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.



Call us: 800-304-5357 Email: btinfo@biotium.com

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 fluoresc amine-reactive SE group.

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- \bullet C₅₆H₇₀N₄O₁₆
- MW: 1055

This datasheet was generated on November 3, 2025 at 08:05:15 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/