

Carboxy-X-rhodamine succinimidyl ester, 5 and 6 isomers

5(6)-ROX, SE (full name: 5-(and-6)-Carboxy-X-rhodamine, succinimidyl ester) is an amine-reactive long wavelength rhodamine dye.

stotum 80

Product attributes

Call us: 800-304-5357

CAS number 114616-32-9 **Excitation/Emission** 568/595 nm

Product Description

5(6)-ROX, SE (full name: 5-(and-6)-Carboxy-X-rhodamine, succinimidyl ester) is an amine-reactive long wavelength rhodamine dye.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 575/600 nm
- Dark red solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- C₃₇H₃₃N₃O₇
- MW: 632
- [114616-32-9]

This datasheet was generated on November 26, 2025 at 11:38:31 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-and-6-carboxy-x-rhodamine-succinimidyl-ester-mixed-isomers-56-rox-se/