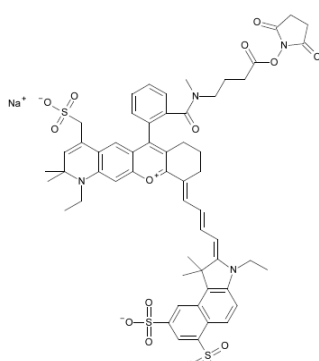
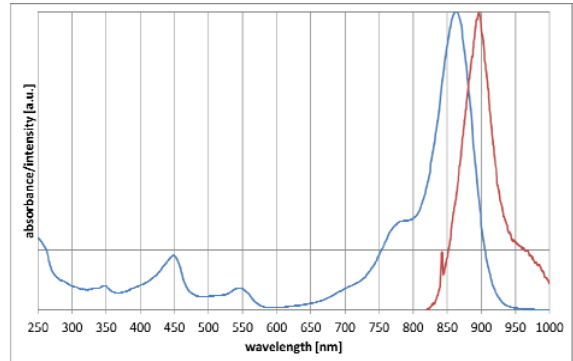


Instruction Manual

Catalog Number	PK-PF865-0-01	PromoFluor-865, carboxylic acid	1 mg
	PK-PF865-0-05	PromoFluor-865, carboxylic acid	5 mg
	PK-PF865-1-01	PromoFluor-865, NHS ester	1 mg
	PK-PF865-1-01A	PromoFluor-865, NHS ester	5 x 0.2 mg
	PK-PF865-1-05	PromoFluor-865, NHS ester	5 mg
	PK-PF865-2-01	PromoFluor-865, amino-modified	1 mg
	PK-PF865-3-01	PromoFluor-865, maleimide	1 mg
	PK-PF865-3-05	PromoFluor-865, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max. 863 nm / 896 nm (in ethanol)

Molar Absorbance 190.000 M⁻¹ cm⁻¹

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	1084.19 g·mol ⁻¹	C ₅₃ H ₅₅ N ₃ O ₁₃ S ₃ *2Na
	NHS-ester	1181.26 g·mol ⁻¹	C ₅₇ H ₅₈ N ₄ O ₁₅ S ₃ *2Na
	Maleimide	1104.29 g·mol ⁻¹	C ₅₅ H ₆₂ N ₅ O ₁₂ S ₃ Na
	Amino-derivative	1206.32 g·mol ⁻¹	C ₅₉ H ₆₁ N ₅ O ₁₄ S ₃ *2Na

Formulation Solid

Solubility Soluble in water, methanol, DMF or DMSO. Slightly soluble in water.

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.promokine.info.

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