TAMRA, 5-isomer

Free acid form of 5-TAMRA (5-Carboxytetramethylrhodamine, single isomer).

nicotum 80

Product attributes

 CAS number
 91809-66-4

 Excitation/Emission
 540/565 nm

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

Product Description

Free acid form of 5-TAMRA (5-Carboxytetramethylrhodamine, single isomer).

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 540/565 nm
- Dark red solid soluble in DMSO, DMF, or MeOH
- C₃₁H₃₇N₅O₅
- MW: 531.5
- [91809-66-4]

$$(H_3C)_2N \longrightarrow 0 \\ CO_2^-$$

$$CO_2H$$

This datasheet was generated on November 4, 2025 at 10:50:47 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-carboxytetramethylrhodamine-single-isomer-5-tamra/