

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.



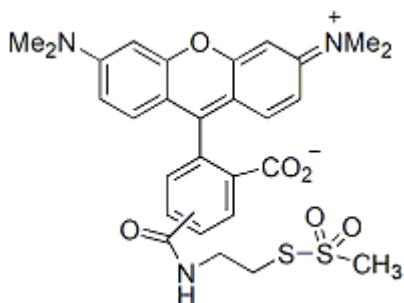
Product attributes

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
- $C_{28}H_{29}N_3O_6S_2$
- MW: 567



References

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

This datasheet was generated on June 3, 2026 at 11:34:23 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/mts-tamra-mixed-isomers-mts-56-carboxytetramethylrhodamine/>