

## 5-TRITC [Tetramethylrhodamine-5-isothiocyanate] \*CAS 80724-19-2\*

Catalog Number: 415, 416

Unit Size: 1 mg, 5 mg

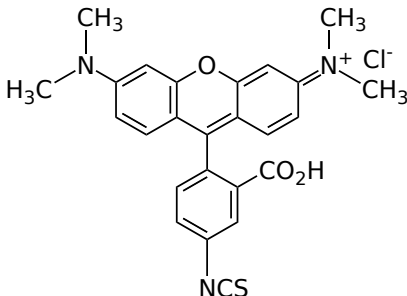
### Product Details

---

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

### Chemical Properties

---

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

### Spectral Properties

---

Excitation Wavelength	544 nm
Emission Wavelength	570 nm

### Applications

---

5-TRITC (also called G isomer) is a purified single isomer of the 5(6)-TRITC mixed isomers. This labeling reagent is predominantly used in labeling peptides and proteins. Cautions must be exercised for the storage of TRITC conjugates.