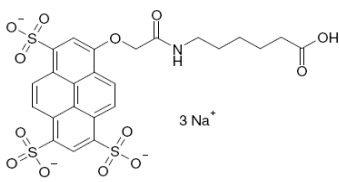
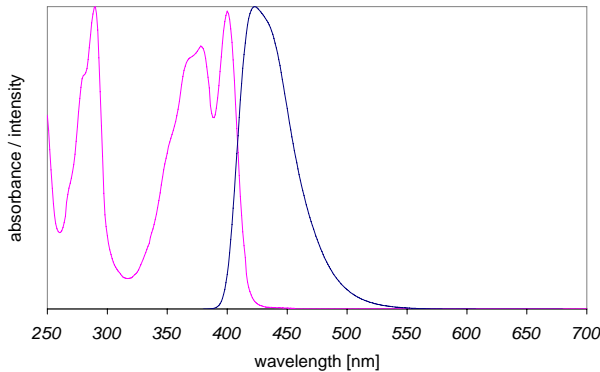


Instruction Manual

Catalog Number	PK-PF405-0-01	PromoFluor-405, carboxylic acid	1 mg
	PK-PF405-0-05	PromoFluor-405, carboxylic acid	5 mg
	PK-PF405-1-01	PromoFluor-405, NHS-ester	1 mg
	PK-PF405-1-01A	PromoFluor-405, NHS-ester	5 x 0.2 mg
	PK-PF405-1-05	PromoFluor-405, NHS-ester	5 mg
	PK-PF405-2-01	PromoFluor-405, amino-modified	1 mg
	PK-PF405-3-01	PromoFluor-405, maleimide	1 mg
	PK-PF405-3-05	PromoFluor-405, maleimide	5 mg
Structure & Spectra	Chemical Structure 		Absorption and Emission Spectra 
	Absorption / Emission max. 400 nm / 420 nm (in ethanol); 400 nm / 423 nm (in PBS)		
Molar Absorbance	32.000 M ⁻¹ cm ⁻¹		
Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	695.59 g·mol ⁻¹	C ₂₄ H ₂₀ NO ₁₃ S ₃ Na ₃
	NHS-ester	792.66 g·mol ⁻¹	C ₂₈ H ₂₃ N ₂ O ₁₅ S ₃ Na ₃
	Maleimide	817.72 g·mol ⁻¹	C ₃₀ H ₂₆ N ₃ O ₁₄ S ₃ Na ₃
Amino-derivative	715.69 g·mol ⁻¹	C ₂₆ H ₂₇ N ₃ O ₁₂ S ₃ Na ₂	
Formulation	Solid		
Solubility	Soluble in water or methanol (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, antibodies, nucleic acids). Labeling protocols are available at www.promkine.info .		

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