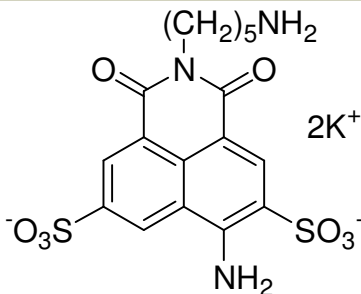


Instruction Manual

Catalog Number	PK-CA707-80018
Description	Like Lucifer Yellow CH (PK-CA707-80015 and PK-CA707-80016), Lucifer Yellow Cadaverine can be used as a fixable fluorescent tracer. In addition, the dye can be used to conjugate to carboxylic acid groups.
Quantity	25 mg
Excitation / Emission Maxima	$\lambda_{ex} \setminus \lambda_{em}$: 428/532 nm; Extinction coefficient = 11,000
Molecular Structure	
Molecular Weight / Molecular Formula	534 Da; C ₁₇ H ₁₇ K ₂ N ₃ O ₈ S ₂
Purity	>95% (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.
Applications	See Description
References	Anal. Biochem. 211, 210(1993)
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.