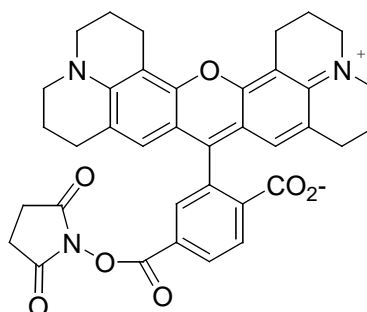


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **6-Carboxy-X-rhodamine, succinimidyl ester, single isomer (6-ROX, SE)**

CATALOG # 90037

MOLECULAR INFORMATION: C₃₇H₃₃N₃O₇
Mwt: 631.7
[216699-36-4]

**PROPERTIES:**

Color & Form Dark purple solid
Solubility Soluble in DMSO or DMF
Absorption/Emission 576nm/601nm (MeOH)
Extinction Coefficient 80,000

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

APPLICATION: 6-ROX is widely used in DNA sequencing and DNA probes based on fluorescent energy transfer (FRET). The succinimidyl ester reacts readily with primary or secondary amines under mild conditions. A stock solution of the dye can be made in anhydrous DMF or DMSO. The coupling reaction is usually carried out in a pH ~8.5 buffer (for example, 0.1-0.2 M sodium bicarbonate buffer). Alternatively, the coupling reaction may also be carried out in an organic solvent for amino compounds that are insoluble in aqueous solution.

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: <i>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.</i>
--