

MTS-5-TAMRA

MTS-5-TAMRA (5-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-5-TAMRA (5-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

