

Carboxyrhodamine 110-C6-Maleimide, 5 and 6 isomers

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



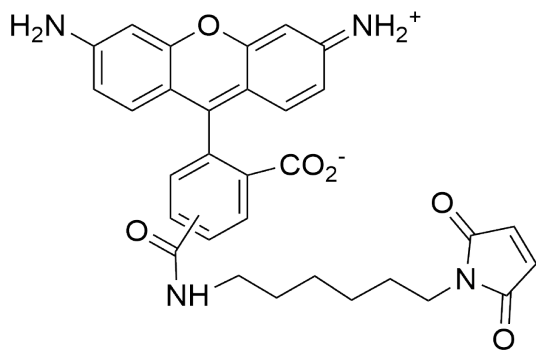
Product attributes

Excitation/Emission	502/524 nm
---------------------	------------

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
- $\text{C}_{31}\text{H}_{28}\text{N}_4\text{O}_6$
- MW: 552.59



This datasheet was generated on January 22, 2026 at 05:50:53 PM. 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/>