

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352 www.biotium.com

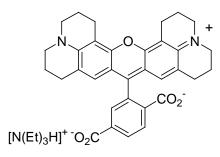
PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:

6-Carboxy-X-rhodamine, triethylammonium salt, single isomer (6-ROX)

CATALOG #

90007



PROPERTIES: Color & Form Purity Solubility Absorption/Emission	Dark red solid ≥ 95% by TLC Soluble in DMSO, DMF or MeOH 568/595nm (MeOH)
Extinction	100,000
Coefficient STORAGE AND HANDLING:	Store at 4°C. For long-term storage, store at -20 °C and protect from light.
APPLICATION:	6-Carboxy-X-rhodamine is a single isomer, long wavelength rhodamine dye.
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.*