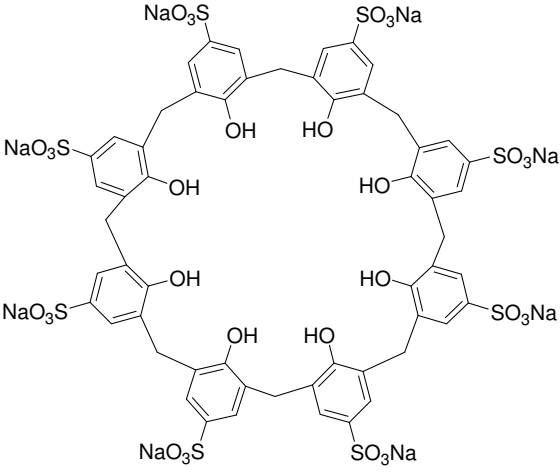


4-sulfonato calix[8]arene, sodium salt

Instruction Manual

Catalog Number	PK-CA707-70037
Description	SCAS is a quencher developed to reduce background fluorescence when using our fixable AM dyes, SynapseGreen and SynapseRed dyes. Unlike ADVASEP-7, SCAS dramatically lowers background fluorescence as soon as it is added to the preparation without the need for repeated washings.
Quantity	100 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	1665.34 Da; C ₅₆ H ₄₀ Na ₈ O ₃₂ S ₈
Purity	>95%
Appearance / Formulation / Solubility	Off-white solid. Soluble in water. Typical concentration of working solution is 0.5 mM
Storage & Stability	Store desiccated at ≤4°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.