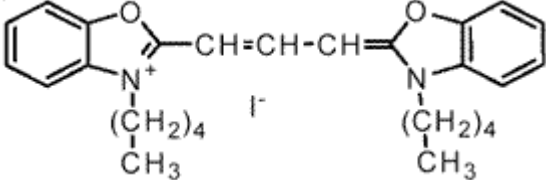


Instruction Manual

Catalog Number	PK-CA707-70007
Description	DiOC ₅ (3) and DiOC ₆ (3) are two of the most widely used carbocyanine dyes used for membrane potential measurements.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 482/497 nm
Molecular Structure	
Molecular Weight / Molecular Formula	544.47 Da; C ₂₇ H ₃₃ IN ₂ O ₂
Purity	NA
Appearance / Formulation / Solubility	Orange solid; soluble in DMSO or DMF.
Storage & Stability	Store at -20°C in a dry environment. Protect from light.
Applications	See Description
References	J. Biol. Chem. 270, 3788 (1995)
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.