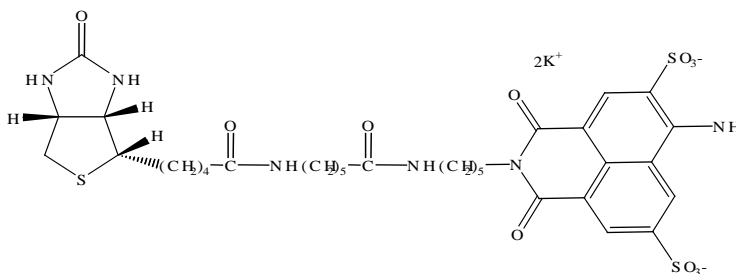


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Lucifer Yellow Cadaverine Biotin-X, dipotaasium salt**CATALOG #:** 80017**MOLECULAR INFORMATION:** C₃₃H₄₂K₂N₆O₁₁S₃
MW: 873.**PROPERTIES:**

Color & Form	Yellow solid
Purity	≥ 95% by HPLC
Solubility	Soluble in water
Absorption/Emission	428/532 nm
Extinction Coefficient	11,000

STORAGE AND HANDLING: Store desiccated at ≤ 4°C. Protect from light.**APPLICATION:** 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.Ref.: 1) *J. Neurosci. Meth.* **53**, 23(1994); 2) *J. Neurosci. Meth.* **46**, 59(1993)**TOXICITY:** Unknown.

FIRST AID:	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.
-------------------	---

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.
