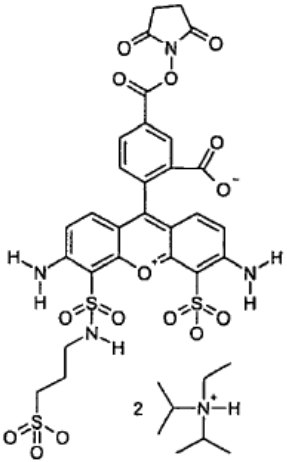
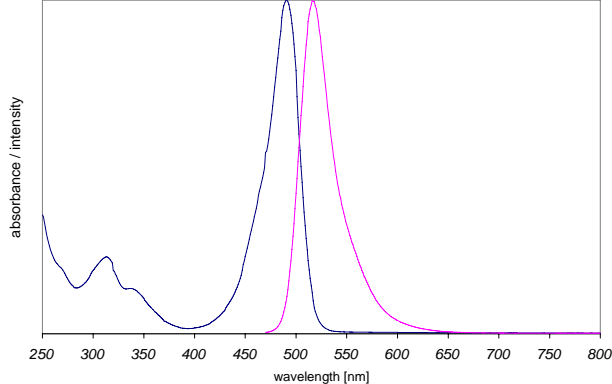


Instruction Manual

Catalog Number	PK-PF488P-0-01	PromoFluor-488 Premium, carboxylic acid	1 mg
	PK-PF488P-0-05	PromoFluor-488 Premium, carboxylic acid	5 mg
	PK-PF488P-1-01	PromoFluor-488 Premium, NHS ester	1 mg
	PK-PF488P-1-01A	PromoFluor-488 Premium, NHS ester	5 x 0.2 mg
	PK-PF488P-1-05	PromoFluor-488 Premium, NHS ester	5 mg
	PK-PF488P-2-01	PromoFluor-488 Premium amino-modified	1 mg
	PK-PF488P-3-01	PromoFluor-488 Premium, maleimide	1 mg
	PK-PF488P-3-05	PromoFluor-488 Premium, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	490 nm / 516 nm (in PBS)
----------------------------	--------------------------

Molar Absorbance	73.000 M ⁻¹ cm ⁻¹
------------------	---

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	769.67 g·mol ⁻¹	C ₂₄ H ₂₂ N ₃ O ₁₃ S ₃ x CF ₃ COOH
	NHS-ester	1011.20 g·mol ⁻¹	C ₂₈ H ₂₂ N ₄ O ₁₅ S ₃ x 2C ₈ H ₂₀ N
	Maleimide	799.74 g·mol ⁻¹	C ₃₀ H ₂₆ N ₅ O ₁₄ S ₃ x Na
	Amino-derivative	925.78 g·mol ⁻¹	C ₂₆ H ₂₉ N ₅ O ₁₂ S ₃ x CF ₃ COO

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.