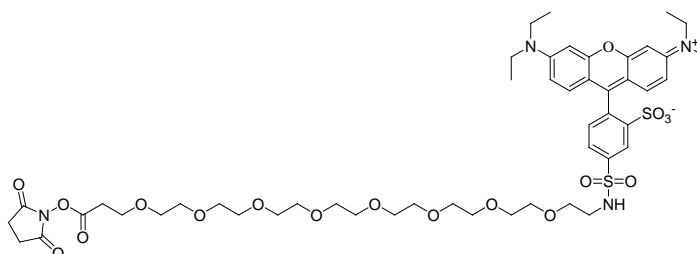


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Sulforhodamine B-PEO₈-propionic acid succinimidy ester**
(SRB-PEO₈SE)

CATALOG # **90101**

MOLECULAR INFORMATION: C₅₀H₇₀N₄O₁₈S₂
Mwt: 1079.24

**PROPERTIES:**

Color & Form Dark purple solid
Purity ≥ 95% by TLC
Solubility Soluble in DMSO or DMF
Absorption/Emission 560/580nm (MeOH)

STORAGE AND HANDLING: Store at -20 °C and protect from light, especially in solution.

APPLICATION: Sulforhodamine B-PEO₈ SE has an approximately 32 Angstrom water-soluble PEG spacer that separates the dye from the amine-reactive SE group

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.*