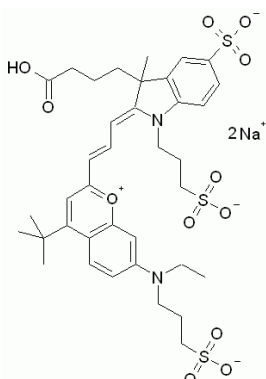
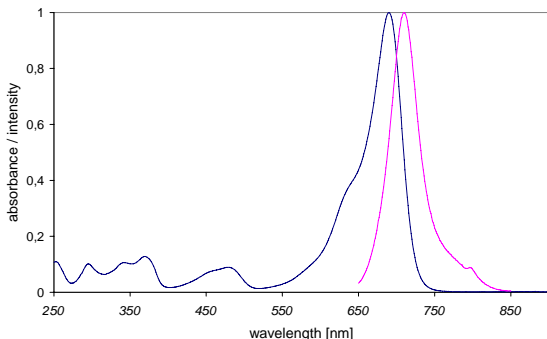


Instruction Manual

Catalog Number	PK-PF680-0-01	PromoFluor-680, carboxylic acid	1 mg
	PK-PF680-0-05	PromoFluor-680, carboxylic acid	5 mg
	PK-PF680-1-01	PromoFluor-680, NHS ester	1 mg
	PK-PF680-1-01A	PromoFluor-680, NHS ester	5 x 0.2 mg
	PK-PF680-1-05	PromoFluor-680, NHS ester	5 mg
	PK-PF680-2-01	PromoFluor-680, amino-modified	1 mg
	PK-PF680-3-01	PromoFluor-680, maleimide	1 mg
	PK-PF680-3-05	PromoFluor-680, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	690 nm / 709 nm (in ethanol)
----------------------------	------------------------------

Molar Absorbance	140.000 M ⁻¹ cm ⁻¹
------------------	--

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	852.96 g·mol ⁻¹	C ₃₇ H ₄₆ N ₂ O ₁₂ S ₃ Na ₂
	NHS-ester	950.03 g·mol ⁻¹	C ₄₁ H ₄₉ N ₃ O ₁₄ S ₃ Na ₂
	Maleimide	975.08 g·mol ⁻¹	C ₄₃ H ₅₂ N ₄ O ₁₃ S ₃ Na ₂
	Amino-derivative	873.05 g·mol ⁻¹	C ₃₉ H ₅₃ N ₄ O ₁₁ 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.promokine.info .
--------------	--

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