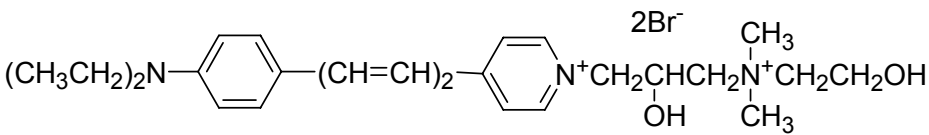


Fast-responding potentiometric probe

Instruction Manual

Catalog Number	PK-CA707-61019
Description	<p>The styryl dye RH795 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. The dye is spectrally similar to RH414 (Cat.No. PK-CA707-61016), but exhibits different physiological effects during staining. For example, RH414 causes arterial constriction during cortex staining, while RH795 does not.</p> <p>In cell membranes, spectra of styryl dyes are typically blue-shifted by as much as 20 nm for absorption or excitation and 80 nm for emission.</p>
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 530/712 nm ; Extinction coefficient = 58,000 (in MeOH)
Molecular Structure	
Molecular Weight / Molecular Formula	585.42 Da; C ₂₆ H ₃₉ Br ₂ N ₃ O ₂
Purity	NA
Appearance / Formulation / Solubility	Orange red solid; soluble water.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	Potentiometric probe
References	<p>1) Nature J. 324, 361(1986)</p> <p>2) J. Neurosci. 14, 2545(1994)</p>
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.