

Carboxyrhodamine 110-X succinimidyl ester, 5-isomer

5-CR110-X, SE (full name: 5-Carboxyrhodamine 110-X, succinimidyl ester) is the amine-reactive form of 5-carboxyrhodamine 110 with a 6-carbon spacer.



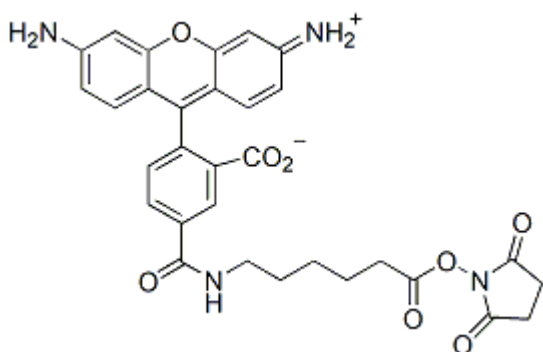
Product attributes

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

Product Description

5-CR110-X, SE (full name: 5-Carboxyrhodamine 110-X, succinimidyl ester) is the amine-reactive form of 5-carboxyrhodamine 110 with a 6-carbon spacer.

- $\lambda_{Ex}/\lambda_{Em}$ (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 January 14, 2026 at 02:50:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-carboxyrhodamine-110-x-succinimidyl-ester-5-cr110-xse/>