

Tetramethylrhodamine maleimide, 5 and 6 isomers

5(6)-Carboxytetramethylrhodamine maleimide has C6-linker that enhances the fluorescence after the dye is conjugated to proteins. The product is a direct replacement for 5(6)-carboxytetramethylrhodamine C5-maleimide (91040).



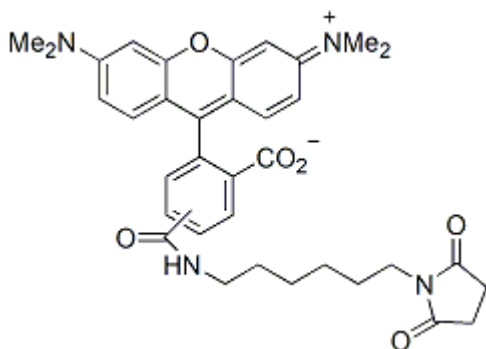
Product attributes

Excitation/Emission	540/565 nm

Product Description

5(6)-Carboxytetramethylrhodamine maleimide has C6-linker that enhances the fluorescence after the dye is conjugated to proteins. The product is a direct replacement for 5(6)-carboxytetramethylrhodamine C5-maleimide (91040)

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 540/565 nm
- Red solid soluble in DMF or DMSO
- Store at -20 °C and protect from light, especially in solution
- $C_{35}H_{36}N_4O_6$
- MW: 608.66



This datasheet was generated on December 5, 2025 at 02:40:23 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/56-tetramethylrhodamine-maleimide/>