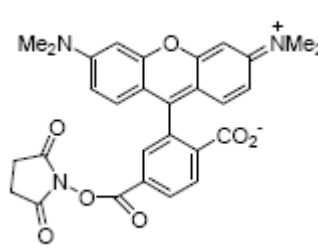


**6-Carboxytetramethylrhodamine succinimidyl ester;
Solution in DMSO**
Instruction Manual

Catalog Number	PK-CA707-90097
Description	Amine-reactive form of 6-carboxytetramethylrhodamine single isomer. Can be used for nucleotide labelling and automated DNA sequencing.
Quantity	100 µl
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	≥98% (as determined by TLC)
Appearance / Formulation / Solubility	Dark red liquid. 90 mg/ml (0.17M) in anhydrous DMSO. Dilute in DMF or DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
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.