

TRITC, 5 and 6 isomers

This item is discontinued and no longer available. We continue to offer TRITC, 5 isomer ([90018](#)) and TRITC, 6-isomer ([90019](#)).



Product attributes

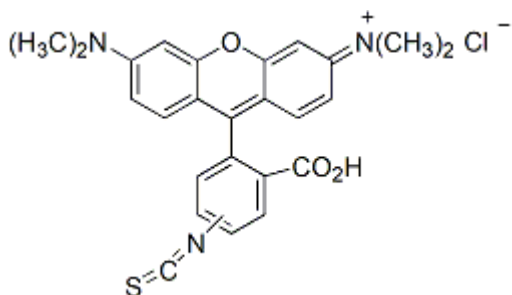
Excitation/Emission	544/572 nm
---------------------	------------

Product Description

This item is discontinued and no longer available. We continue to offer TRITC, 5 isomer ([90018](#)) and TRITC, 6-isomer ([90019](#)).

5(6)-TRITC (Tetramethylrhodamine isothiocyanate chloride) is an amine-reactive rhodamine derivative that 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 02:58:24 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tetramethylrhodamine-isothiocyanate-chloride-56-tritcmixed-isomers/>