

Instruction Manual

Catalog Number	PK-CA707-90055
Description	Biocytin is a very useful cellular tracer that can be introduced into cells by microinjection.
Quantity	100 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	372.48 Da; C ₁₆ H ₂₈ N ₄ O ₄ S
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	White solid; soluble in water or DMSO.
Storage & Stability	Store desiccated at ≤4°C.
Applications	See Description
References	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)
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.