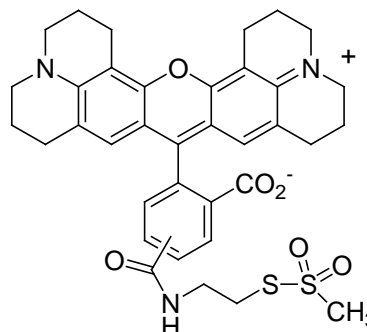


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: MTS-ROX, mixed isomers (MTS-5(6)-carboxy-X-rhodamine)**CATALOG #** 91037**MOLECULAR INFORMATION:** C₃₆H₃₇N₃O₆S₂
MW: 671.8**PROPERTIES:**

Color & Form	Dark red solid
Purity	≥ 95% by TLC
Solubility	Soluble in DMF or DMSO
Absorption/Emission	568/595 nm (MeOH)
Extinction Coefficient	≥55,000

STORAGE AND HANDLING: Store at -20 °C and protect from light.**APPLICATION:** MTS-ROX is a fast-reacting and highly selective thiol-reactive rhodamine dye.**TOXICITY:** Unknown

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--