

Instruction Manual

Catalog Number	PK-CA707-70041
Description	SynapseRed C1 is similar to the widely used red fluorescent nerve terminal dye SynapseRed C2 (also called FM4-64, a trademark of Molecular Probes, Inc.) except that the lipophilic tail of the latter is one carbon longer than that of the former. Thus, SynapseRed C1 is expected to have a greater "off-rate" than SynapseRed C2 does.
Quantity	5 x 1 mg
Excitation / Emission Maxima	$\lambda_{exc}/\lambda_{em}$ (in MeOH) = 543/too weak to detect. Note: The excitation/emission of styryl dyes will undergo a blue shift and the dyes will become significantly more fluorescent in membranes.
Molecular Structure	<p>The chemical structure shows a central chain of three vinyl groups (-CH=CH-) connecting two benzene rings. The left benzene ring has a dimethylamino group (-N(CH₃)₂) at the para position. The right benzene ring has a quaternary ammonium group (-N⁺(CH₂)₃-N⁺(CH₂CH₃)₃) at the para position. Two bromide ions (2Br⁻) are shown as counterions.</p>
Molecular Weight / Molecular Formula	580 Da; C ₂₈ H ₄₁ Br ₂ N ₃
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Deep 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	NA
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.