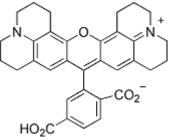
Carboxy-X-rhodamine free acid, 6-isomer

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

Product Description

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

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO, DMF or MeOH
- Store at -20°C and protect from light
- C₃₃H₃₀N₂O₅
- MW: 534.6





Product attributes

Excitation/Emission

568/595 nm



This datasheet was generated on November 3, 2025 at 10:11:36 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/6-carboxy-x-rhodamine-free-acid-6-rox-free