

ROX-PEO8 SE, 5 isomer

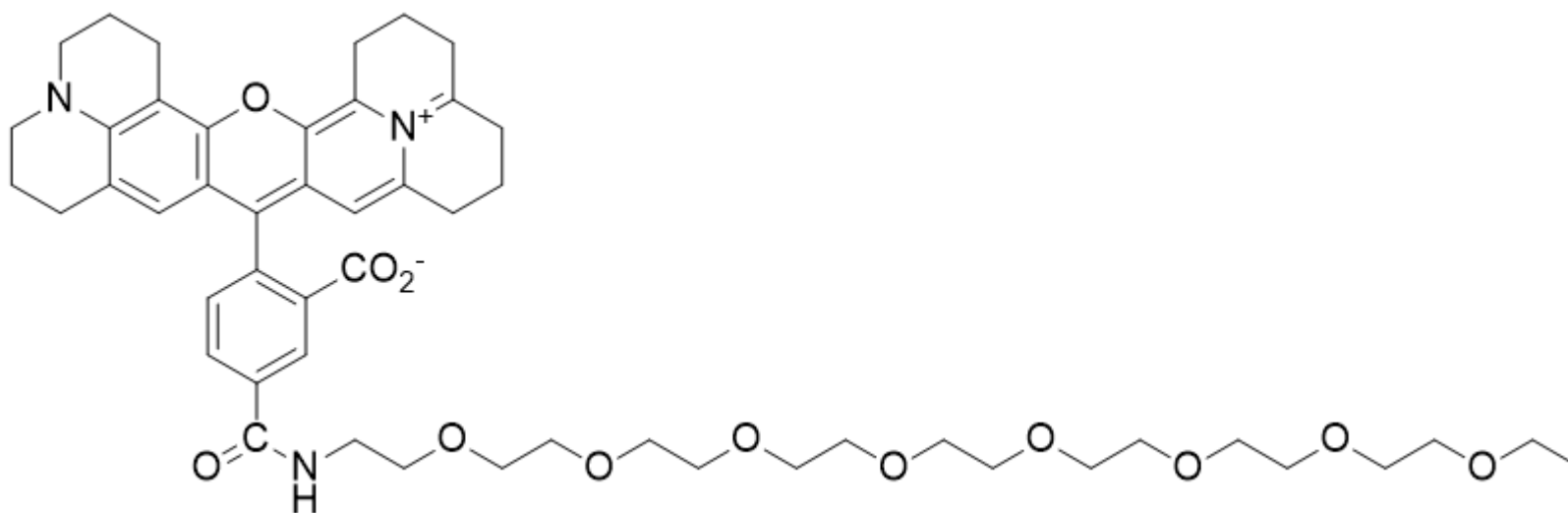
5-ROX-PEO₈ SE (full name: 5-Carboxy-X-rhodamine-PEO₈-propionate, succinimidyl ester) has an approximately 32 Angstrom water-soluble PEG spacer that separates the red fluorescent ROX dye and the amine-reactive SE group.



Product Description

5-ROX-PEO₈ SE (full name: 5-Carboxy-X-rhodamine-PEO₈-propionate, succinimidyl ester) has an approximately 32 Angstrom water-soluble PEG spacer that separates the red fluorescent dye and the amine-reactive SE group.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- $\text{C}_{56}\text{H}_{70}\text{N}_4\text{O}_{16}$
- MW: 1055



This datasheet was generated on June 3, 2026 at 09:31:19 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-carboxy-x-rhodamine-peo8-propionate-succinimidyl-ester-5-rox-peo8-se/>