

TRITC, 6 isomer

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



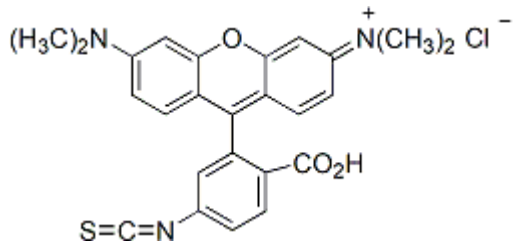
Product attributes

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

Product Description

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

- $\lambda_{Ex}/\lambda_{Em}$: 544/572 nm (MeOH)
- Red solid soluble in DMF
- MW: 479.98



This datasheet was generated on June 16, 2026 at 09:00:37 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/tetramethylrhodamine-6-isothiocyanate-chloride-6-tritc-isomer/>