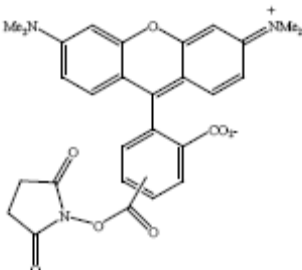


**5-(and-6)-Carboxytetramethylrhodamine succinimidyl ester
(mixed isomers)**
Instruction Manual

Catalog Number	PK-CA707-90022
Description	Carboxytetramethylrhodamine is one of the most commonly used red fluorescent dyes, which can be optimally excited by the 546 Hg line. The succinimidyl ester reacts readily with primary or secondary amines under mild conditions. PromoKine offers a very high grade TAMRA SE.
Quantity	25 mg
Excitation / Emission Maxima	$\lambda_{\text{ex}} / \lambda_{\text{em}} = 540/565 \text{ nm (MeOH)}$; Extinction Coefficient : 95,000
Molecular Structure	
Molecular Weight / Molecular Formula	527.5 Da; C ₂₉ H ₂₅ N ₃ O ₇
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Red solid. Soluble in DMF or DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	see Description
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.