

Product Information

Carboxyrhodamine 110 Succinimidyl Ester, 6-Isomer

Catalog Number: 90031

Unit Size: 5 mg

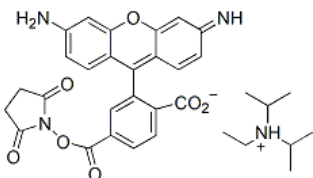
Storage and Handling

Store desiccated at -20°C, protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended. We recommend using anhydrous DMSO (see related products) to prepare stock solutions of succinimidyl ester dyes immediately before use.

Molecular Information

$C_{33}H_{36}N_4O_7$

MW: 600.67



Properties

Color & Form: Orange red solid

Solubility: Soluble in DMF or DMSO

Ex/Em (MeOH) = 502/524 nm

Product Description

Amine-reactive form of single isomer carboxyrhodamine 110, useful in DNA sequencing. For many single-dye-labeling applications, the dye is preferred over other green fluorescent dyes including fluorescein and Alexa Fluor® 488 because of its exceptional photostability and fluorescence insensitivity to pH between 4 and 9.

Related Products

Catalog number	Product
22004	Ultrafiltration vial, 10K MWCO (5 per pack)
22018	Ultrafiltration vial, 3K MWCO (5 per pack)
90082	DMSO, anhydrous
22013	Bovine Serum Albumin, Fraction V
22014	Bovine Serum Albumin, 30% solution
22020	10X Phosphate Buffered Saline
41024-4L	Water, Ultrapure Molecular Biology Grade

Please visit www.biotium.com to learn about our next-generation CF® dyes with superior brightness and photostability. We offer a wide selection of reactive dye chemistries, labeling kits, labeled antibodies, other conjugates, and more.

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