



Product Information

Carboxyrhodamine 6G, Hydrochloride

Unit Size: 5 mg

Catalog no.	Product
90011	6-CR6G (6-Carboxyrhodamine 6G, hydrochloride, single isomer)
90012	5-CR6G (5-Carboxyrhodamine 6G, hydrochloride, single isomer)
90120	5(6)-CR6G (5-(and-6)-Carboxyrhodamine 6G, hydrochloride, mixed isomers)

Storage and Handling

Store at 4°C desiccated and protected from light. Product is stable for at least 5 years from date of receipt when stored as recommended.

Stock solutions may be prepared in DMF or DMSO. Solutions can be stored with desiccant and protected from light at 4° C, for up to 4 years.

Molecular Information: C₂₇H₂₇CIN₂O₅

Molecular Weight: 495

Color and Form: Orange red solid **Solubility:** Soluble in DMF or DMSO

Absorption/Emission: 520/546 nm (MeOH)

Figure 1. 5(6)-CR6G (5-(and-6)-Carboxyrhodamine 6G, hydrochloride, mixed isomers)

Product Description

Free acid forms of orange fluorescent 5-CR6G (5-carboxyrhodamine 6G, hydrochloride, single isomer), 6-CR6G (6-carboxyrhodamine 6G, hydrochloride, single isomer), and 5(6)-CR6G (5-(and-6)-carboxyrhodamine 6G hydrochloride, mixed isomers).

Related Products

Catalog number	Product
90082	DMSO, Anhydrous
90032	Carboxyrhodamine 6G Succinimidyl Ester, 5-isomer
90033	Carboxyrhodamine 6G Succinimidyl Ester, 6-isomer
90000	Carboxyrhodamine 110, Hydrochloride, 5 and 6 isomers
90030	Carboxyrhodamine 110 Succinimidyl Ester, 5-isomer
90031	Carboxyrhodamine 110 Succinimidyl Ester, 6-isomer
90001	Carboxyrhodamine 110 Succinimidyl Ester, 5 and 6 isomers
92103	CF®514 Succinimidyl Ester
80500	CFQ520 Succinimidyl Ester
59002	EDC (EDAC)
29052	ROX Reference Dye, 25 uM in TE Buffer

Please visit our website at www.biotium.com to view our wide selection of fluorescent dyes including our bright and photostable CF® dyes. Also learn more about our other life science research products, including environmentally friendly EvaGreen® qPCR master mixes, antibody conjugates, apoptosis reagents, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.