

## Carboxyrhodamine 110-PEO3-amine, TFA salt, 5 isomer

5-CR110-PEO<sub>3</sub>-amine has a long water-soluble PEG spacer that separates the green fluorescent CR110 and the amine functional group.



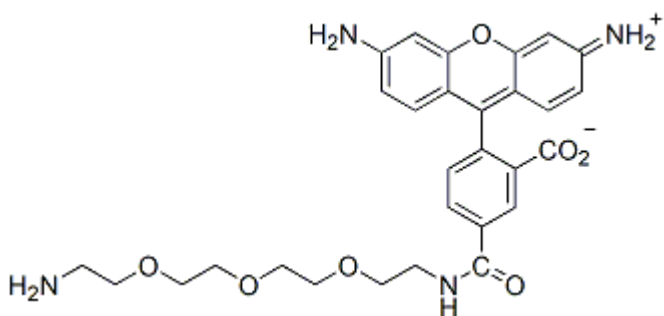
### Product attributes

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

### Product Description

5-CR110-PEO<sub>3</sub>-amine has a long water-soluble PEG spacer that separates the green fluorescent CR110 and the amine functional group.

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 502/524 nm
- Orange yellow solid soluble in DMSO, DMF, or H<sub>2</sub>O
- Store at -20 °C and protect from light
- C<sub>33</sub>H<sub>34</sub>F<sub>6</sub>N<sub>4</sub>O<sub>11</sub>
- MW: 776.63



This datasheet was generated on June 22, 2026 at 02:02:41 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/5-cr110-peo3-amine-tfa-salt/>