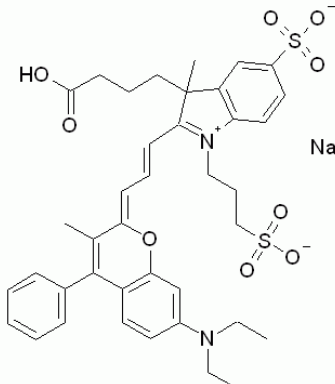
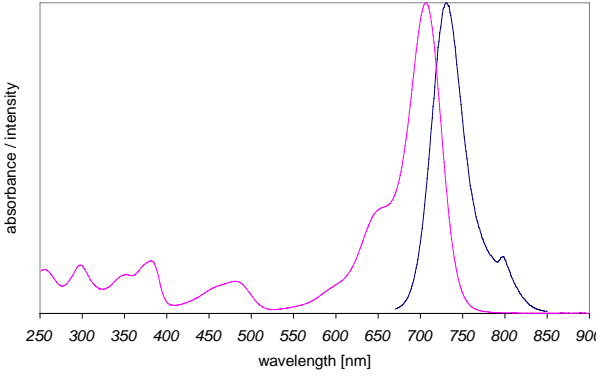


Instruction Manual

Catalog Number	PK-PF700P-0-01	PromoFluor-700P, carboxylic acid	1 mg
	PK-PF700P-0-05	PromoFluor-700P, carboxylic acid	5 mg
	PK-PF700P-1-01	PromoFluor-700P, NHS ester	1 mg
	PK-PF700P-1-01A	PromoFluor-700P, NHS ester	5 x 0.2 mg
	PK-PF700P-1-05	PromoFluor-700P, NHS ester	5 mg
	PK-PF700P-2-01	PromoFluor-700P, amino-modified	1 mg
	PK-PF700P-3-01	PromoFluor-700P, maleimide	1 mg
	PK-PF700P-3-05	PromoFluor-700P, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	706 nm / 731 nm (in ethanol)		
Molar Absorbance	140.000 M ⁻¹ cm ⁻¹		
Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	770.90 g·mol ⁻¹	C ₃₉ H ₄₃ N ₂ O ₉ S ₂ Na
	NHS-ester	867.97 g·mol ⁻¹	C ₄₃ H ₄₆ N ₃ O ₁₁ S ₂ Na
	Maleimide	893.03 g·mol ⁻¹	C ₄₅ H ₄₉ N ₄ O ₁₀ S ₂ Na
	Amino-derivative	791.01 g·mol ⁻¹	C ₄₁ H ₅₀ N ₄ O ₈ S
Formulation	Solid		
Solubility	Soluble in methanol, DMF or DMSO (up to 10 mg/ml).		
Storage & Stability	Store at -20°C and protect from light and humidity.		
Applications	Fluorescent labeling of biomolecules (e.g. proteins/peptides, nucleic acids). Labeling protocols are available at www.promkine.info .		

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.