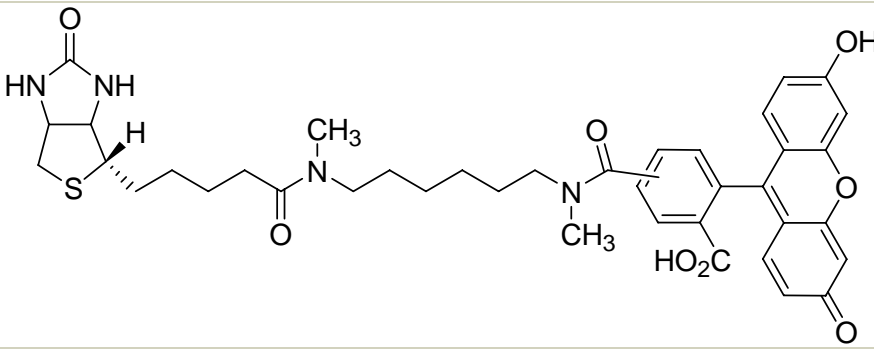


Instruction Manual

Catalog Number	PK-CA707-51024
Description	Flubi-2 is the hydrolyzed product of the site-directed fluorogenic probe Flubida-2 (Cat.No. PK-CA707-51022) and can be used as a standard for calibration. The dye can also be microinjected into cells to probe the local pH where avidin-chimera proteins are located.
Quantity	5 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	728.85 Da; C ₃₉ H ₄₄ N ₄ O ₈ S
Purity	NA
Appearance / Formulation / Solubility	Yellow solid; soluble in DMSO.
Storage & Stability	Store at -20°C and protect from light.
Applications	Calibration standard for Flubida-2
References	Chemistry & Biology 7, 197(2000)
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.