

ROX maleimide, 5 and 6 isomers

5(6)-ROX maleimide has C6-linker that enhances the fluorescence after the dye is conjugated to proteins. The product is a direct replacement of 5(6)-carboxy-X-rhodamine C5-maleimide (91041).



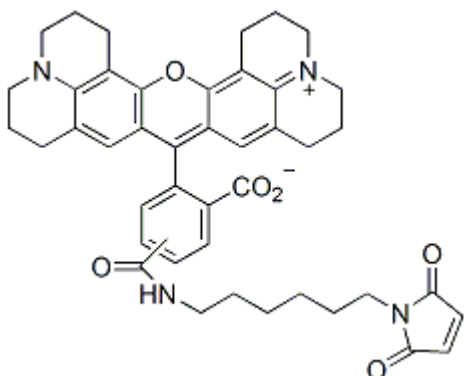
Product attributes

Excitation/Emission	568/595 nm
---------------------	------------

Product Description

5(6)-ROX maleimide has C6-linker that enhances the fluorescence after the dye is conjugated to proteins. The product is a direct replacement of 5(6)-carboxy-X-rhodamine C5-maleimide (91041).

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 568/595 nm
- Red solid soluble in DMF or DMSO
- Store at -20°C and protect from light, especially in solution
- $C_{43}H_{44}N_4O_6$
- MW: 712.81



This datasheet was generated on June 22, 2026 at 12:19:41 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/56-rox-maleimide/>