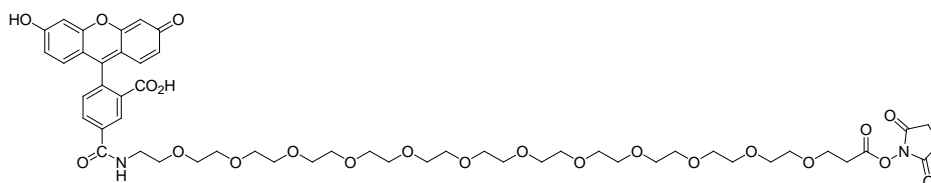


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-FAM-PEO₁₂ SE (5-carboxyfluorescein-PEO₁₂ propionate succinimidyl ester)**CATALOG #** 90093**MOLECULAR INFORMATION:** C₅₂H₆₈N₂O₂₂
MWt: 1073.1**PROPERTIES:**

Color & Form Yellow orange solid
Solubility Soluble in DMF, DMSO or pH 6.5 buffer
Absorption/Emission 494/521 nm
Extinction Coefficient 90,000

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

APPLICATION: 5-FAM-PEO₁₂ SE has an approximately 47 Angstrom water-soluble PEG spacer that separates fluorescein 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.