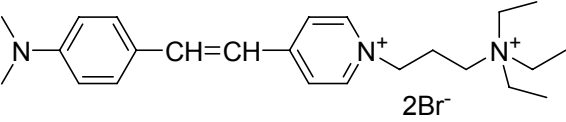


## Instruction Manual

Catalog Number	PK-CA707-70042
Description	SynapseGreen C1 is a fluorescent membrane probe that can be used to follow synaptic vesicle recycling. The dye is nonfluorescent in water but becomes highly fluorescent once incorporated into membranes. The dye's maximal emission is in the orange red spectral region. However, because of its relatively broad spectrum, it can also be seen in the green region. Among the series of nerve terminal dyes we offer, SynaptoGreen C1 has the shortest lipophilic tail (C1) and is the most water soluble of the nerve terminal dye series. Thus, it has the fastest de-staining rate, i.e. it is expected to have the slowest "on-rate" but the fastest "off-rate".
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex}$ \ $\lambda_{em}$ : 505/620 nm (in MeOH); 480/598 nm (in membranes); Extinction coefficient: ~60,000
Molecular Structure	
Molecular Weight / Molecular Formula	528 Da; C <sub>24</sub> H <sub>37</sub> Br <sub>2</sub> N <sub>3</sub>
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) Proc. R. Soc. Lond. B 255, 61(1994) 2) J Biol Chem 275, 15279-15286 (2000); 3) J Neurochem 75, 1645-1653 (2000) 4) Neuron 28, 221-231 (2000) 5) Proc Natl Acad Sci U S A 97, 12828-12833 (2000) 6) Neurosci Lett 253, 1-4 (1998) 7) Nature 394, 581-585 (1998)
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.