

Carboxy-X-rhodamine triethylammonium salt, 6-isomer

6-ROX (6-Carboxy-X-rhodamine) is a single isomer, long wavelength rhodamine dye.



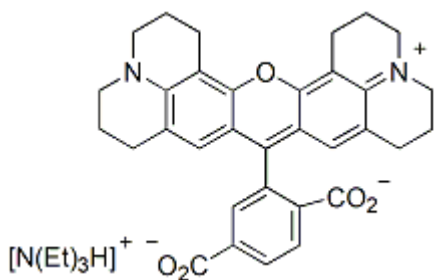
Product attributes

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

Product Description

6-ROX (6-Carboxy-X-rhodamine) 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 January 31, 2026 at 09:38:00 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/6-carboxy-x-rhodamine-triethylammonium-salt-single-isomer6-rox/>