

## Carboxy-X-rhodamine triethylammonium salt, 5-isomer

otichan 680

Product attributes

Call us: 800-304-5357

Excitation/Emission 568/595 nm

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

## **Product Description**

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

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 568/595 \text{ nm}$
- Dark red solid soluble in DMSO, DMF or MeOH
- Store at 4°C and protect from light
- C<sub>39</sub>H<sub>45</sub>N<sub>3</sub>O<sub>5</sub>
- MW: 635.6

This datasheet was generated on November 19, 2025 at 04:14:29 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/5-carboxy-x-rhodamine-triethylammonium-salt-single-isomer5-rox/">https://biotium.com/product/5-carboxy-x-rhodamine-triethylammonium-salt-single-isomer5-rox/</a>