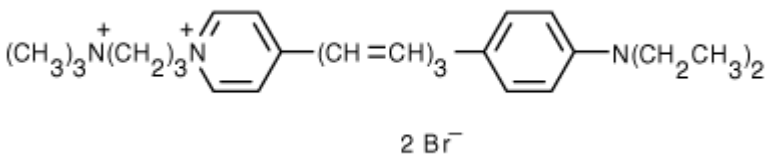


(also known as FM5-95*)

Instruction Manual

Catalog Number	PK-CA707-70019
Description	SynapseRed C2M (also known as FM5-95, a trademark of Molecular Probes, Inc.) is a red fluorescent probe used to follow synaptic vesicles. SynapseRed C2M is slightly more water soluble than SynapseRed C2 (also known as FM4-64, a trademark of Molecular Probes, Inc.).
Quantity	5 x 1 mg
Excitation / Emission Maxima	$\lambda_{exc}/\lambda_{em}$ (in MeOH) = 560/734 nm; Extinction coefficient (in MeOH) = 43,000
Molecular Structure	 <chem>C[N+](C)(C)C1=CC=C(C=C1)C=CC2=CC=C(C=C2)N(CC)CC.[Br-].[Br-]</chem>
Molecular Weight / Molecular Formula	565.43 Da; C ₂₇ H ₃₉ Br ₂ N ₃
Purity	>99% (as determined by TLC)
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) Science 297, 1525-1531 (2002) 2) Cell 107, 605-616 (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.

* FM5-95 is a trademark of Molecular Probes, Inc.

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