

Sulfonated β -cyclodextrin derivative

Instruction Manual

Catalog Number	PK-CA707-70029
Description	ADVASEP-7 is a sulfonated β -cyclodextrin derivative that has been reported to reduce background fluorescence when using SynapseGreen C4 (also known as FM1-43, a trademark of Molecular Probes, Inc. ; PK-CA707-70020) to stain brain slices.
Quantity	250 mg
Excitation / Emission Maxima	NA
Molecular Structure	NA
Molecular Weight / Molecular Formula	~ 2163 Da; $C_{42}H_{70-n}O_{35}(C_4H_8SO_3Na)_n \cdot \text{hydrate}$, where n = Average Degree of Substitution = 6.5
Purity	>96%
Appearance / Formulation / Solubility	White solid; soluble in water.
Storage & Stability	Store desiccated at 4°C. Highly hygroscopic!
Applications	See Description
References	Kay, A. R., A. Alfonso, et al. "Imaging synaptic activity in intact brain and slices with FM1-43 in <i>C. elegans</i> , Lamprey and Rat." <i>Neuron</i> 24: 809-817 (1999).
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.