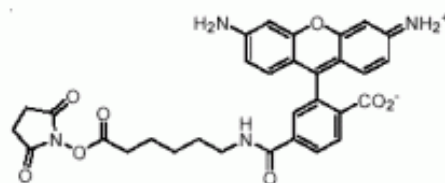


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Carboxyrhodamine 110-X succinimidyl ester, 6-isomer
(6-CR110-X, SE)

CATALOG # 90015

MOLECULAR INFORMATION: C₃₁H₂₈N₄O₈
MW: 584.58

**PROPERTIES:**

Color & Form Orange red solid soluble in DMSO or DMF
Purity ≥ 95% by TLC
Solubility Soluble DMF or DMSO
Absorption/Emission 502 nm/524 nm (MeOH)

STORAGE AND HANDLING:

Store desiccated at 4 °C. Store at -20°C when in solution and protect from light.

APPLICATION:

6-Carboxyrhodamine 110-X succinimidyl ester is the amine-reactive form of 6-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.