Carboxyrhodamine 6G hydrochloride, 6-isomer

atictum 8

Excitation/Emission 520/546 nm

Product attributes

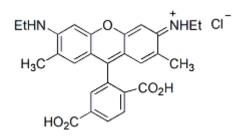
Call us: 800-304-5357 Email: btinfo@biotium.com

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

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
- C₂₇H₂₇CIN₂O₅
- MW: 494.97



This datasheet was generated on November 5, 2025 at 01:34:02 AM. Visit product page to check for updated information before use. Product link: $\frac{\text{https://biotium.com/product/6-carboxyrhodamine-6g-hydrochloride-single-isomer-6-cr6g/}{\text{https://biotium.com/product/6-carboxyrhodamine-6g-hydrochloride-single-isomer-6-cr6g/}}$