

Sulfo-Cyanine 3 carboxylic acid

Catalog number: 260, 261 Unit size: 25 mg, 100 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Solid
Molecular Weight	638.73
Soluble In	DMSO
Chemical Structure	
Spectral Properties	
Excitation Wavelength	555 nm
Emission Wavelength	569 nm

Applications

A variety of cyanine 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. Cy3[®] dyes are one type of the most common red fluorophores. These versatile fluorophores can tolerate a pH range of 3-10 for use in a variety of applications at biologically relevant pHs. The dyes are also DMSO tolerant and photostable to enable transfer from storage to assay without loss of performance. The aqueous solubility eliminates the need for organic solvents in the assay buffers. Sulfo-Cyanine 3 is an alternative dye primarily used for labeling peptides and oligos in replacing Cy3 that was originally developed by GE Healthcare. Cy3[®] is the trademark of GE Healthcare.