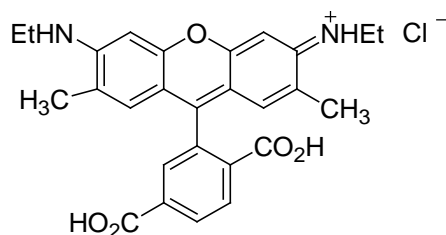


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 6-Carboxyrhodamine 6G, hydrochloride, single isomer (6-CR 6G)**CATALOG #** 90011**MOLECULAR INFORMATION:** C₂₇H₂₇ClN₂O₅
Mwt:494.97**PROPERTIES:**

Color & Form Orange red solid
Purity ≥ 95% by HPLC
Solubility Soluble in water (pH 7), DMF or DMSO
Absorption/Emission 520nm/546nm (MeOH)
Extinction Coefficient 100,000

STORAGE AND HANDLING: Store desiccated at 4 °C. Protect from light, especially when in solution.**APPLICATION:** 6-Carboxyrhodamine 6G, hydrochloride, single isomer is the free acid form of 6-carboxyrhodamine 6G single isomer.**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--