

Product Information Sheet

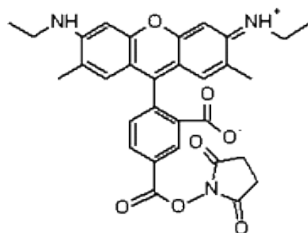
Ordering Information

Product Number: 345
Product Name: 5-CR6G, SE [5-Carboxyrhodamine 6G, succinimidyl ester]
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 555.59

Chemical Structure:



Soluble in: MeOH
Excitation Wavelength: 524 nm
Emission Wavelength: 556 nm

Application Notes

5-CR6G, SE is amine-reactive derivative of 5-CR6G. It is preferred for some complicated biological applications where reproducibility is more critical than material cost since the minor positional difference between 5-CR6G and 6-CR6G might significantly affect some biological properties of the underlying conjugates.