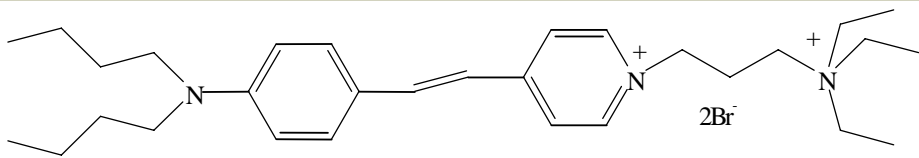


(also known as FM1-43*)

Instruction Manual

Catalog Number	PK-CA707-70022
Description	SynapseGreen C4 (also known as FM1-43, a trademark of Molecular Probes, Inc.) is a widely used green fluorescent dye for following synaptic activities by staining synaptic vesicles at the synapse or neuromuscular junctions. When used in combination with the red fluorescent dye SynapseRed C2 (PK-CA707-70021), synapses or neuromuscular junctions can be imaged independently in two colors. Recent research results from Tsien's and Kay's labs showed that brain slices can be stained with the styryl dye SynapseGreen when SR101 (Sulforhodamine 101, PK-CA707-80101) or ADVASEP-7 (PK-CA707-70029) is used to reduce the background staining. PromoKine offers several Nerve Terminal Staining Kits for staining brain slices.
Quantity	5 x 1 mg
Excitation / Emission Maxima	$\lambda_{ex}\backslash\lambda_{em}$ (in MeOH) = 510/625 nm; $\lambda_{ex}\backslash\lambda_{em}$ (in membranes) = 480/598 nm Extinction coefficient = 56,000
Molecular Structure	
Molecular Weight / Molecular Formula	612 Da; C ₃₀ H ₄₉ Br ₂ N ₃
Purity	>98% (as determined by HPLC)
Appearance / Formulation / Solubility	Red solid; soluble in water or DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	See Description
References	1) J. Neurosci. 12, 363(1992) 2) Science. 255, 200(1992) 2) Neuron 24, 803(1999) 3) Neuron 24, 809(1999)
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.

* FM1-43 is a trademark of Molecular Probes, Inc.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.