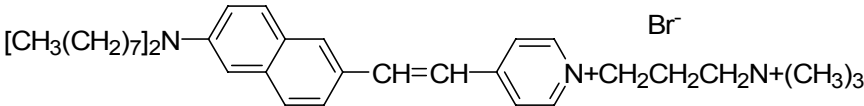


## Fast-responding membrane-potential dye

### Instruction Manual

Catalog Number	PK-CA707-61014
Description	Di-8-ANEPPQ is a fast-responding membrane potential dye. It is more hydrophobic than Di-2-ANEPEQ and is useful for potential-sensitive retrograde labeling of neurons.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 498/713 nm; Extinction coefficient = 38,000 (in MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	731 Da; C <sub>39</sub> H <sub>61</sub> Br <sub>2</sub> N <sub>3</sub>
Purity	>99% (as determined by HPLC)
Appearance / Formulation / Solubility	Dark-brown solid; soluble in DMF or DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	Membrane-potential dye
References	1) J. Neurosci. Methods. 70, 121(1996) 2) J. Neurosci. Methods 70, 111(1996)
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.