

Carboxy-X-rhodamine triethylammonium salt, 5 and 6 isomers (5(6)-ROX)



Product attributes

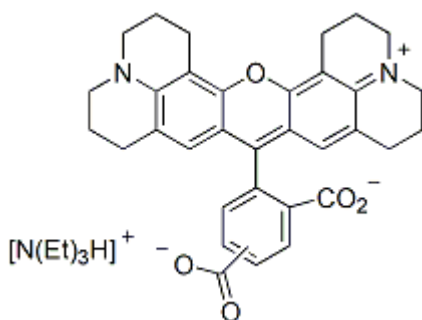
Excitation/Emission	568/595 nm
CAS number	198978-94-8

5(6)-ROX [full name: 5-(and-6)-Carboxy-X-rhodamine, triethylammonium salt (mixed isomers)] is a long wavelength rhodamine dye.

Product Description

5(6)-ROX (full name: 5-(and-6)-Carboxy-X-rhodamine, triethylammonium salt (mixed isomers)) is a long wavelength rhodamine dye.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 568/595 nm
- Dark red solid soluble in DMSO, DMF or MeOH
- Store at -20 °C and protect from light
- C₃₉H₄₅N₃O₅
- MW: 635.6
- [198978-94-8]



This datasheet was generated on June 18, 2026 at 07:21:20 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/5-and-6-carboxy-x-rhodamine-triethylammonium-salt-mixed-isomers-56-rox/>