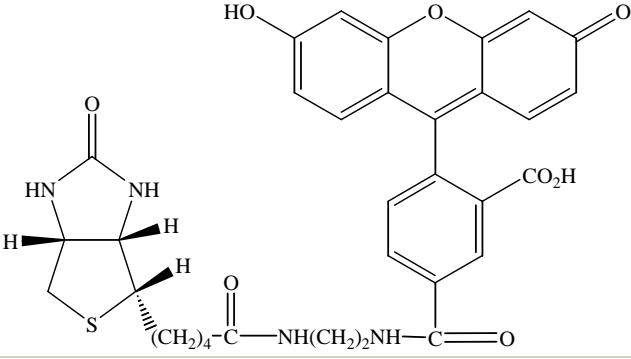


### Instruction Manual

Catalog Number	PK-CA707-90062
Description	Biotin-4-fluorescein has been reported to be a superior fluorescent biotin derivative for accurate measurement of avidin and streptavidin in crude biofluids. The biotin binding to avidin and streptavidin is high affinity, fast and noncooperative. Also see Fluorescein biotin (PK-CA707-80019). Please see also biotin-rhodamine 110 (PK-CA707-80022) for superior photostability.
Quantity	10 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em} = 494/523$ nm
Molecular Structure	
Molecular Weight / Molecular Formula	644.7 Da; C <sub>33</sub> H <sub>32</sub> N <sub>4</sub> O <sub>8</sub> S
Purity	>99% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow solid; soluble in DMSO or buffers with pH>7.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	See Description
References	1) PNAS 102, 24, pp 8507 (2005) 2) Biochim Biophys Acta 1427, 33 (1999)
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.