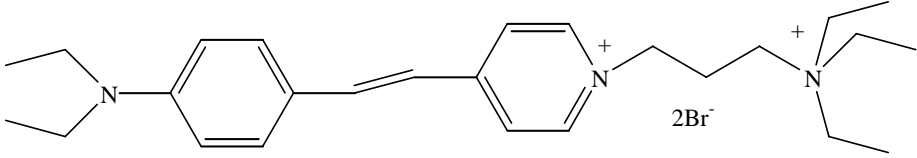


Nerve terminal probe (also known as FM2-10*)

Instruction Manual

Catalog Number	PK-CA707-70045
Description	SynapseGreen C2 (also known as FM2-10) is a green fluorescent probe used to follow synaptic vesicles. SynaptoGreen C2 is more water soluble than SynaptoGreen C4 and thus has a faster de-staining rate than the latter.
Quantity	5 x 1 mg
Excitation / Emission Maxima	λ_{exc} λ_{em} : 505/620 nm (in MeOH); 480/598 nm (in membranes); Extinction coefficient: ~60,000
Molecular Structure	
Molecular Weight / Molecular Formula	556 Da; C ₂₆ H ₄₁ Br ₂ N ₃
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.

* FM2-10 is a trademark of Molecular Probes, Inc.

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