

Carboxy-X-rhodamine free acid, 6-isomer

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



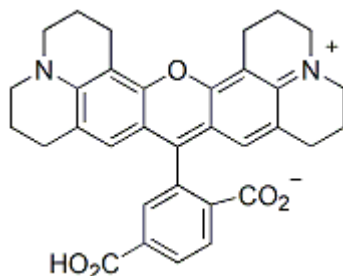
Product attributes

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

Product Description

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 -20°C and protect from light
- $\text{C}_{33}\text{H}_{30}\text{N}_2\text{O}_5$
- MW: 534.6



This datasheet was generated on February 2, 2026 at 10:01:19 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/6-carboxy-x-rhodamine-free-acid-6-rox-free-acid/>