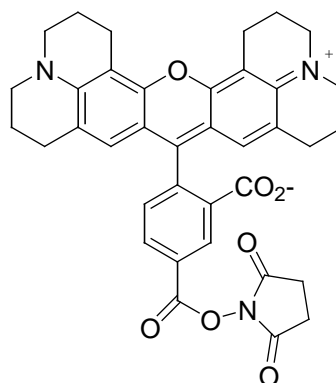


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Carboxy-X-rhodamine, succinimidyl ester, single isomer (5-ROX, SE)

CATALOG # 90036

MOLECULAR INFORMATION: C₃₇H₃₃N₃O₇
Mwt: 631.7

**PROPERTIES:**

Color & Form Dark red solid
Solubility Soluble in DMSO or DMF
Absorption/Emission 568/595nm (MeOH)
Extinction Coefficient

STORAGE AND HANDLING: We recommend using anhydrous DMSO to make stock solution. The solution is good for at least one month if stored at -20°C.
For long-term storage, store the solid desiccated at -20 °C and protect from light.

APPLICATION: Amine-reactive form of 5-carboxy-X-rhodamine single isomer.

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.
