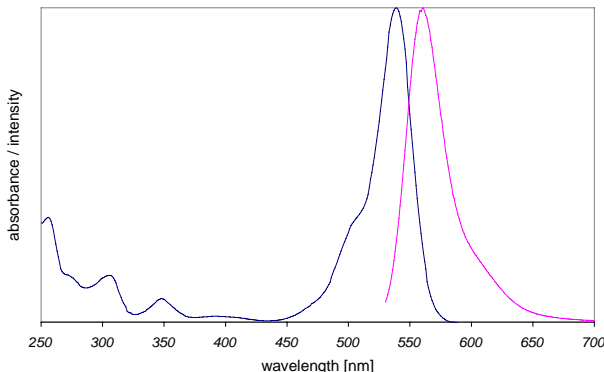


Instruction Manual

Catalog Number	PK-PF532-0-01	PromoFluor-532, carboxylic acid	1 mg
	PK-PF532-0-05	PromoFluor-532, carboxylic acid	5 mg
	PK-PF532-1-01	PromoFluor-532, NHS ester	1 mg
	PK-PF532-1-01A	PromoFluor-532, NHS ester	5 x 0.2 mg
	PK-PF532-1-05	PromoFluor-532, NHS ester	5 mg
	PK-PF532-2-01	PromoFluor-532, amino-modified	1 mg
	PK-PF532-3-01	PromoFluor-532, maleimide	1 mg
	PK-PF532-3-05	PromoFluor-532, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
	Structure available on request.	

Absorption / Emission max.	539 nm / 561 nm (in ethanol)
----------------------------	------------------------------

Molar Absorbance	100.000 M ⁻¹ cm ⁻¹
------------------	--

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	695.75 g·mol ⁻¹	C ₃₁ H ₄₃ N ₃ O ₁₀ S ₂ *Na
	NHS-ester	792.82 g·mol ⁻¹	C ₃₅ H ₃₇ N ₄ O ₁₂ S ₂ *Na
	Maleimide	817.88 g·mol ⁻¹	C ₃₇ H ₄₀ N ₅ O ₁₁ S ₂ *Na
	Amino-derivative	715.78 g·mol ⁻¹	C ₃₃ H ₄₁ N ₅ O ₉ S ₂

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

Solubility	Soluble in water, methanol, ethanol, DMF, 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.promkine.info .
--------------	--

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