

## 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).



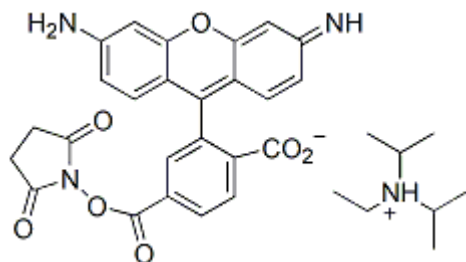
### Product attributes

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  $-20^{\circ}\text{C}$  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.