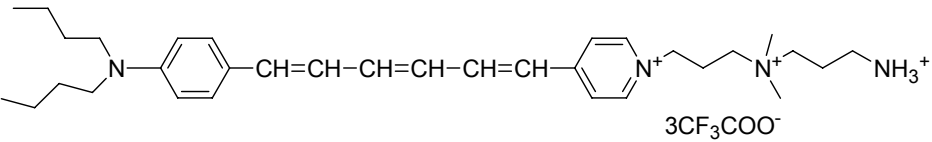


Instruction Manual

Catalog Number	PK-CA707-70039
Description	AM4-65 has been developed as a fixable, activity-dependent fluorescent nerve terminal probe. It functions similarly as SynapseRed C2 (also called FM4-64, PK-CA707-70021) does, except that AM4-65 has a fixable amine group. AM4-65 and FM4-64 have identical excitation and emission wavelengths.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in CHCl_3) = 558/734 nm; Extinction coefficient: ~46,000; Note: The excitation/emission of styryl dyes will undergo a blue shift and the dyes will become significantly more fluorescent in membranes.
Molecular Structure	
Molecular Weight / Molecular Formula	844.85 Da; $\text{C}_{39}\text{H}_{53}\text{F}_9\text{N}_4\text{O}_6$
Purity	NA
Appearance / Formulation / Solubility	Dark purple solid; soluble in water (1 mM).
Storage & Stability	Store desiccated at -20°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.