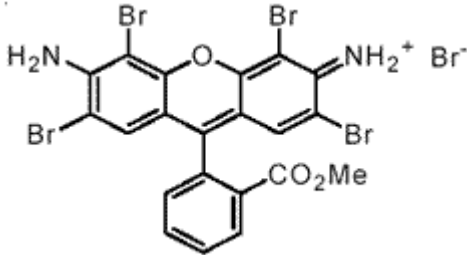


Instruction Manual

Catalog Number	PK-CA707-70013
Description	Tetrabromorhodamine 123 can be used to generate singlet oxygen (1O_2) in mitochondrial membranes. The dye has a quantum yield for singlet oxygen generation of 0.65-0.7 and is particularly toxic to carcinoma cells.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 524/550 nm Extinction coefficient (in MeOH) > 85,000
Molecular Structure	
Molecular Weight / Molecular Formula	740.86 Da; $C_{21}H_{13}Br_5N_2O_3$
Purity	≥95% as determined by HPLC analysis.
Appearance / Formulation / Solubility	Orange solid; soluble in DMSO or DMF.
Storage & Stability	Store at 4°C and protect from light, especially when in solution.
Applications	See Description
References	1) Photochem. Photobiol. 55, 81(1992) 2) Proc. SPIE 997, 48 (1988) 3) Cancer Res. 49, 3961(1989)
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.