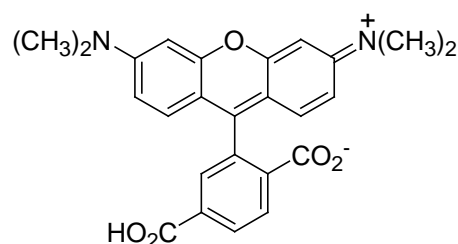


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 6-Carboxytetramethylrhodamine, single isomer (6-TAMRA)**CATALOG #:** 90008**MOLECULAR INFORMATION:** Mwt: 430.5  
C<sub>25</sub>H<sub>22</sub>N<sub>2</sub>O<sub>5</sub>**PROPERTIES:**

**Color & Form** Dark red powder  
**Isomer purity**  $\geq 98\%$   
**Solubility** Soluble in DMSO, DMF, MeOH, or H<sub>2</sub>O (pH > 6)  
**Absorption/Emission** 540/565 nm (MeOH)

**STORAGE AND HANDLING:**Store at  $\leq 4$  °C and protect from light, especially in solution.**APPLICATION:**

6-Carboxytetramethylrhodamine, triethylammonium salt is a single isomer rhodamine dye. The 6-carboxy group can be activated to succinimidyl ester or mixed anhydride for coupling with amines using standard methods.

**TOXICITY:**

Not established.

**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:** 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.