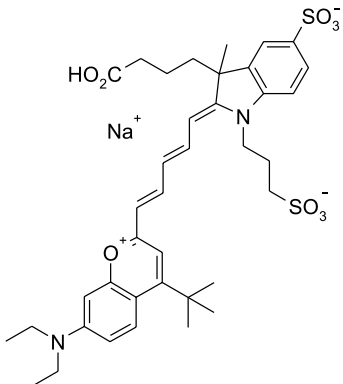
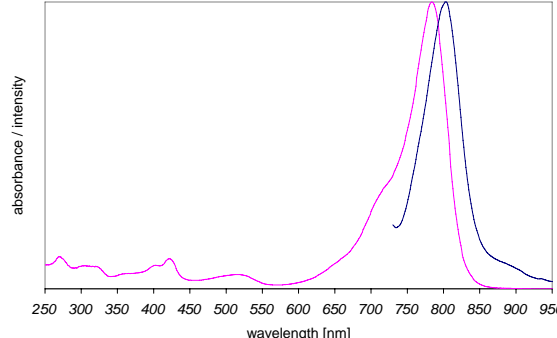


Instruction Manual

Catalog Number	PK-PF780-0-01	PromoFluor-780, carboxylic acid	1 mg
	PK-PF780-0-05	PromoFluor-780, carboxylic acid	5 mg
	PK-PF780-1-01	PromoFluor-780, NHS ester	1 mg
	PK-PF780-1-01A	PromoFluor-780, NHS ester	5 x 0.2 mg
	PK-PF780-1-05	PromoFluor-780, NHS ester	5 mg
	PK-PF780-2-01	PromoFluor-780, amino-modified	1 mg
	PK-PF780-3-01	PromoFluor-780, maleimide	1 mg
	PK-PF780-3-05	PromoFluor-780, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	783 nm / 800 nm (in ethanol)
----------------------------	------------------------------

Molar Absorbance	170.000 M ⁻¹ cm ⁻¹
------------------	--

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	762.92 g·mol ⁻¹	C ₃₈ H ₄₇ N ₂ O ₉ S ₂ Na
	NHS-ester	860.00 g·mol ⁻¹	C ₄₂ H ₅₀ N ₃ O ₁₁ S ₂ Na
	Maleimide	885.05 g·mol ⁻¹	C ₄₄ H ₅₃ N ₄ O ₁₀ S ₂ Na
	Amino-derivative	783.03 g·mol ⁻¹	C ₄₀ H ₅₄ N ₄ O ₈ S ₂

Formulation	Solid
-------------	-------

Solubility	Soluble in water, methanol, DMF and 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.promokine.info .
--------------	--

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