## TAMRA SE, 6 isomer in DMSO

6-TAMRA SE (full name: 6-Carboxytetramethylrhodamine succinimidyl ester) is the amine-reactive form of 6-TAMRA single isomer (90035).



## **Product attributes**

Excitation/Emission

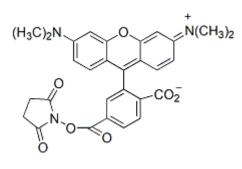
540/565 nm

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

## **Product Description**

6-TAMRA SE (full name: 6-Carboxytetramethylrhodamine succinimidyl ester) is the amine-reactive form of 6-TAMRA single isomer (90035). Supplied at 90 mg/mL (0.17 M) in anhydrous DMSO. The solution is suitable for labeling oligonucleotides.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 540/565 nm
- Dark red solution
- Store desiccated at -20°C and protect from light
- C29H25N3O7
- MW: 527.5



This datasheet was generated on November 4, 2025 at 10:30:51 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/6-carboxytetramethylrhodamine-succinimidyl-ester-90mgml-0-17m-in-anhydrous-dmso-6-tamra-se-in-dmso/">https://biotium.com/product/6-carboxytetramethylrhodamine-succinimidyl-ester-90mgml-0-17m-in-anhydrous-dmso-6-tamra-se-in-dmso/</a>