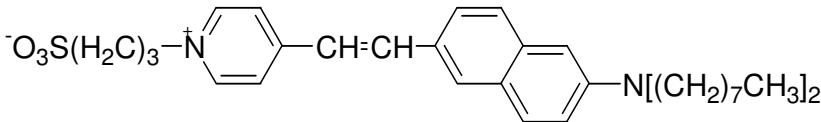


Fast-responding membrane-potential dye

Instruction Manual

Catalog Number	PK-CA707-61012
Description	ANEPPS dyes are fast responding membrane potential dyes. PromoKine offers both Di-4-ANEPPS and Di-8-ANEPPS. Di-4-ANEPPS has the problem of being internalized in cells over long-term, so it is mainly used for short-term experiments. However, because Di-8-ANEPPS is more lipophilic, the dye is better retained in the outer leaflet of cell plasma membranes and thus more suitable for long-term membrane potential studies. Di-8-ANEPPS is also more photostable and less phototoxic than Di-4-ANEPPS.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 498/713 nm; Extinction coefficient = 44,000 (in MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	593 kDa; C ₃₆ H ₅₂ N ₂ O ₃ S
Purity	>99% (as determined by HPLC)
Appearance / Formulation / Solubility	Orange red solid; soluble in EtOH or DMSO.
Storage & Stability	Store desiccated at $\leq 4^{\circ}\text{C}$. Protect from light, especially when in solution.
Applications	Membrane-potential dye
References	1) J. Neurosci. 15, 1392(1995) 2) Neuron 13,1187(1994) 3) Neuron,9,393(1992)
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.