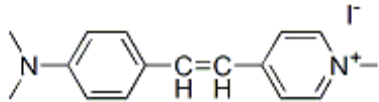


4-(4-(Dimethylamino)-styryl)-N-methylpyridinium iodide; stains mitochondria in live cells

Instruction Manual

Catalog Number	PK-CA707-70006
Description	4-Di-1-ASP is a styryl dye that can be used to stain mitochondria of live cells. It has also been used to stain glioma cells in living brain tissues. ¹
Quantity	500 mg
Excitation / Emission Maxima	$\lambda_{ex} / \lambda_{em} = 474/606$ nm (MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	366.2 Da; C ₁₆ H ₁₉ IN ₂
Purity	NA
Appearance / Formulation / Solubility	Red solid. Soluble in DMF and DMSO.
Storage & Stability	Store desiccated at 4°C. Protect from light, especially when in solution.
Applications	see Description
References	1) BioTechniques, doi:10.2144/000113940
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.