

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:	Biocytin (ε-biotinoyl –L-lysine)
CATALOG #:	90055
MOLECULAR INFORMATION:	C ₁₆ H ₂₈ N ₄ O ₄ S MW: 372.48
	$HN^{\prime} \qquad NH \\ H \qquad H \qquad O \qquad NH_{3}+ \\ H \qquad H \qquad H \qquad H \qquad H^{\prime} \qquad $
PROPERTIES: Color & Form Purity Solubility	White solid ≥ 95% by TLC Soluble in water or DMSO
STORAGE AND HANDLING:	Store desiccated at \leq 4 °C
APPLICATION:	Biocytin is a very useful cellular tracer that can be introduced into cells by microinjection.
	Ref.: 1) Anat Embryol 199, 149 (1998); 2)Brain Res 782, 113 (1998); 3)Brain Res Devel Brain Res 106, 79 (1998); 4) Exp Brain Res 119, 375 (1998).
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.
Discloiment Magnials for	seek medical advice immediately. n Biotium are sold for research use only, and are not intended for food, drug, household, or

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.