

MTS-6-TAMRA

MTS-6-TAMRA (6-carboxytetramethylrhodamine methanethiosulfonate) is a fast-reacting and highly selective thiol-reactive label that permits one to attach the popular fluorescent dye TAMRA to a thiol group.



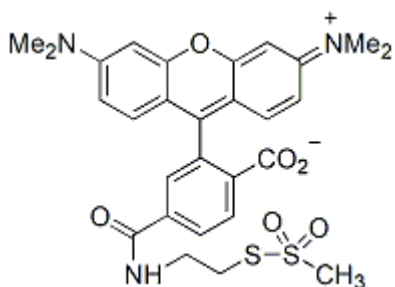
Product attributes

Excitation/Emission	540/565 nm
---------------------	------------

Product Description

MTS-6-TAMRA (6-carboxytetramethylrhodamine methanethiosulfonate) is a fast-reacting and highly selective thiol-reactive label that permits one to attach the popular fluorescent dye TAMRA to a thiol group.

- $\lambda_{Ex}/\lambda_{Em}$: 540/565 nm (MeOH)
- Red solid soluble in DMF or DMSO
- MW: 567



This datasheet was generated on December 18, 2025 at 04:55:39 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mts-6-tetramethylrhodamine-mts-6-tamra/>