

Sulforhodamine 101-PEO8 SE

SR101-PEO₈ SE (Sulforhodamine 101-PEO₈-propionic acid succinimydyl ester; Texas Red®-PEO₈ SE) has an approximately 32 Angstrom water-soluble PEG spacer that separates the dye from the amine-reactive SE group.



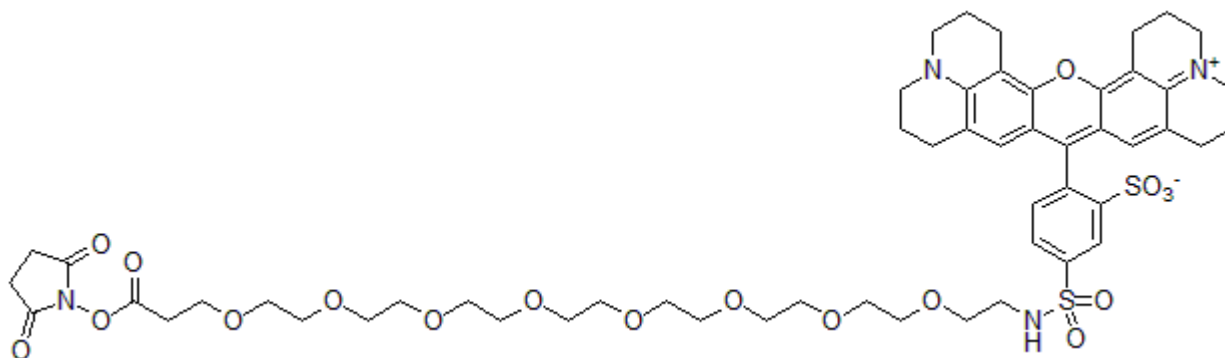
Product attributes

Excitation/Emission	583/603 nm
---------------------	------------

Product Description

SR101-PEO₈ SE (Sulforhodamine 101-PEO₈-propionic acid succinimydyl ester; Texas Red®-PEO₈ SE) has an approximately 32 Angstrom water-soluble PEG spacer that separates the dye from the amine-reactive SE group.

- $\lambda_{Ex}/\lambda_{Em}$ = 583/603 nm
- Dark purple solid soluble in DMF or DMSO
- Store at -20 °C and protect from light, especially in solution
- C₅₄H₇₀N₄O₁₈S₂
- MW: 1127.28



Texas Red is a registered trademark of Thermo Fisher Scientific.

This datasheet was generated on June 13, 2026 at 11:04:50 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/sulforhodamine-101-peo8-propionic-acid-succinimydyl-ester-sr101-peo8-se-or-texas-red-peo8-se/>