

## 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.



### Product attributes

Excitation/Emission	568/595 nm
---------------------	------------

### 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.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO or DMF
- Store at  $-20^{\circ}\text{C}$  and protect from light
- $\text{C}_{37}\text{H}_{33}\text{N}_3\text{O}_7$
- MW: 632
- [344402-35-3]

