Carboxy-X-rhodamine succinimidyl ester, 5 isomer

5-ROX, SE (5-Carboxy-X-rhodamine, succinimidyl ester, single isomer) is the amine-reactive form of 5-carboxy-X-rhodamine single isomer.

ation and a second

Product attributes

Excitation/Emission

568/595 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

5-ROX, SE (5-Carboxy-X-rhodamine, succinimidyl ester, single isomer) is the amine-reactive form of 5-carboxy-X-rhodamine single isomer.

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- C₃₇H₃₃N₃O₇
- MW: 632
- [344402-35-3]

This datasheet was generated on November 3, 2025 at 09:13:22 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-carboxy-x-rhodamine-succinimidyl-ester-single-isomer5-rox-se/