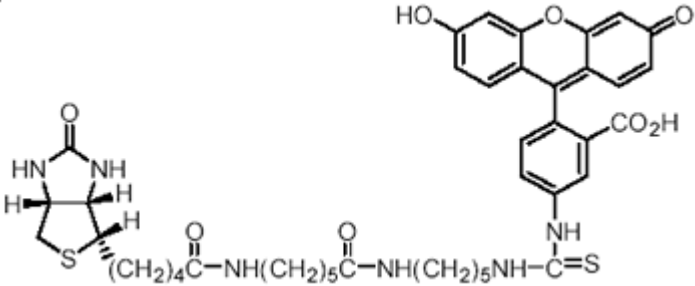


5-((N-(5-(N-(6-(biotinoyl)amino)hexanoyl)amino)pentyl)thioureidyl) fluorescein

Instruction Manual

Catalog Number	PK-CA707-80019
Description	Fluorescein biotin can be used to detect and quantify biotin binding sites and the degree of biotinylation of proteins. Also see biotin-rhodamine 110 (Cat.No. PK-CA707-80022) and biotin-4-fluorescein (Cat.No. PK-CA707-90062).
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \setminus \lambda_{em}$: 494/518 nm; Extinction coefficient = 75,000
Molecular Structure	
Molecular Weight / Molecular Formula	831 Da; C ₄₂ H ₅₀ N ₆ O ₈ S ₂
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow solid; soluble in DMF or DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	See Description
References	1) Biochim Biophys Acta 1381, 203 (1998) 2) Clin.Chem. 40, 2112(1994)
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.