Carboxyrhodamine 6G succinimidyl ester, 6-isomer

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

siotum &

Product attributes

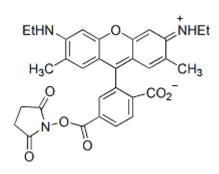
Excitation/Emission 520/546 nm

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

Product Description

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

- λεx/λεm (MeOH) = 520/546 nm
- Orange red solid moderately soluble in DMSO
- Store desiccated at -20°C and protect from light
- C₃₁H₂₉N₃O₇
- MW: 555.6



This datasheet was generated on November 3, 2025 at 02:14:54 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/