

## FAM-PEO8 SE, 5 isomer

5-FAM-PEO<sub>8</sub> SE (full name: 5-Carboxyfluorescein-PEO<sub>8</sub> propionate succinimidyl ester) has an approximately 32 Angstrom water-soluble PEG spacer that separates fluorescein and the amine-reactive SE group.



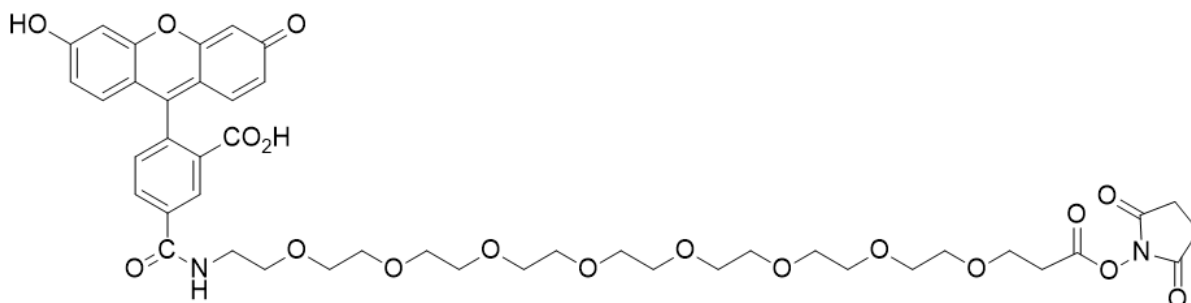
### Product attributes

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

## Product Description

5-FAM-PEO<sub>8</sub> SE (full name: 5-Carboxyfluorescein-PEO<sub>8</sub> propionate succinimidyl ester) has an approximately 32 Angstrom water-soluble PEG spacer that separates fluorescein and the amine-reactive SE group.

- $\lambda_{Ex}/\lambda_{Em}$  (pH 9) = 494/521 nm
- Yellow orange solid soluble in DMF, DMSO, or pH 6.5 buffer
- Store at -20 °C and protect from light
- C<sub>44</sub>H<sub>52</sub>N<sub>2</sub>O<sub>18</sub>
- MW: 896.89



This datasheet was generated on June 22, 2026 at 06:33:25 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-fam-peo8-se-5-carboxyfluorescein-peo8-propionate-succinimidyl-ester/>