## Carboxyrhodamine 110 succinimidyl ester, 5 and 6 isomers

5-(and-6)-Carboxyrhodamine 110, SE (5-(and-6)-Carboxyrhodamine 110, succinimidyl ester, mixed isomers), also known as Carboxyrhodamine 110-NHS, is a superior green fluorescent dye that readily reacts with primary or secondary amines

## alcitum

## Product attributes

Excitation/Emission

502/524 nm

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

## **Product Description**

5-(and-6)-Carboxyrhodamine 110, SE (5-(and-6)-Carboxyrhodamine 110, succinimidyl ester, mixed isomers), also known as Carboxyrhodamine 110-NHS, is a superior green fluorescent dye that readily reacts with primary or secondary amines. For many applications, the dye is preferred over 5-(and-6)-carboxyfluorescein succinimidyl ester (5(6)-FAM) or FITC because of its exceptional photostability and fluorescence insensitivity to pH between pH 4 and pH 9.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 502/524 nm
- Orange red solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C25H17N3O7
- MW:471.42

This datasheet was generated on November 10, 2025 at 07:51:38 AM. Visit product page to check for updated information before use.

Product link: <a href="https://biotium.com/product/5-and-6-carboxyrhodamine-110-succinimidyl-ester-mixed-isomers-56-cr-110-se-or-carboxyrhodamine-110-nhs/">https://biotium.com/product/5-and-6-carboxyrhodamine-110-nhs/</a>