

5(6)-CR6G, SE [5-(and-6)-Carboxy-rhodamine 6G, succinimidyl ester]

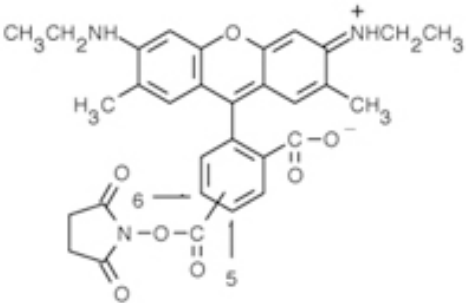
Catalog Number	Packaging Size
C130	5 mg

Storage upon receipt: -20°C, protect from light

Introduction

5(6)-CR6G, SE is a amine-reactive fluorescent probe for preparation of bioconjugates. The excitation and emission spectra of carboxyrhodamine 6G (CR 6G) fall between those of fluorescein and tetramethylrhodamine and with higher fluorescence quantum yield than tetramethylrhodamine conjugates. With a peak absorption at 525 nm, conjugated probes of carboxyrhodamine 6G are an excellent match to the 514 nm spectral line of the argon-ion laser.

Specifications

Label:	Rhodamine 6G	
Ex/Em:	525/550 nm	
Detection Method:	Fluorescent	
Solubility:	DMSO, DMF	
Molecular Formula:	C ₃₁ H ₂₉ N ₃ O ₇	
Molecular Weight:	555.59	
CAS Number:	349672-89-5	
Storage Conditions:	-20°C, protect from light	
Shipping Condition:	Room Temperature	

Applications

Fluorescent labeling

References:

- [Multiplex detection of single-nucleotide variations using molecular beacons.](#)
Marras SA, Kramer FR, Tyagi S
Genet Anal (1999) 14:151-156
- Intramolecular Excitonic Dimers in Protease Substrates: Modification of the Backbone Moiety to Probe the H-Dimer Structure.
Packard BZ, et al.
J Phys Chem B (1998) 102:1820-1820
- [Single-molecule detection and identification of multiple species by multiparameter fluorescence detection.](#)
Widengren J, Kudryavtsev V, Antonik M, Berger S, Gerken M, Seidel CA
Anal Chem (2006) 78:2039-2050