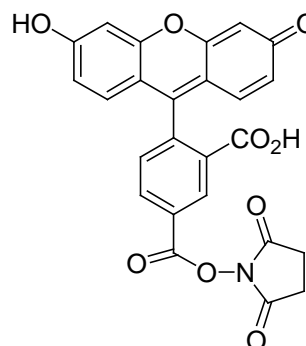


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Carboxyfluorescein, succinimidyl ester (5-FAM, SE)**CATALOG #** 90029**MOLECULAR INFORMATION:** C₂₅H₁₅NO₉
MWt: 473
[92557-80-7]**PROPERTIES:**

Color & Form	Yellow solid
Purity	99% by TLC
Solubility	Soluble in DMF, DMSO
Absorption/Emission	495nm/519nm (pH 9)
Extinction Coefficient	>70,000

STORAGE AND HANDLING: Store desiccated at -20 °C. Protect from light, especially when in solution.**APPLICATION:** 5-FAM SE is an amine-reactive form of 5-carboxyfluorescein single isomer.Ref: *Bioconjugate Chem.* 6, 447(1995)**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: <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>
--