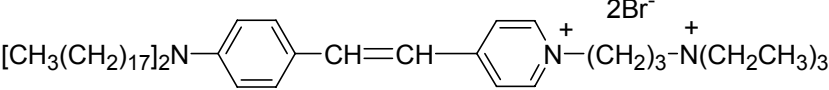


Nerve terminal probe (also known as FM3-25\*)

## Instruction Manual

Catalog Number	PK-CA707-70049
Description	SynapseGreen C18 (also known as FM3-25) is a membrane stain similar to other nerve terminal dyes (see our website), except that it has the longest lipophilic tail (C18).
Quantity	5 x 1 mg
Excitation / Emission Maxima	$\lambda_{exc}$ $\lambda_{em}$ : 510/625 nm (in MeOH); 480/598 nm (in membranes); Extinction coefficient = 56,000
Molecular Structure	 <p>The chemical structure shows a central trans-stilbene core. The left phenyl ring is substituted with a long-chain dialkylammonium group, <math>[CH_3(CH_2)_{17}]_2N</math>. The right phenyl ring is substituted with a trimethylammonium group, <math>N^+(CH_2)_3-N^+(CH_2CH_3)_3</math>. Two bromide ions, <math>2Br^-</math>, are shown as counterions.</p>
Molecular Weight / Molecular Formula	1004 Da; $C_{58}H_{105}Br_2N_3$
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Red solid; soluble in DMF and DMSO.
Storage & Stability	Store desiccated at $-20^\circ C$ . Protect from light, especially when in solution. To dissolve the material, the solution mixture can be safely sonicated and heated up to $80^\circ C$ .
Applications	See Description
References	Meyers, J.R. et al. J Neurosci., 23, 4054(2003)
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.

\* FM3-25 is a trademark of Molecular Probes, Inc.

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