

5(6)-TRITC [Tetramethylrhodamine-5-(and-6)-isothiocyanate] *CAS 95197-95-8*

Catalog Number: 409, 410

Unit Size: 25 mg, 5 mg

Product Details

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

Chemical Properties

Appearance	Solid dark purple
Molecular Weight	479.98
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	544 nm
Emission Wavelength	570 nm

Applications

5(6)-TRITC is an amino-reactive labeling reagent that is widely used in preparing bioconjugates of proteins and nucleic acids. The resultant conjugates have similar spectral properties to those prepared from 5(6)-TAMRA, SE. However, the latter conjugates are much more stable. Cautions must be exercised for the storage of TRITC conjugates.