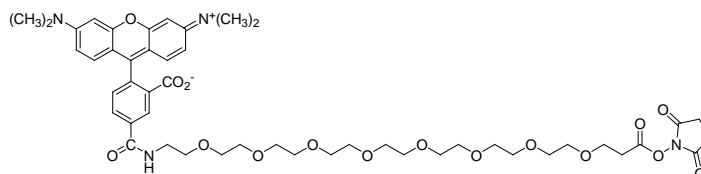


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Carboxytetramethylrhodamine-PEO₈-propionate succinimidyl ester
(5-TAMRA-PEO₈ SE)

CATALOG # 90087

MOLECULAR INFORMATION: C₄₈H₆₂N₄O₁₆
MWt: 951.02

**PROPERTIES:**

Color & Form Dark red solid
Purity ≥ 90% by TLC (remaining amount is free acid)
Solubility Soluble in DMSO or DMF
Absorption/Emission 540/565 nm (MeOH)

STORAGE AND HANDLING: Store at -20 °C and protect from moisture and light

APPLICATION: 5-TAMRA-PEO₈ SE has a long water-soluble PEG spacer between tetramethylrhodamine and the amine-reactive SE group.

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.
