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.

siètum 8

Product attributes

Excitation/Emission

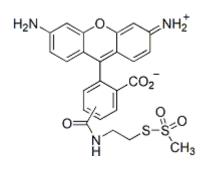
502/524 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

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
- C₂₄H₂₁N₃O₆S₂
- MW: 511



This datasheet was generated on November 4, 2025 at 04:28:19 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/