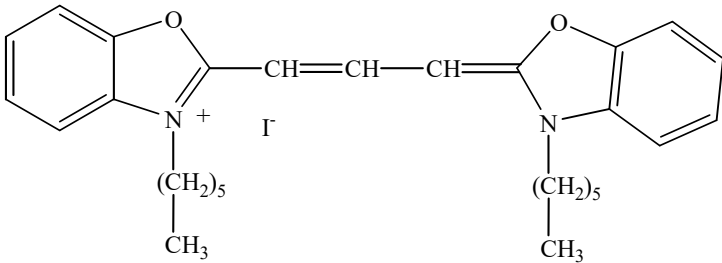


Instruction Manual

Catalog Number	PK-CA707-70009
Description	DiOC ₆ (3) is a green fluorescent membrane dye that has been used to stain the ER in both live and fixed cells. The dye has also been used to study structural interactions and dynamics of the ER in neurons and yeast. In addition, the dye is widely used as a membrane potential dye.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 484/501 nm Extinction coefficient = 150,000 (in MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	572.53 Da; C ₂₉ H ₃₇ IN ₂ O ₂
Purity	>98% (as determined by HPLC)
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	<ol style="list-style-type: none"> 1) J. Cell Biol. 103, 1557(1986) 2) Cell 38, 101(1984) 3) J. Cell Biol. 127, 1021(1994); 6050(1980) 4) Nature, 310, 53(1984) 5) Cell Motil. Cytoskeleton, 25, 111(1993)
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.