Carboxyrhodamine 110-X hydrochloride, 5 and 6 isomers

5-(and-6)-Carboxyrhodamine 110-X (full name:

5-(and-6)-Carboxyrhodamine 110-X, hydrochloride) is similar to

5-(and-6)-carboxyrhodamine 110 (#90000).

nicotum 80

Product attributes

Excitation/Emission

502/524 nm

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

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
- C₂₈H₃₀CIN₃O₆
- MW:524

This datasheet was generated on November 11, 2025 at 06:53:24 AM. 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/