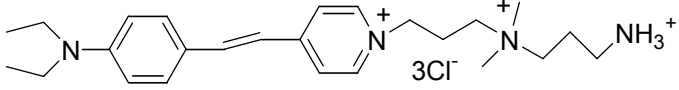


Instruction Manual

Catalog Number	PK-CA707-70036
Description	AM2-10 has been synthesized as a fixable analog of SynapseGreen C2 (FM2-10). It is spectrally and structurally similar to SynapseGreen C2 (FM2-10). However, AM2-10 has an amino group, which makes the dye fixable with the standard fixation agent. The utility of this probe is still under evaluation.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 502/625 nm; Extinction coefficient (in MeOH): ~50,000
Molecular Structure	
Molecular Weight / Molecular Formula	503.05 Da; C ₂₅ H ₃₇ Cl ₃ N ₄
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Dark red solid; soluble in water.
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.