

Fast-responding potentiometric probe

Instruction Manual

Catalog Number	PK-CA707-61018
Description	The styryl dye RH237 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. 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.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 520/712 nm ; Extinction coefficient = 58,000 (in MeOH)
Molecular Structure	$[\text{CH}_3(\text{CH}_2)_3]_2\text{N}-\text{C}_6\text{H}_4-(\text{CH}=\text{CH})_3-\text{C}_6\text{H}_4-\text{N}^+(\text{CH}_2)_4\text{SO}_3^-$
Molecular Weight / Molecular Formula	496.71 Da; C ₂₉ H ₄₀ N ₂ O ₃ S
Purity	NA
Appearance / Formulation / Solubility	Dark solid; soluble in DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	Potentiometric probe
References	1) Neuroscience 31, 613 (1989) 2) The data is for methanol solution of the dye.
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.