

Sulforhodamine 101-PEO12 SE

SR101-PEO₁₂ SE (Sulforhodamine 101-PEO₁₂-propionic acid succinimidylyl ester; Texas Red®-PEO₁₂ SE) has an approximately 47 Angstrom water-soluble PEG spacer that separates the dye from the amine-reactive SE group.



Product attributes

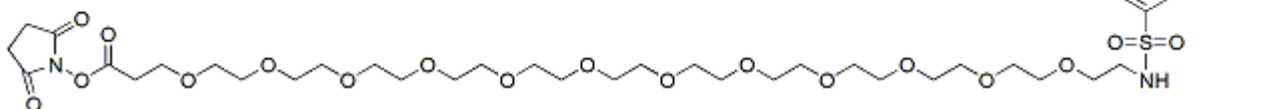
Excitation/Emission

583/603 nm

Product Description

SR101-PEO₁₂ SE (Sulforhodamine 101-PEO₁₂-propionic acid succinimidylyl ester; Texas Red®-PEO₁₂ SE) has an approximately 47 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
- C₆₂H₈₆N₄O₂₂S₂
- MW: 1303.49



Texas Red is a registered trademark of Thermo Fisher Scientific.

This datasheet was generated on June 3, 2026 at 09:30:16 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/sulforhodamine-101-peo12-propionic-acid-succinimidylyl-ester-sr101-peo12-se-or-texas-red-peo12-se/>