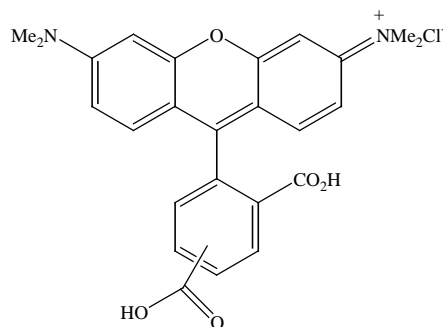


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 5-(and-6)-carboxytetramethylrhodamine, mixed isomers (5(6)-TAMRA)

**CATALOG #** 90027

**MOLECULAR INFORMATION:** C<sub>25</sub>H<sub>23</sub>ClN<sub>2</sub>O<sub>5</sub>  
MWt: 466.92

**PROPERTIES:**

<b>Color &amp; Form</b>	Dark red solid
<b>Purity</b>	97% by HPLC
<b>Solubility</b>	Soluble in DMF, DMSO or MeOH
<b>Absorption/Emission</b>	540nm/565nm (MeOH)
<b>Extinction Coefficient</b>	95,000

**STORAGE AND HANDLING:** Store desiccated at ≤ 4 °C. Protect from light, especially when in solution. When stored as recommended, the product is stable for at least one year from the date it is received.

**APPLICATION:** 5-(and-6)-carboxytetramethylrhodamine is a widely used red fluorescent dye, which can be optimally excited by the 546 Hg line. Biotium offers carboxytetramethylrhodamine as 5- and 6- mixed isomers. The dye can be coupled to primary or secondary amines via standard peptide chemistry.

**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> <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>
--