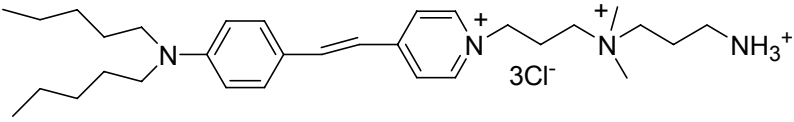


Instruction Manual

Catalog Number	PK-CA707-70038
Description	AM1-43 is a fixable, activity-dependent fluorescent nerve terminal probe and functions similarly as our SynapseGreen (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. The newly improved version, AM1-44 probe functions similarly as AM1-43, except that AM1-44 has one more lipophilic tail longer than that of AM1-43, resulting in better fixability. The dye can also be used as a general probe for following endocytosis.
Quantity	1 mg
Excitation / Emission Maxima	λ_{ex} \ λ_{em} : 510/625 nm (in MeOH); 480/598 nm (in membranes); $\epsilon = 56,000$
Molecular Structure	
Molecular Weight / Molecular Formula	587.5 Da; C ₃₁ H ₅₂ Cl ₃ N ₄
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Dark red solid; soluble in water and methanol.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	See Description
References	1) <i>Neuron</i> , 29 , 469 (2001) 2) <i>The Journal of Neuroscience</i> , 23 (10), 4054(2003)
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.