

TRITC, 6 isomer

6-TRITC (Tetramethylrhodamine 6-isothiocyanate chloride) is spectrally similar to TAMRA ([90022](#)).



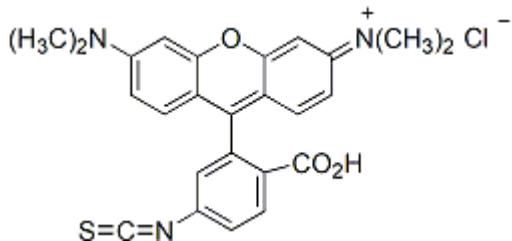
Product attributes

Excitation/Emission	544/572 nm
---------------------	------------

Product Description

6-TRITC (Tetramethylrhodamine 6-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



This datasheet was generated on January 21, 2026 at 04:28:49 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tetramethylrhodamine-6-isothiocyanate-chloride-6-tritc-risomer/>