

Carboxyrhodamine 110-X succinimidyl ester, 5 and 6 isomers

5-(and-6)-Carboxyrhodamine 110-X, succinimidyl ester (5(6)-CR 110-X, SE) readily couples to a primary or secondary amine. The dye is a superior alternative to 5(6)-SFX because of its exceptional photostability and fluorescence insensitivity to pH (4-9).



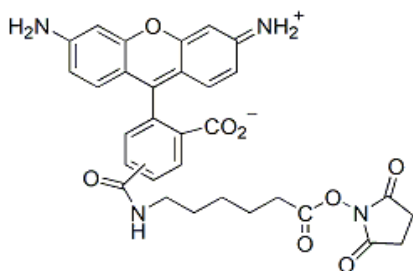
Product attributes

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

Product Description

5-(and-6)-Carboxyrhodamine 110-X, succinimidyl ester (5(6)-CR 110-X, SE) readily couples to a primary or secondary amine. The dye is a superior alternative to 5(6)-SFX because of its exceptional photostability and fluorescence insensitivity to pH (4-9).

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 502/524 nm
- Orange red solid soluble in DMSO or DMF.
- Store at -20°C and protect from light
- $C_{31}H_{28}N_4O_8$
- MW: 584.58



This datasheet was generated on June 23, 2026 at 07:42:51 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/5-and-6-carboxyrhodamine-110-x-succinimidyl-ester-56-cr-110-x-se/>