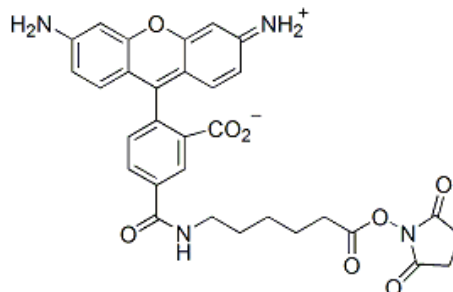


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Carboxyrhodamine 110-X succinimidyl ester, 5-isomer

CATALOG # 90014

MOLECULAR INFORMATION: C₃₁H₂₈N₄O₈
Mwt: 584.58



PROPERTIES:

Color & Form Orange red solid
Solubility Soluble in DMSO or DMF
Absorption/Emission 502/524 nm

STORAGE AND HANDLING: Store at -20°C and protect from light.

APPLICATION: 5-CR110-X, SE (full name: 5-Carboxyrhodamine 110-X, succinimidyl ester) is the amine-reactive form of 5-carboxyrhodamine 110 with a 6-carbon spacer.

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.