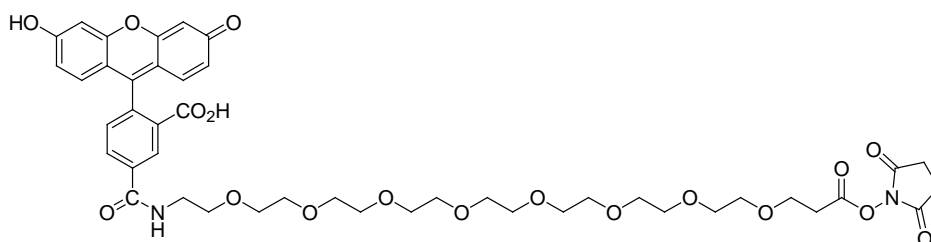


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 5-FAM-PEO<sub>8</sub> SE (5-carboxyfluorescein-PEO<sub>8</sub> propionate succinimidyl ester)**CATALOG #** 90085**MOLECULAR INFORMATION:** C<sub>44</sub>H<sub>52</sub>N<sub>2</sub>O<sub>18</sub>  
MWt: 896.89**PROPERTIES:**

**Color & Form** Yellow orange solid  
**Solubility** Soluble in DMF, DMSO or pH 6.5 buffer  
**Absorption/Emission** 494/521 nm  
**Extinction Coefficient** 94,000

**STORAGE AND HANDLING:** Store at -20 °C and protect from light**APPLICATION:** 5-FAM-PEO<sub>8</sub> SE has an approximately 32 Angstrom water-soluble PEG spacer that separates fluorescein and the amine-reactive SE group.

**RELATED PRODUCTS** 5-FAM-PEO<sub>12</sub> SE(#90093); 5-CR110-PEO<sub>4</sub> SE(#90081); 5-CR110-PEO<sub>8</sub> SE(#90084); 5-CR110-PEO<sub>12</sub> SE(#90086); 5-TAMRA-PEO<sub>8</sub> SE(#90087); 5-TAMRA-PEO<sub>12</sub> SE(#90088); 5-ROX-PEO<sub>8</sub> SE(#90089); 5-ROX-PEO<sub>12</sub> SE(#90090); Texas Red-PEO<sub>2</sub> SE(#90039); Texas Red-PEO<sub>8</sub> SE(#90099); Texas Red-PEO<sub>12</sub> SE(#90100); Sulforhodamine B-PEO<sub>2</sub> SE(#90043); Sulforhodamine B-PEO<sub>8</sub> SE(#90101); Sulforhodamine B-PEO<sub>12</sub> SE(#90102)

**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> 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.
---