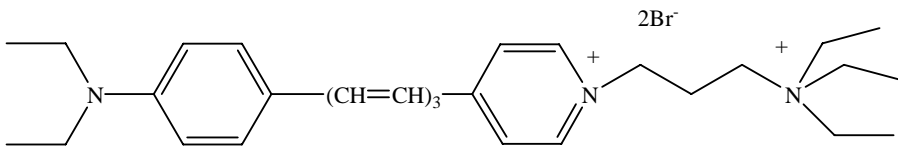


(equivalent to FM4-64\*)

## Instruction Manual

Catalog Number	PK-CA707-70021
Description	SynapseRed C2 (equivalent to FM4-64, a trademark of Molecular Probes, Inc.) is a widely used red fluorescent dye for following synaptic activities by staining synaptic vesicles at the synapse or neuromuscular junctions. When used in combination with the green fluorescent dye SynapseGreen C4 (PK-CA707-70020), synapses or neuromuscular junctions can be imaged independently in two colors. SynapseRed C2 has also been used to study endocytosis in Dictyostelia.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{exc}$ (in MeOH) = 543/too weak to detect; Extinction coefficient (in MeOH) = 43,000
Molecular Structure	
Molecular Weight / Molecular Formula	608 Da; C <sub>30</sub> H <sub>45</sub> Br <sub>2</sub> N <sub>3</sub>
Purity	>96% (as determined by HPLC)
Appearance / Formulation / Solubility	Dark purple 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) 3) J. Cell Biol. 121, 1311(1993)
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.

\* FM4-64 is a trademark of Molecular Probes, Inc.

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