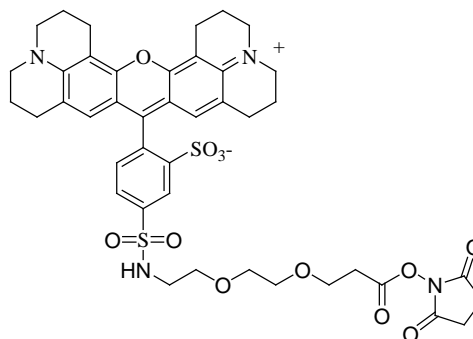


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** **Sulforhodamine 101-PEO<sub>2</sub>-propionic acid succinimydyl ester** (Texas Red-PEO<sub>2</sub>-P SE)

**CATALOG #** 90039

**MOLECULAR INFORMATION:** C<sub>42</sub>H<sub>46</sub>N<sub>4</sub>O<sub>12</sub>S<sub>2</sub>  
MWt: 862.26

**PROPERTIES:**

**Color & Form** Dark purple solid  
**Solubility** Soluble in DMF or DMSO  
**Absorption/Emission** 583/603 nm  
**Extinction Coefficient** 104,000

**STORAGE AND HANDLING:**

Store desiccated at -20 °C. Protect from light, especially in solution.

**APPLICATION:**

Biotium developed Sulforhodamine101-EO<sub>2</sub>-P SE as a superior alternative to the popular amine-reactive dye sulforhodamine 101 sulfonyl chloride (also called Texas Red sulfonyl chloride). Sulforhodamine 101 sulfonyl chloride undergoes hydrolysis fairly rapidly under normal conditions used for labeling amines in aqueous solutions, making it difficult to store and use. On the other hand, SR101-EO<sub>2</sub>-P SE reacts with amines readily but with a much slower hydrolysis rate. In addition, the long spacer between the fluorophore and the carboxylic acid group enhances the fluorescence of the dye and minimizes the interaction between the dye and the biomolecule to be labeled. SR101-EO<sub>2</sub>-P SE is also a better alternative to Texas Red-X SE for protein labeling because the former has a more water-soluble ethylene oxide linker that reduces dye aggregation, a common cause for fluorescence quenching.

**TOXICITY:**

Not established. Not listed by NTP, IARC or OSHA.

**FIRST AID:**

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.