

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

5-CR110-PEO₃-amine has a long water-soluble PEG spacer that separates the green fluorescent CR110 and the amine functional group.



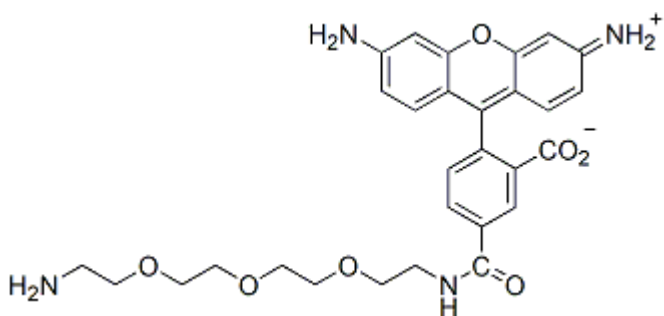
Product attributes

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

Product Description

5-CR110-PEO₃-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 H₂O
- Store at -20 °C and protect from light
- C₃₃H₃₄F₆N₄O₁₁
- MW: 776.63



This datasheet was generated on December 19, 2025 at 06:33:40 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/5-cr110-peo3-amine-tfa-salt/>