

MTS-Rhodamine 6G

MTS-Rhodamine 6G is a fast-reacting and highly selective thiol-reactive rhodamine dye.



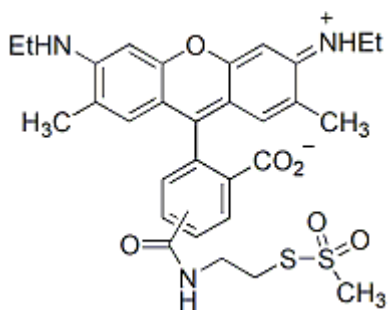
Product attributes

Excitation/Emission	520/546 nm
---------------------	------------

Product Description

MTS-Rhodamine 6G is a fast-reacting and highly selective thiol-reactive rhodamine dye.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 520/546 nm
- Orange red solid soluble in DMF or DMSO
- Store at -20 °C and protect from light
- $C_{30}H_{33}N_3O_6S_2$
- MW: 595



This datasheet was generated on December 16, 2025 at 09:09:58 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mts-rhodamine-6g/>