

## Carboxyrhodamine 6G hydrochloride, 6-isomer

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



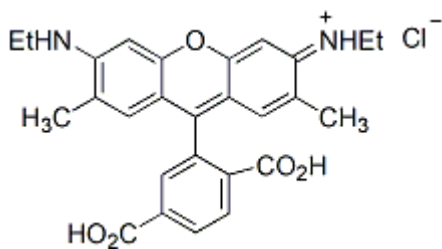
### Product attributes

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

### Product Description

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

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 520/546 nm
- Orange red solid soluble in water (pH 7), DMF or DMSO
- Store at 4°C and protect from light
- $C_{27}H_{27}ClN_2O_5$
- MW: 494.97



This datasheet was generated on June 3, 2026 at 05:58:06 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/6-carboxyrhodamine-6g-hydrochloride-single-isomer-6-cr6g/>