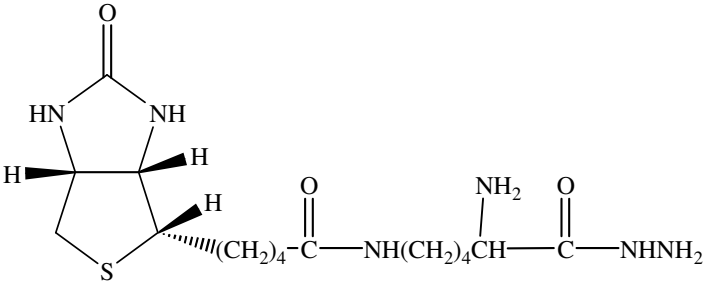


### Instruction Manual

Catalog Number	PK-CA707-90060
Description	Biocytin-Hydrazide is an aldehyde-fixable cellular tracer that can be introduced into cells by microinjection. Biocytin hydrazide can also be used to label glycoproteins or other carbohydrate molecules.
Quantity	25 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	386.51 Da; C <sub>16</sub> H <sub>30</sub> N <sub>6</sub> O <sub>3</sub> S
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	White solid; soluble in DMSO.
Storage & Stability	Store at -20°C.
Applications	See Description
References	<ol style="list-style-type: none"> <li>1) Glycoconjugate J. 9, 21(1992)</li> <li>2) Anal Biochem 171, 1(1988)</li> <li>3) Anal. Biochem. 170, 271(1988)</li> <li>4) Biochem. Biophys. Res. Comm. 136, 80(1986)</li> </ol>
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.