

Carboxyrhodamine 110 succinimidyl ester, 5-isomer

Dicitum Pictum

Product attributes

Excitation/Emission 502/524 nm

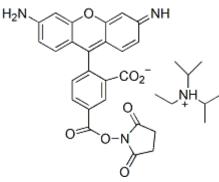
Call us: 800-304-5357

Amine-reactive form of single isomer carboxyrhodamine 110, useful in DNA sequencing. For many single-dye-labeling applications, the dye is preferred over other green fluorescent dyes including fluorescein and Alexa Fluor® 488 because of its exceptional photostability and fluorescence insensitivity to pH.

Product Description

Amine-reactive form of single isomer carboxyrhodamine 110, useful in DNA sequencing. For many single-dye-labeling applications, the dye is preferred over other green fluorescent dyes including fluorescein and Alexa Fluor® 488 because of its exceptional photostability and fluorescence insensitivity to pH between 4 and 9.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 502/524 nm
- Orange red solid soluble in DMF or DMSO
- Store desiccated at -20°C and protect from light
- C33H36N4O7
- MW: 600.67



Alexa Fluor is a registered trademark of Thermo Fisher Scientific.

This datasheet was generated on November 30, 2025 at 10:56:33 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-carboxyrhodamine-110-succinimidyl-ester-single-isomer-5-cr110-se/