

TAMRA, 5-isomer

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



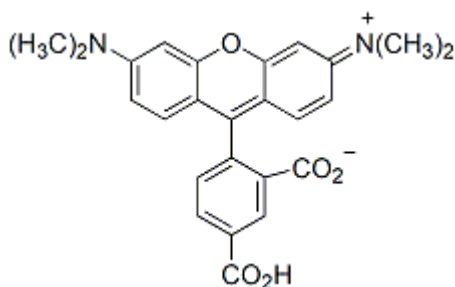
Product attributes

CAS number	91809-66-4
Excitation/Emission	540/565 nm

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
- $C_{31}H_{37}N_5O_5$
- MW: 531.5
- [91809-66-4]



This datasheet was generated on December 25, 2025 at 09:02:00 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/5-carboxytetramethylrhodamine-single-isomer-5-tamra/>