

## Sulforhodamine 101 (Texas Red®)-PEO3-Amine

nfortum 680

Excitation/Emission

Call us: 800-304-5357

**Product attributes** 

583/603 nm

Sulforhodamine 101-PEO<sub>3</sub>-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 Description**

Sulforhodamine  $101-PEO_3$ -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
- $\bullet$  C<sub>41</sub>H<sub>52</sub>N<sub>4</sub>O<sub>9</sub>S<sub>2</sub>
- MW: 809

This datasheet was generated on November 23, 2025 at 12:37:25 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/sulforhodamine-101-peo3-amine-texas-redtm-peo3-amine/">https://biotium.com/product/sulforhodamine-101-peo3-amine-texas-redtm-peo3-amine/</a>