

SFX succinimidyl ester, 5 and 6 isomers

5(6)-SFX has a 6-atom aminohexanoyl spacer group that separates the fluorescein dye and the amine-reactive group.



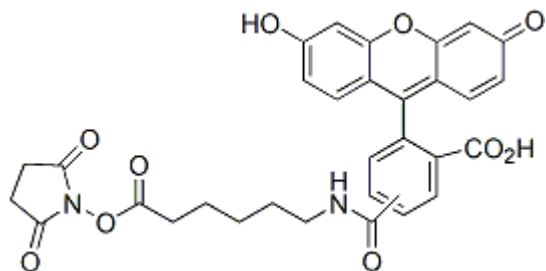
Product attributes

Excitation/Emission	494/521 nm (pH 9)
---------------------	-------------------

Product Description

5(6)-SFX (full name: 6-(Fluorescein-5-(and-6)-carboxamido)hexanoic acid,succinimidyl ester) has a 6-atom aminohexanoyl spacer group that separates the fluorescein dye and the amine-reactive group.

- $\lambda_{Ex}/\lambda_{Em}$ (pH 9) = 494/521 nm
- Yellow solid soluble in DMF, DMSO or pH>6.5 buffers
- Store at -20 °C and protect from light
- $C_{31}H_{26}N_2O_{10}$
- MW:586.55



This datasheet was generated on February 2, 2026 at 08:42:50 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/56-sfx-6-fluorescein-5-and-6-carboxamido-hexanoic-acid-succinimidyl-ester/>