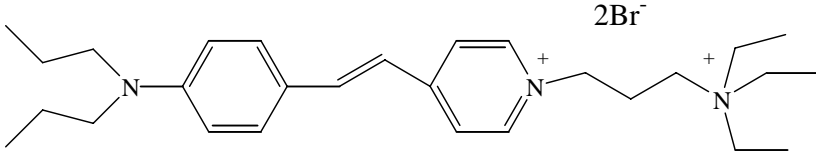


Instruction Manual

Catalog Number	PK-CA707-70026
Description	SynapseGreenC3 is similar to SynapseGreen C4 (Cat.No. PK-CA707-70020) except that the lipophilic tail is one carbon shorter and also the hydrophilic end is a trimethylammonium group. Thus, SynapseGreen C3 is slightly more water soluble than SynapseGreen.
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: NA
Molecular Structure	
Molecular Weight / Molecular Formula	542 Da; C ₂₅ H ₃₉ Br ₂ N ₃
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Dark red solid; soluble in water.
Storage & Stability	Store desiccated at -20°C.
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.

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