Carboxyrhodamine 110 succinimidyl ester, 6-isomer

Amine-reactive form of single isomer carboxyrhodamine 110. For certain applications, the dye is preferred over the corresponding 6-carboxyfluorescein because of its exceptional photostability and fluorescence insensitivity to pH change (4-9).

nicitum

Product attributes

Excitation/Emission

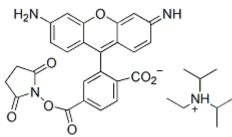
502/524 nm

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

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.

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 502/524 \text{ nm}$
- 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 5, 2025 at 12:37:49 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/6-carboxyrhodamine-110-succinimidyl-ester-single-isomer-6-cr110-se/