

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](#)).



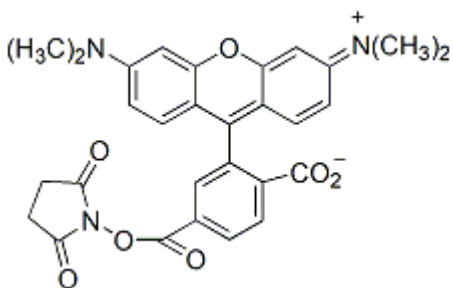
Product attributes

Excitation/Emission	540/565 nm
---------------------	------------

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
- $C_{29}H_{25}N_3O_7$
- MW: 527.5



This datasheet was generated on June 6, 2026 at 09:56:52 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/6-carboxytetramethylrhodamine-succinimidyl-ester-90mgml-0-17m-in-anhydrous-dms0-6-tamra-se-in-dms0/>