Cy3B PEG8 methyltetrazine can be used for conjugations with TCO-containing biological molecules under copper-free conditions. Compared to the other click reactions, the TCO-tetrazine reaction gives much higher yields under mild reaction conditions. Cy3B has poor water solubility. This Cy3B PEG8 methyltetrazine contains a PEG8 linker that significantly improves its water solubility. Cy3® is the trademark of GE Healthcare.