

Sulforhodamine 101 (Texas Red®)-PEO3-Amine

Sulforhodamine 101-PEO₃-amine has a long water-soluble spacer separating the dye and the amine group. The dye can be coupled to activated carboxy groups or sulfonyl chlorides.



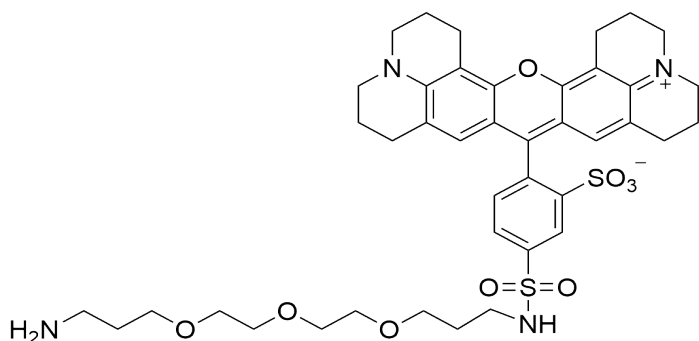
Product attributes

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

Product Description

Sulforhodamine 101-PEO₃-amine has a long water-soluble spacer separating the dye and the amine group. The dye can be coupled to activated carboxy groups or sulfonyl chlorides.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 583/603 nm
- Dark purple solid soluble in DMSO or DMF
- Store at 4°C and protect from light
- C₄₁H₅₂N₄O₉S₂
- MW: 809



This datasheet was generated on June 14, 2026 at 09:07:17 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/sulforhodamine-101-peo3-amine-texas-redtm-peo3-amine/>