

Carboxy-X-rhodamine succinimidyl ester, 5 and 6 isomers

5(6)-ROX, SE (full name: 5-(and-6)-Carboxy-X-rhodamine, succinimidyl ester) is an amine-reactive long wavelength rhodamine dye.



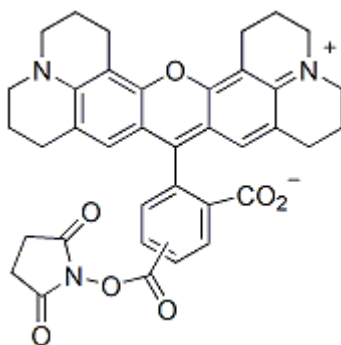
Product attributes

| | |
|---------------------|-------------|
| CAS number | 114616-32-9 |
| Excitation/Emission | 568/595 nm |

Product Description

5(6)-ROX, SE (full name: 5-(and-6)-Carboxy-X-rhodamine, succinimidyl ester) is an amine-reactive long wavelength rhodamine dye.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 575/600 nm
- Dark red solid soluble in DMSO or DMF
- Store at -20 °C and protect from light
- $C_{37}H_{33}N_3O_7$
- MW: 632
- [114616-32-9]



This datasheet was generated on June 3, 2026 at 06:06:15 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/5-and-6-carboxy-x-rhodamine-succinimidyl-ester-mixed-isomers-56-rox-se/>