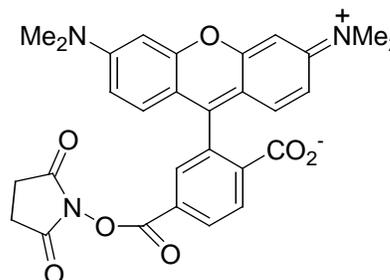


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 6-Carboxytetramethylrhodamine succinimidyl ester, 90mg/mL (0.17M) in anhydrous DMSO

**CATALOG #** 90097

**MOLECULAR INFORMATION:** C<sub>29</sub>H<sub>25</sub>N<sub>3</sub>O<sub>7</sub>  
Mwt:527.5

**PROPERTIES:**

**Color & Form** Dark red liquid  
**Purity** ≥ 98% by TLC  
**Solubility** Soluble in DMSO or DMF  
**Absorption/Emission** 540nm/565nm (MeOH)  
**Extinction Coefficient** 90,000

**STORAGE AND HANDLING:** The solution is good for at least one month if stored at -20°C and keep away from moisture.

**APPLICATION:** Amine-reactive form of 6-carboxytetramethylrhodamine single isomer.

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

**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.