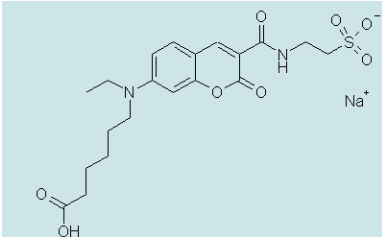
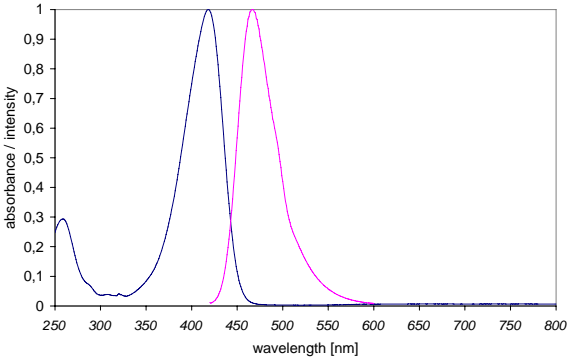


Instruction Manual

Catalog Number	PK-PF415-0-01	PromoFluor-415, carboxylic acid	1 mg
	PK-PF415-0-05	PromoFluor-415, carboxylic acid	5 mg
	PK-PF415-1-01	PromoFluor-415, NHS ester	1 mg
	PK-PF415-1-01A	PromoFluor-415, NHS ester	5 x 0.2 mg
	PK-PF415-1-05	PromoFluor-415, NHS ester	5 mg
	PK-PF415-2-01	PromoFluor-415, amino-modified	1 mg
	PK-PF415-3-01	PromoFluor-415, maleimide	1 mg
	PK-PF415-3-05	PromoFluor-415, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	418 nm / 467 nm (in ethanol)
----------------------------	------------------------------

Molar Absorbance	34.000 M ⁻¹ cm ⁻¹
------------------	---

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	476.48 g·mol ⁻¹	C ₂₀ H ₂₅ N ₂ O ₈ SNa
	NHS-ester	573.56 g·mol ⁻¹	C ₂₄ H ₂₈ N ₃ O ₁₀ SNa
	Maleimide	598.61 g·mol ⁻¹	C ₂₆ H ₃₁ N ₄ O ₉ SNa
	Amino-derivative	496.59 g·mol ⁻¹	C ₂₂ H ₃₂ N ₄ O ₇ S

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

Solubility	Soluble in water, methanol, ethanol, 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.promokine.info .
--------------	--

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