

TAMRA SE, 5 isomer

Full name: 5-Carboxytetramethylrhodamine succinimidyl ester (single isomer). Amine-reactive form of carboxytetramethylrhodamine (TAMRA) single isomer.



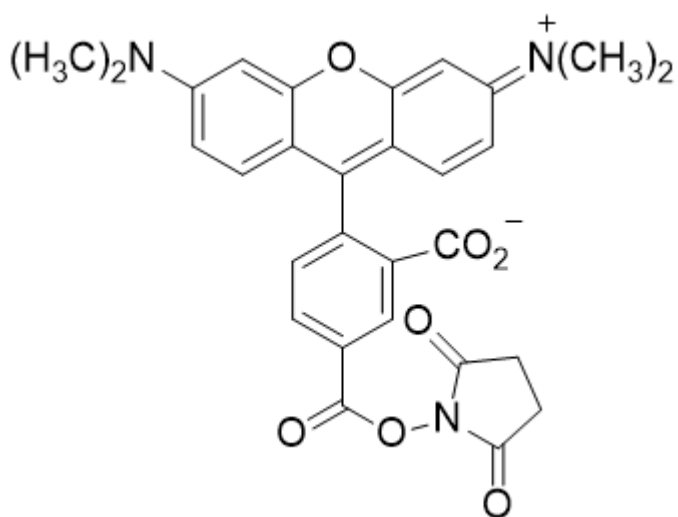
Product attributes

CAS number	150810-68-7
Excitation/Emission	540/565 nm

Product Description

Full name: 5-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
- [150810-68-7]



This datasheet was generated on June 3, 2026 at 06:07:07 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-carboxytetramethylrhodamine-succinimidyl-ester-single-isomer-5-tamra-se/>