

Cy3B Tetrazine

Catalog number: 948

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 906.07

Soluble In DMSO

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 tetrazine 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. Cy3® is the trademark of GE Healthcare.