

TAMRA SE, 6-isomer

Full name: 6-Carboxytetramethylrhodamine succinimidyl ester single isomer. Amine-reactive form of carboxytetramethylrhodamine (TAMRA) single isomer.



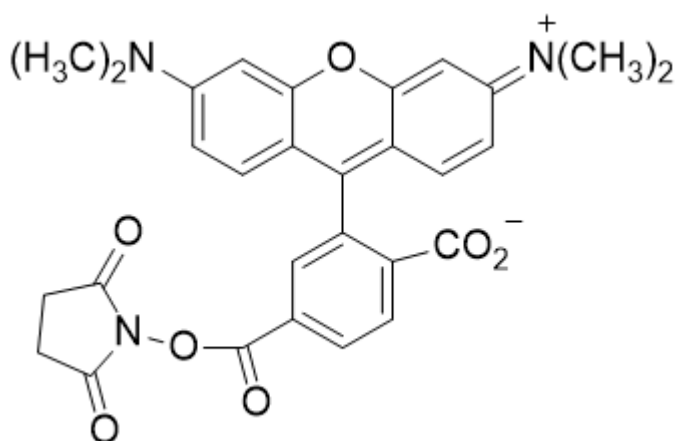
Product attributes

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

Product Description

Full name: 6-Carboxytetramethylrhodamine succinimidyl ester single isomer. Amine-reactive form of carboxytetramethylrhodamine (TAMRA) single isomer.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 540/565 nm
- Dark red solid soluble in DMSO or DMF
- Store at -20 °C and protect from light
- $C_{29}H_{25}N_3O_7$
- MW: 527.5
- [150810-69-8]



This datasheet was generated on January 9, 2026 at 04:01:01 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/6-carboxytetramethylrhodamine-succinimidyl-ester-single-isomer-6-tamra-se/>