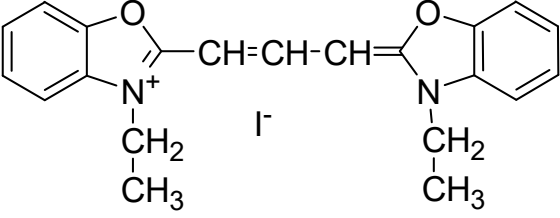


3,3'-Diethyloxocarboyanine, iodide

Instruction Manual

Catalog Number	PK-CA707-70008
Description	DiOC ₂ (3) has been used for measuring membrane potentials in bacteria. The green fluorescent dye forms red fluorescent aggregates with increasing membrane potential. The spectral shifting phenomenon has been exploited to measure membrane potential ratiometrically.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 482/497 nm Extinction coefficient = 150,000 (in MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	460.31 Da; C ₂₁ H ₂₁ N ₂ O ₂
Purity	≥98% (as determined by TLC)
Appearance / Formulation / Solubility	Orange solid; soluble in DMSO or DMF.
Storage & Stability	Store at ≤4°C. Protect from light, especially when in solution.
Applications	See Description
References	Biophysics J. 56, 979 (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.