

MTS-CR110

MTS-CR110 (MTS-5-(and-6)-carboxyrhodamine 110, mixed isomers) is a green fluorescent thiol-reactive dye. CR110 (5-(and-6)-carboxyrhodamine 110) has absorption and emission spectra (502/524 nm) similar to those of fluorescein (495/519 nm) but its photostability is far better than that of the latter.



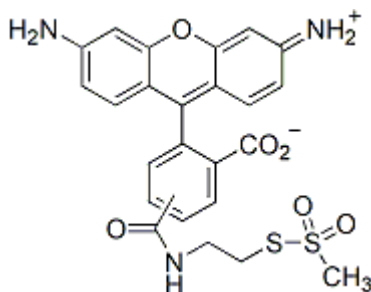
Product attributes

Excitation/Emission	502/524 nm
---------------------	------------

Product Description

MTS-CR110 (MTS-5-(and-6)-carboxyrhodamine 110, mixed isomers) is a green fluorescent thiol-reactive dye. CR110 (5-(and-6)-carboxyrhodamine 110) has absorption and emission spectra (502/524 nm) similar to those of fluorescein (495/519 nm) but its photostability is far better than that of the latter. Thus, MTS-CR110 is a superior green fluorescent dye for labeling thiols.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 502/524 nm
- Orange solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- $C_{24}H_{21}N_3O_8S_2$
- MW: 511



This datasheet was generated on June 14, 2026 at 01:26:32 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/mts-cr110-mts-5-and-6-carboxyrhodamine-110-mixed-isomers/>