

## 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.



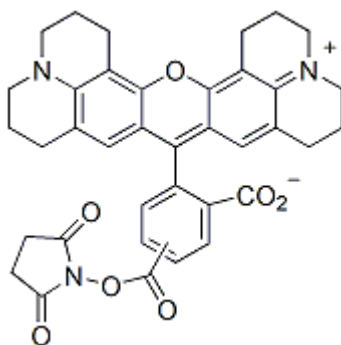
### Product attributes

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 January 28, 2026 at 12:50:25 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/>