# 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

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

## **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<sub>31</sub>H<sub>28</sub>N<sub>4</sub>O<sub>8</sub>
- MW: 584.58

$$H_2N$$
 $CO_2$ 
 $CO_2$ 

This datasheet was generated on November 10, 2025 at 05:31:38 AM. 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/