

## TAMRA, 6-isomer

Free acid form of 6-TAMRA (6-carboxytetramethylrhodamine) single isomer.



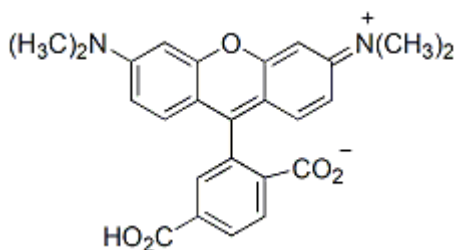
### Product attributes

CAS number	91809-67-5
Excitation/Emission	540/565 nm

## Product Description

Free acid form of 6-TAMRA (6-carboxytetramethylrhodamine) single isomer.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 540/565 nm
- Dark red solid soluble in DMSO, DMF, MeOH, or pH>6.5 buffers
- $C_{25}H_{22}N_2O_5$
- MW: 430.5
- [91809-67-5]



This datasheet was generated on June 10, 2026 at 04:52:50 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/6-carboxytetramethylrhodamine-single-isomer-6-tamra/>