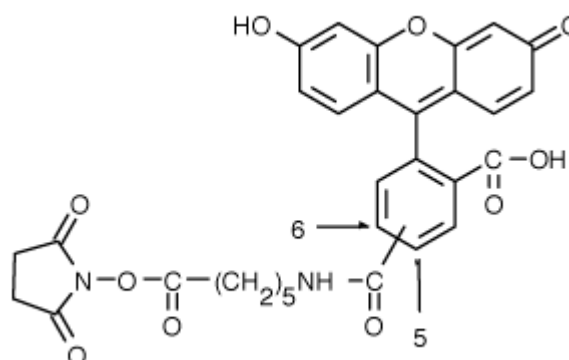


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 6-(fluorescein-5-(and-6)-carboxamido) hexanoic acid, succinimidyl ester
(5(6)-SFX), mixed isomers

CATALOG # 90044

MOLECULAR INFORMATION: C₃₁H₂₆N₂O₁₀
MWt: 586.55
[114116-31-8]

**PROPERTIES:**

Color & Form Orange solid
Purity 98% by TLC
Solubility Soluble in DMF, DMSO
Absorption/Emission 492nm\520nm
Extinction Coefficient 75,000

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

APPLICATION: 5(6)-SFX, mixed isomer is an amine-reactive fluorescein with a spacer group.

Ref.: 1) Jamieson GP, et al. *Cytometry* **14**, 32 (1993); 2) Prussin C., et al. *J Immunol Methods* **188**, 117 (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: 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.