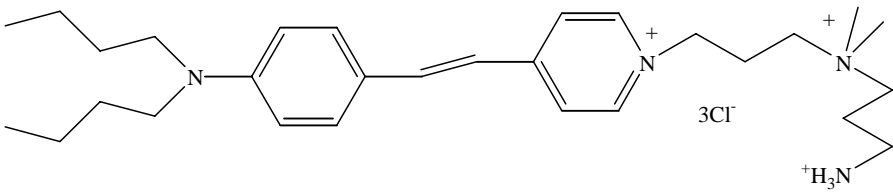


Instruction Manual

Catalog Number	PK-CA707-70024
Description	AM1-43 is a fixable, activity-dependent fluorescent nerve terminal probe and functions similarly as our SynapseGreen C4 (PK-CA707-70020; equivalent to FM1-43, a trademark of Molecular Probes, Inc.) does, but with the additional property of being fixable. Thus, AM1-43 is a useful tool for synapse studies where subsequent fluorescent immunochemistry is desired. AM1-43 and FM1-43 have identical excitation and emission wavelengths.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 510/625 nm; $\lambda_{ex}/\lambda_{em}$ (in membranes) = 480/598 nm Extinction coefficient = 56,000
Molecular Structure	
Molecular Weight / Molecular Formula	560.09 Da; C ₂₉ H ₄₉ Cl ₃ N ₄
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Dark-red solid; soluble in water.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	See Description
References	Renger, J.J., et al., A developmental switch in neurotransmitter flux enhances synaptic efficacy by affecting AMPA receptor activation. <i>Neuron</i> 29, 469, (2001).
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.