Carboxyrhodamine 110-C6-Maleimide, 5 and 6 isomers

Excitation/Emission

Product attributes

502/524 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

CR110-C6-maleimide (full name: 5(6)-Carboxyrhodamine 110-C6-maleimide) is a maleimide derivative of the green fluorescent dye carboxyrhodamine 110.

Product Description

CR110-C6-maleimide (full name: 5(6)-Carboxyrhodamine 110-C6-maleimide) is a maleimide derivative of the green fluorescent dye carboxyrhodamine 110. The dye has spectral properties ($\lambda_{\text{Ex}}/\lambda_{\text{Em}}$ (MeOH) = 502/524 nm) similar to those of fluorescein (fluorescein-5-maleimide, for example), but the photostability of the dye is among the best of all green fluorescent dyes. Thus, the dye is a superior green fluorescent dye for labeling thiols.

- Orange red solid soluble in DMF or DMSO
- Store at -20°C and protect from light, especially in solution
- C₃₁H₂₈N₄O₆
- MW: 552.59

This datasheet was generated on November 5, 2025 at 01:34:06 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/56-carboxyrhodamine-110-c5-maleimide-56-cr110-c5-maleimide/