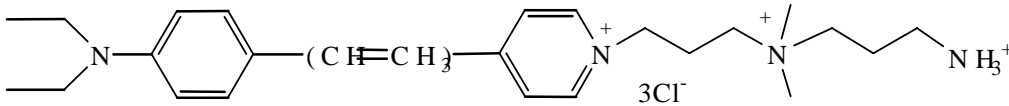


Instruction Manual

Catalog Number	PK-CA707-70025
Description	AM4-64 has been developed to be a fixable nerve terminal dye. However, its application has not been fully tested.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 543/- (too weak to detect); Extinction coefficient = 43,000
Molecular Structure	
Molecular Weight / Molecular Formula	555.5 Da; C ₂₉ H ₄₅ Cl ₃ N ₄
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.