

6-TRITC [Tetramethylrhodamine-6-isothiocyanate] *CAS 80724-20-5*

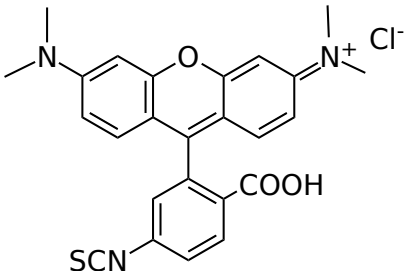
Catalog Number: 417, 418

Unit Size: 1 mg, 5 mg

Product Details

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

Chemical Properties

Appearance	Solid purple
Molecular Weight	479.98
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	544 nm
Emission Wavelength	570 nm

Applications

6-TRITC (also called R isomer) is the other isomer of the TRITC labeling reagent that is widely used in preparing bioconjugates of proteins and nucleic acids. Complimentary to 5-TRITC, the 6-isomer is predominantly used in labeling nucleotides and nucleic acids. Cautions must be exercised for the storage of TRITC conjugates.