TAMRA SE, 6-isomer

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



Product attributes

Excitation/Emission

540/565 nm

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

Product Description

Full name: 6-Carboxytetramethylrhodamine succinimidyl ester single isomer. Amine-reactive form ofcarboxytetramethylrhodamine (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
- C₂₉H₂₅N₃O₇
- MW: 527.5
- [150810-69-8]

This datasheet was generated on November 6, 2025 at 07:30:00 AM. 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/