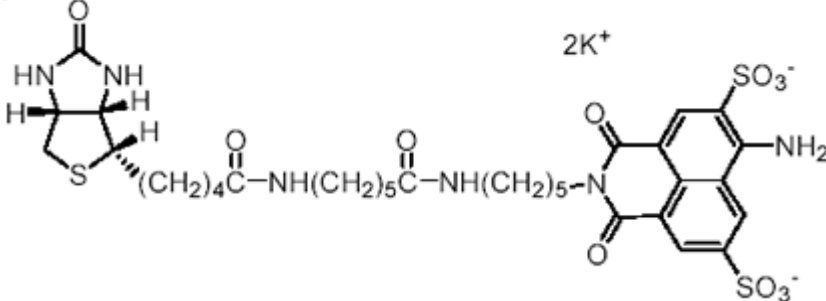


Lucifer Yellow Cadaverine Biotin-X, dipotassium salt

Tracer for studying neuronal morphology

Instruction Manual

Catalog Number	PK-CA707-80017
Description	This fluorescent biotin derivative is a very useful tracer for studying neuronal morphology. The biotin portion of the molecule allows researchers to amplify the fluorescent signal using standard biotin-avidin or biotin-streptavidin biochemistry. The staining can tolerate conditions of aldehyde fixation and sectioning, and washing with detergent.
Quantity	10 mg
Excitation / Emission Maxima	λ_{ex} \ λ_{em} : 428/532 nm; Extinction coefficient = 11,000
Molecular Structure	
Molecular Weight / Molecular Formula	873 Da; C ₃₃ H ₄₂ K ₂ N ₆ O ₁₁ S ₃
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow solid; soluble in water.
Storage & Stability	Store desiccated at -20°C. Protect from light, especially when in solution.
Applications	See Description
References	1) J. Neurosci. Meth. 53, 23(1994) 2) J. Neurosci. Meth. 46, 59(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.