

Carboxyrhodamine 6G succinimidyl ester, 6-isomer

Amine-reactive form of 6-carboxyrhodamine 6G single isomer.



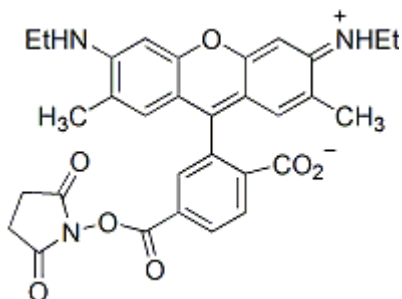
Product attributes

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

Product Description

Amine-reactive form of 6-carboxyrhodamine 6G single isomer.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 520/546 nm
- Orange red solid moderately soluble in DMSO
- Store desiccated at -20 °C and protect from light
- $C_{31}H_{29}N_3O_7$
- MW: 555.6



This datasheet was generated on June 3, 2026 at 06:06:14 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/6-carboxyrhodamine-6g-succinimidyl-ester-single-isomer6-cr6g-se/>