

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:	Lucifer Yellow Cadaverine	
CATALOG #:	80018	
MOLECULAR INFORMATION:	C ₁₇ H ₁₇ K ₂ N ₃ O ₈ S ₂ MW: 534	(CH ₂) ₅ NH ₂ O N O 2K ⁺ SO ₃ ⁻ NH ₂

PROPERTIES: Color & Form Purity Solubility Absorption/Emission Extinction Coefficient	Yellow solid ≥ 95% by HPLC Soluble in water 428/532 nm 11,000
STORAGE AND HANDLING:	Store desiccated at 4°C. Protect from light.
APPLICATION:	Like Lucifer Yellow CH (80015 and 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.
	Ref.: Anal. Biochem. 211, 210(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.*