

SFX (6-(Fluorescein-6-carboxamido)hexanoic acid) Succinimidyl Ester, 6 Isomer



Product attributes

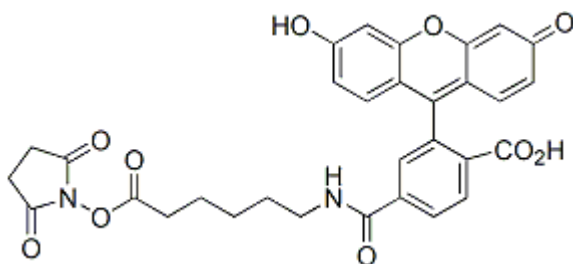
Excitation/Emission	494/521 nm (pH 9)
---------------------	-------------------

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

Product Description

6-SFX (full name: 6-(Fluorescein-6-carboxamido)hexanoic acid, succinimidyl ester, single isomer) 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
- C₃₁H₂₆N₂O₁₀
- MW:586.55



This datasheet was generated on June 22, 2026 at 03:30:37 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/6-sfx-6-fluorescein-6-carboxamidohexanoic-acid-succinimidyl-estersingle-isomer/>