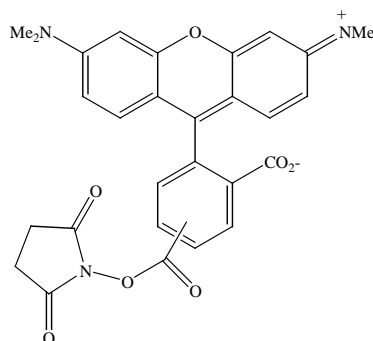


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** **5-(and-6)-carboxytetramethylrhodamine succinimidyl ester (5(6)-TAMRA SE)**

**CATALOG #** **90022**

**MOLECULAR INFORMATION:** C<sub>29</sub>H<sub>25</sub>N<sub>3</sub>O<sub>7</sub>  
MWt: 528  
[246256-50-8]

**PROPERTIES:**

<b>Color &amp; Form</b>	Red solid
<b>Purity</b>	≥ 98% by TLC
<b>Solubility</b>	Soluble in DMF, DMSO
<b>Absorption/Emission</b>	545nm/575nm (MeOH)
<b>Extinction Coefficient</b>	60,000

**STORAGE AND HANDLING:** Store desiccated at -20 °C. Protect from light, especially when in solution.

**APPLICATION:** Carboxytetramethylrhodamine is one of the most commonly used red fluorescent dye, which can be optimally excited by 546 Hg line. The succinimidyl ester reacts readily with primary or secondary amines under mild conditions. Biotium offers a very high grade TAMRA SE.

**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>
--