

## СуЗВ ТСО

Catalog number: 946 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	1093.34
Soluble In	DMSO
Spectral Properties	
Excitation Wavelength	560 nm
Emission Wavelength	571 nm

## Applications

Cy3<sup>®</sup> dye is one of the most common cyanine dyes used for labeling proteins, nucleic acids and other biological molecules. A variety of Cy3<sup>®</sup> 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 TCO can be used for conjugations with tetrazine-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<sup>®</sup> is the trademark of GE Healthcare.