

Carboxyrhodamine 110-X hydrochloride, 5 and 6 isomers



Product attributes

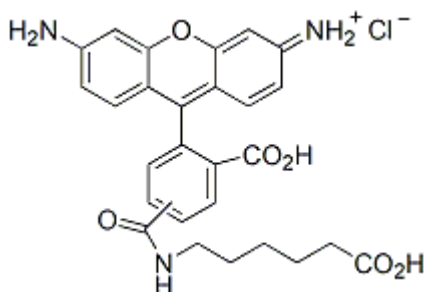
Excitation/Emission	502/524 nm
---------------------	------------

5-(and-6)-Carboxyrhodamine 110-X (full name: 5-(and-6)-Carboxyrhodamine 110-X, hydrochloride) is similar to 5-(and-6)-carboxyrhodamine 110 ([#90000](#)).

Product Description

5-(and-6)-Carboxyrhodamine 110-X (full name: 5-(and-6)-Carboxyrhodamine 110-X, hydrochloride) is similar to 5-(and-6)-carboxyrhodamine 110 ([#90000](#)). The X linker molecule is useful to reduce the interaction between the dye and the molecule to be labeled.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 502/524 nm
- Orange red solid soluble in water, DMF or DMSO
- Store at 4 °C and protect from light
- $C_{28}H_{30}ClN_3O_6$
- MW:524



This datasheet was generated on June 17, 2026 at 03:14:39 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/5-and-6-carboxyrhodamine-110-x-hydrochloride-56-cr-110-x/>