

Instruction Manual

Catalog Number	PK-CA707-70051
Description	AM3-25 has been developed as a potentially fixable, activity-dependent fluorescent nerve terminal probe. AM3-25 is a lipophilic styryl dye that has an amine group at the hydrophilic head. The amine group can react with formaldehyde or glutaldehyde to make the dye fixable. Preliminary results have indicated that AM3-25 functions similarly as SynapseGreen C18 (also called FM3-25, PK-CA707-70049), except that AM3-25 has a fixable amine group. However, its full utility has not been thoroughly evaluated. AM3-25 and FM3-25 have identical excitation and emission wavelengths.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex} \setminus \lambda_{em}$: 510/625 nm (in MeOH); 480/598 nm (in membranes)
Molecular Structure	<p>The chemical structure shows a lipophilic styryl dye. It consists of a long-chain dialkylammonium group, $[CH_3(CH_2)_{17}]_2N$, attached to a para-substituted phenyl ring. This phenyl ring is connected via a vinyl bridge ($-CH=CH-$) to another para-substituted phenyl ring. The second phenyl ring is attached to a trimethylammonium group, $N^+(CH_2)_3N^+(CH_3)_2$, which is further connected to a primary ammonium group, $-CH_2CH_2CH_2NH_3^+$. The counterion is $3CF_3CO_2^-$.</p>
Molecular Weight / Molecular Formula	952.83 Da; $C_{57}H_{105}N_4Cl_3$
Purity	NA
Appearance / Formulation / Solubility	Dark red solid; soluble in MeOH, DMF or DMSO.
Storage & Stability	Store desiccated at $-20^\circ C$. Protect from light, especially when in solution.
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.

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