

MTS-CR110-X

MTS-CR110-X (MTS-5-(and-6)-carboxyrhodamine 110-X, mixed isomers) is a green fluorescent thiol-reactive dye. It is similar to MTS-CR110 (catalog no. 91035) except that the former has a X spacer molecule between the dye and the MTS reactive group.



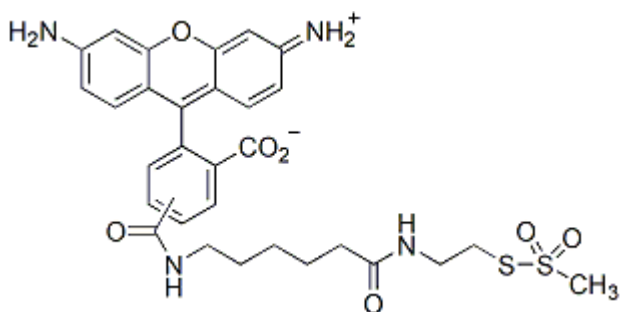
Product attributes

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

Product Description

MTS-CR110-X (MTS-5-(and-6)-carboxyrhodamine 110-X, mixed isomers) is a green fluorescent thiol-reactive dye. It is similar to MTS-CR110 (catalog no. [91035](#)) except that the former has a X spacer molecule between the dye and the MTS reactive group. The linker molecule may enhance fluorescence when the dye is attached to proteins.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 502/524 nm
- Orange solid soluble in DMF or DMSO
- Store at -20 °C and protect from light
- $C_{30}H_{32}N_4O_7S_2$
- MW: 624



This datasheet was generated on June 22, 2026 at 12:29:38 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/mts-cr110-x-mts-5-and-6-carboxyrhodamine-110-x-mixed-isomers/>