

## TAMRA free acid, 5 and 6-isomer

Mixed isomers of carboxytetramethylrhodamine (TAMRA) free acid. The dye can be coupled to primary or secondary amines via standard peptide chemistry.



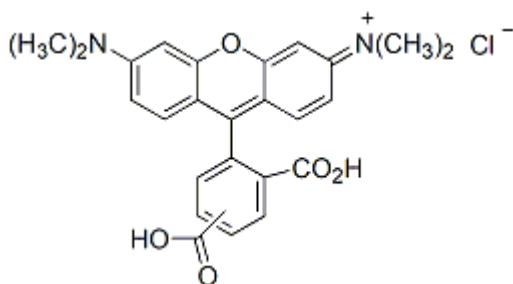
### Product attributes

Excitation/Emission	540/565 nm
---------------------	------------

## Product Description

Mixed isomers of carboxytetramethylrhodamine (TAMRA) free acid. The dye can be coupled to primary or secondary amines via standard peptide chemistry.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 540/565 nm
- Dark red solid soluble in DMSO, DMF, MeOH, or pH >6.5 buffer
- Store at -20 °C and protect from light
- $C_{25}H_{23}ClN_2O_5$
- MW: 466.92



This datasheet was generated on December 15, 2025 at 06:20:23 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/5-and-6-carboxytetramethylrhodamine-mixed-isomers-56-tamra/>