## 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

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

## **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<sub>29</sub>H<sub>25</sub>N<sub>3</sub>O<sub>7</sub>
- MW: 527.5
- [150810-68-7]

This datasheet was generated on November 4, 2025 at 10:50:39 AM. 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/