

Nerve terminal probe (also known as FM1-84*)

Instruction Manual

Catalog Number	PK-CA707-70046
Description	SynapseGreen C5 (also known as FM1-84) is similar to SynapseGreen C4 (Cat.No. PK-CA707-70020) except that the lipophilic tail is one carbon longer. Thus SynapseGreen C5 is less water soluble than SynapseGreen C4 and thus faster staining rate but slower de-staining rate.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \setminus \lambda_{em}$: 510/625 nm (in MeOH); 480/598 nm (in membranes); Extinction coefficient = 56,000
Molecular Structure	$[\text{CH}_3(\text{CH}_2)_4]_2\text{N}-\text{C}_6\text{H}_4-\text{CH}=\text{CH}-\text{C}_6\text{H}_4-\text{N}^+(\text{CH}_2)_3\text{N}^+(\text{CH}_2\text{CH}_3)_3 \cdot 2\text{Br}^-$
Molecular Weight / Molecular Formula	640 Da; C ₃₂ H ₅₃ Br ₂ N ₃
Purity	>99% (as determined by HPLC)
Appearance / Formulation / Solubility	Dark orange solid; soluble in water and DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution. To dissolve the material, the solution mixture can be safely sonicated and heated up to 80°C.
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.

* FM1-84 is a trademark of Molecular Probes, Inc.

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