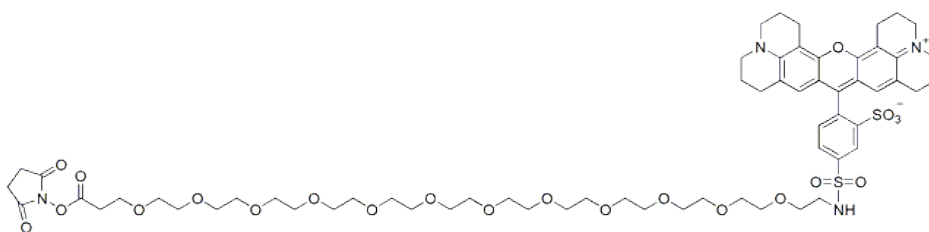


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Sulforhodamine 101-PEO12 SE

CATALOG # 90100

MOLECULAR INFORMATION: C₆₂H₈₆N₄O₂₂S₂
Mwt: 1303.49



PROPERTIES:

Color & Form Dark purple solid
Solubility Soluble in DMF or DMSO
Absorption/Emission 583/603 nm

STORAGE AND HANDLING:

Store at -20°C and protect from light, especially in solution

APPLICATION:

SR101-PEO₁₂ SE (Sulforhodamine 101-PEO₁₂-propionic acid succinimide ester; Texas Red-PEO₁₂ SE) has an approximately 47 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.