ROX, SE, 6 isomer in DMSO

6-ROX, SE is the amine-reactive form of 6-ROX single isomer (90037) dissolved in anhydrous DMSO, suitable for labeling oligonucleotides.



Product attributes

CAS number 216699-36-4 **Excitation/Emission** 576/601 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

6-ROX, SE is the amine-reactive form of 6-ROX single isomer (90037) dissolved in anhydrous DMSO, suitable for labeling oligonucleotides.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 576/601 nm
- Dark red solution
- Store at -20°C and protect from moisture and light
- C37H33N3O7
- MW: 632
- [216699-36-4]

This datasheet was generated on November 4, 2025 at 10:34:19 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/6-carboxy-x-rhodamine-succinimidyl-ester-83mgml-0-13m-in-anhydrous-dmso-6-rox-se-in-dmso/