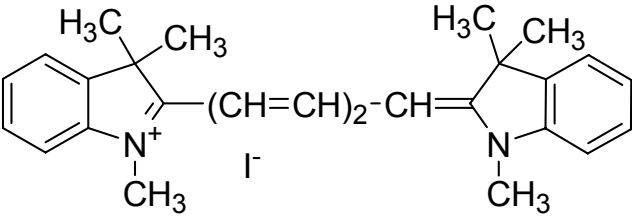


1,1'-dimethyl-3,3,3',3'-tetramethylindodicarbocyanine, iodide

Instruction Manual

Catalog Number	PK-CA707-70015
Description	The carbocyanine dye DiIC ₁ (5) has been used to measure mitochondrial membrane potentials in apoptotic cells. Loss of membrane potential in apoptotic cells is reflected by a decrease of fluorescence signal in the infrared channel.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 646/666 nm; Extinction coefficient = 226,000 (646 nm, in DMSO)
Molecular Structure	
Molecular Weight / Molecular Formula	510.46 Da; C ₂₇ H ₃₁ IN ₂
Purity	>99% (as determined by TLC)
Appearance / Formulation / Solubility	Dark green solid; soluble in DMSO.
Storage & Stability	Store desiccated at $\leq 4^{\circ}\text{C}$. Protect from light, especially when in solution.
Applications	See Description
References	Cytometry 33, 333(1998)
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.