

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.



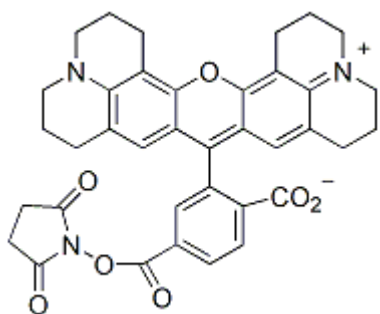
Product attributes

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

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
- $C_{37}H_{33}N_3O_7$
- MW: 632
- [216699-36-4]



This datasheet was generated on June 3, 2026 at 09:29:55 PM. 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-dms-6-rox-se-in-dms->