

## TRITC, 5 isomer

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



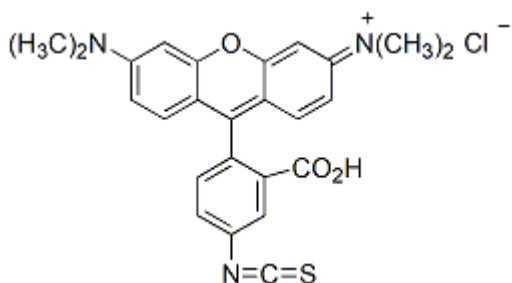
### Product attributes

Excitation/Emission	544/572 nm
CAS number	80724-19-2

## Product Description

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

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



This datasheet was generated on January 16, 2026 at 07:53:13 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/tetramethylrhodamine-5-isothiocyanate-chloride-5-tritc-gisomer/>