

Carboxyrhodamine 110-C6-Maleimide, 5 and 6 isomers



Product attributes

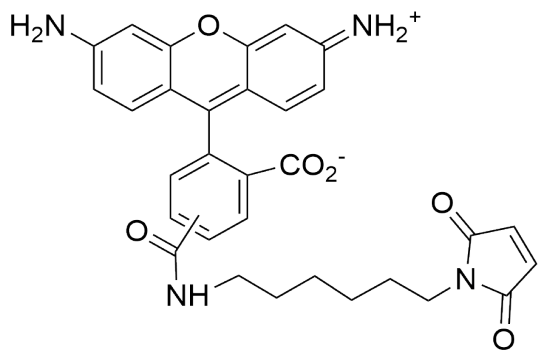
Excitation/Emission	502/524 nm
---------------------	------------

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 ($\lambda_{Ex}/\lambda_{Em}$ (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
- $C_{31}H_{28}N_4O_6$
- MW: 552.59



This datasheet was generated on June 3, 2026 at 07:35:11 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/>