

Carboxyrhodamine 6G hydrochloride, 5-isomer

5-CR6G (full name: 5-Carboxyrhodamine 6G, hydrochloride, single isomer) is the free acid form of 5-carboxyrhodamine 6G single isomer.



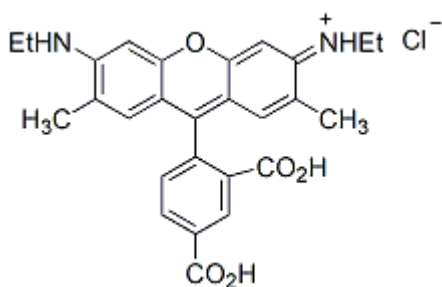
Product attributes

Excitation/Emission	520/546 nm
---------------------	------------

Product Description

5-CR6G (full name: 5-Carboxyrhodamine 6G, hydrochloride, single isomer) is the free acid form of 5-carboxyrhodamine 6G single isomer.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 520/546 \text{ nm}$
- Orange red solid soluble in water (pH 7), DMF or DMSO
- Store at 4°C and protect from light
- $\text{C}_{27}\text{H}_{27}\text{ClN}_2\text{O}_5$
- MW: 494.97



This datasheet was generated on June 23, 2026 at 05:11:30 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-carboxyrhodamine-6g-hydrochloride-single-isomer-5-cr6g/>