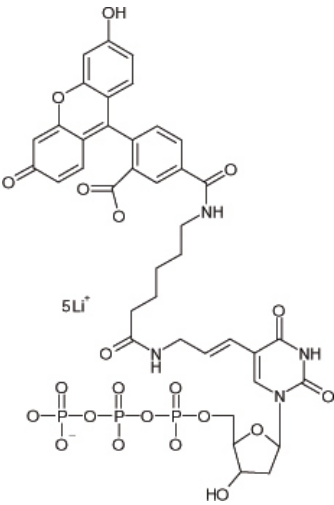
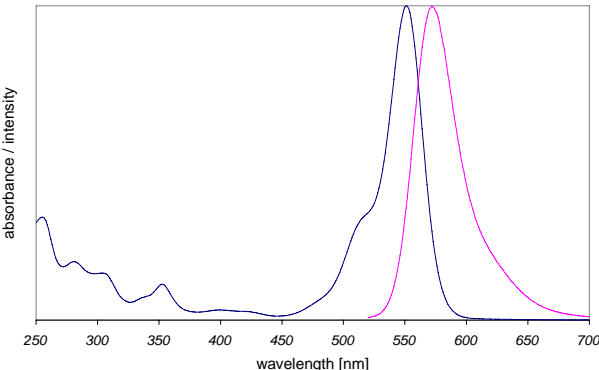


Instruction Manual

Catalog Number	PK-PF546-0-01	PromoFluor-546, carboxylic acid	1 mg
	PK-PF546-0-05	PromoFluor-546, carboxylic acid	5 mg
	PK-PF546-1-01	PromoFluor-546, NHS ester	1 mg
	PK-PF546-1-01A	PromoFluor-546, NHS ester	5 x 0.2 mg
	PK-PF546-1-05	PromoFluor-546, NHS ester	5 mg
	PK-PF546-2-01	PromoFluor-546, amino-modified	1 mg
	PK-PF546-3-01	PromoFluor-546, maleimide	1 mg
	PK-PF546-3-05	PromoFluor-546, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	551 nm / 572 nm (in ethanol)
----------------------------	------------------------------

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

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	580.08 g·mol ⁻¹	C ₃₁ H ₃₄ N ₃ O ₆ Cl
	NHS-ester	677.16 g·mol ⁻¹	C ₃₅ H ₃₇ N ₄ O ₈ Cl
	Maleimide	665.74 g·mol ⁻¹	C ₃₇ H ₃₉ N ₅ O ₇
	Amino-derivative	622.17 g·mol ⁻¹	C ₃₃ H ₄₀ N ₅ O ₅ Cl

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

Solubility	Soluble in 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.