TRITC, 5 isomer

5-TRITC (Tetramethylrhodamine 5-isothiocyanate chloride) is spectrally similar to TAMRA (90022).



Product attributes

 Excitation/Emission
 544/572 nm

 CAS number
 80724-19-2

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

5-TRITC (Tetramethylrhodamine 5-isothiocyanate chloride) is spectrally similar to TAMRA (90022).

- λ_{Ex}/λ_{Em}: 544/572 nm (MeOH)
- Red solid soluble in DMF
- MW: 479.98
- [80724-19-2]

$$(H_3C)_2N$$
 O
 $N(CH_3)_2$
 CI
 CO_2H
 $N=C=S$

This datasheet was generated on November 17, 2025 at 09:49:46 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/tetramethylrhodamine-5-isothiocyanate-chloride-5-tritc-gisomer/