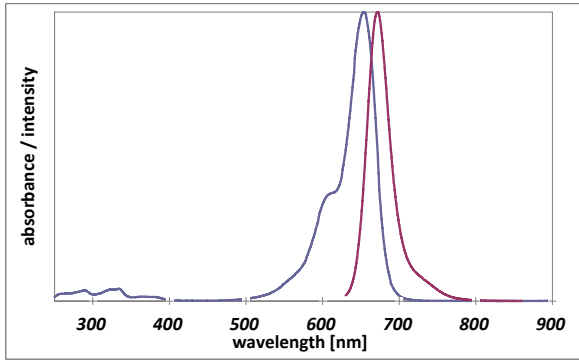


Instruction Manual

Catalog Number	PK-PF647P-0-01	PromoFluor-647P, carboxylic acid	1 mg
	PK-PF647P-0-05	PromoFluor-647P, carboxylic acid	5 mg
	PK-PF647P-1-01	PromoFluor-647P, NHS ester	1 mg
	PK-PF647P-1-01A	PromoFluor-647P, NHS ester	5 x 0.2 mg
	PK-PF647P-1-05	PromoFluor-647P, NHS ester	5 mg
	PK-PF647P-2-01	PromoFluor-647P, amino-modified	1 mg
	PK-PF647P-3-01	PromoFluor-647P, maleimide	1 mg
	PK-PF647P-3-05	PromoFluor-647P, maleimide	5 mg
Structure & Spectra	Chemical Structure		Absorption and Emission Spectra
	Structure available on request.		
Absorption / Emission max.	654 nm / 672 nm (in PBS); 657 nm / 678 nm (in ethanol)		
Molar Absorbance	250.000 M ⁻¹ cm ⁻¹		
Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	969.03 g·mol ⁻¹	C ₃₈ H ₄₇ N ₂ O ₁₅ S ₄ * 3 Na
	NHS-ester	1066.1 g·mol ⁻¹	C ₄₂ H ₅₀ N ₃ O ₁₇ S ₄ * 3 Na
	Maleimide	1091.16 g·mol ⁻¹	C ₄₄ H ₅₃ N ₄ O ₁₆ S ₄ * 3 Na
	Amino-derivative	989.13 g·mol ⁻¹	C ₄₀ H ₅₄ N ₄ O ₁₄ S ₄ * 2 Na
Formulation	Solid		
Solubility	Soluble in water, 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, antibodies, nucleic acids). Labeling protocols are available at www.promkine.info .		

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