

## MTS-ROX , mixed isomers

MTS-ROX (MTS-5(6)-carboxy-X-rhodamine) is a fast-reacting and highly selective thiol-reactive rhodamine dye.



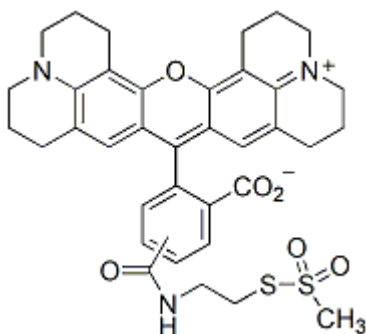
### Product attributes

Excitation/Emission	568/595 nm
---------------------	------------

## Product Description

MTS-ROX (MTS-5(6)-carboxy-X-rhodamine) is a fast-reacting and highly selective thiol-reactive rhodamine dye.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 568/595 nm
- Dark Red solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- $C_{36}H_{37}N_3O_6S_2$
- MW: 671.8



This datasheet was generated on June 22, 2026 at 12:22:17 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mts-rox-mixed-isomers-mts-56-carboxy-x-rhodamine/>