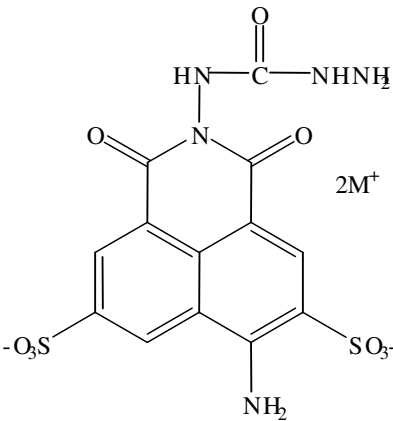


Instruction Manual

Catalog Number	PK-CA707-80015
Description	Lucifer Yellow CH lithium salt (LY CH, lithium salt) is a widely used polar molecular tracer for studying neuronal morphology. The fluorescent molecule contains a carbohydrazide that allow the molecule to be fixable with an aldehyde fixation agent.
Quantity	25 mg
Excitation / Emission Maxima	$\lambda_{ex} \setminus \lambda_{em}$: 428/536 nm; Extinction coefficient = 12,000
Molecular Structure	 <p>80015 M = Li⁺ 80016 M = K⁺</p>
Molecular Weight / Molecular Formula	457 Da; C ₁₃ H ₉ Li ₂ N ₅ O ₉ S ₂
Purity	>99% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow solid; soluble in water.
Storage & Stability	Store desiccated at $\leq 4^{\circ}\text{C}$. Protect from light, especially when in solution.
Applications	See Description
References	1) Nature 292, 17(1981) 2) Cell 14, 741(1978)
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.