

5-(and-6)-Carboxyrhodamine 6G, Hydrochloride

The free acid form of orange fluorescent 5-(and-6)-carboxyrhodamine 6G mixed isomers.



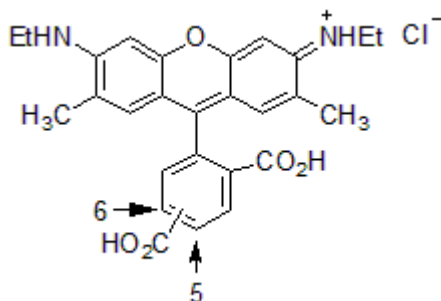
Product attributes

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

Product Description

5-(and-6)-Carboxyrhodamine 6G, hydrochloride, is the free acid form of orange fluorescent 5-(and-6)-carboxyrhodamine 6G mixed isomers. Also called 5(6)-CR6G.

- Ex/Em (MeOH) = 520/546 nm
- Orange red solid moderately soluble in DMSO
- Store at -20 °C, protect from light
- MW: 494.97



This datasheet was generated on December 7, 2025 at 08:07:26 PM. Visit product page to check for updated information before use.

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