

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).



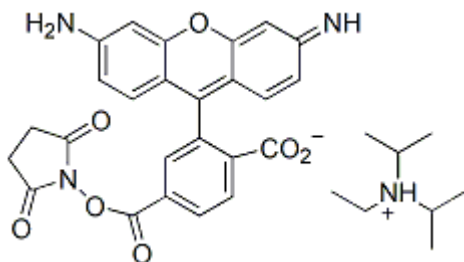
Product attributes

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

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.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 502/524 \text{ nm}$
- Orange red solid soluble in DMF or DMSO
- Store desiccated at -20°C and protect from light
- $\text{C}_{33}\text{H}_{36}\text{N}_4\text{O}_7$
- MW: 600.67



Alexa Fluor is a registered trademark of Thermo Fisher Scientific.

This datasheet was generated on June 6, 2026 at 03:39:52 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/>