

Carboxy-X-rhodamine succinimidyl ester, 6-isomer

6-ROX, SE (6-Carboxy-X-rhodamine, succinimidyl ester, single isomer) is the amine-reactive form of 6-carboxy-X-rhodamine single isomer.



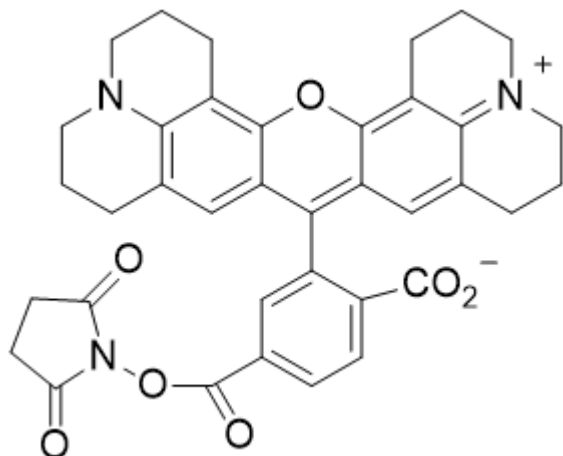
Product attributes

CAS number	216699-36-4
Excitation/Emission	568/595 nm

Product Description

6-ROX, SE (6-Carboxy-X-rhodamine, succinimidyl ester, single isomer) is the amine-reactive form of 6-carboxy-X-rhodamine single isomer.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 576/601 nm
- Dark red solid soluble in DMSO or DMF
- Store desiccated at -20 °C and protect from light
- $C_{37}H_{33}N_3O_7$
- MW: 632
- [344402-35-3]



This datasheet was generated on February 3, 2026 at 10:35:38 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/6-carboxy-x-rhodamine-succinimidyl-ester-single-isomer6-rox-se/>