

Carboxy-X-rhodamine triethylammonium salt, 5-isomer

5-ROX (full name: 5-Carboxy-X-rhodamine, triethylammonium salt, single isomer) is a single isomer, long wavelength rhodamine dye.



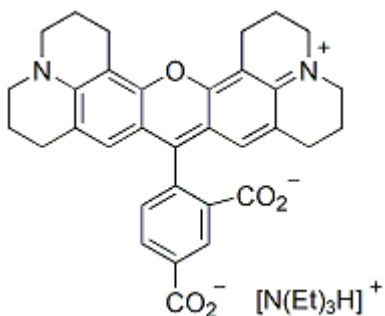
Product attributes

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

Product Description

5-ROX (full name: 5-Carboxy-X-rhodamine, triethylammonium salt, single isomer) is a single isomer, long wavelength rhodamine dye.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO, DMF or MeOH
- Store at 4 °C and protect from light
- $\text{C}_{39}\text{H}_{45}\text{N}_3\text{O}_5$
- MW: 635.6



This datasheet was generated on June 23, 2026 at 11:17:16 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-carboxy-x-rhodamine-triethylammonium-salt-single-isomer5-rox/>