

MTS-TAMRA

MTS-TAMRA (MTS-5(6)-carboxytetramethylrhodamine, mixed isomers) is a fast-reacting and highly selective thiol-reactive label that permits one to attach the popular fluorescent dye carboxytetramethylrhodamine to a thiol group.



Product attributes

Call us: 800-304-5357

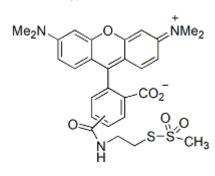
Excitation/Emission

540/565 nm

Product Description

MTS-TAMRA (MTS-5(6)-carboxytetramethylrhodamine, mixed isomers) is a fast-reacting and highly selective thiol-reactive label that permits one to attach the popular fluorescent dye carboxytetramethylrhodamine to a thiol group.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 540/565 nm
- Red solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C₂₈H₂₉N₃O₆S₂
- MW: 567



References

1. J Biol Chem 281(39), 28837 (2006).

This datasheet was generated on November 26, 2025 at 04:04:54 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mts-tamra-mixed-isomers-mts-56-carboxytetramethylrhodamine/