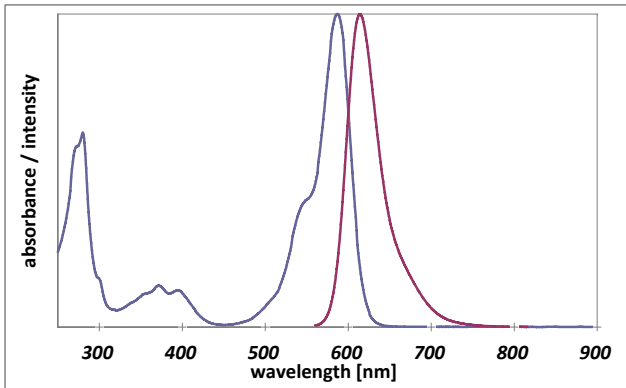


Instruction Manual

Catalog Number	PK-PF594-0-01	PromoFluor-594, carboxylic acid	1 mg
	PK-PF594-0-05	PromoFluor-594, carboxylic acid	5 mg
	PK-PF594-1-01	PromoFluor-594, NHS ester	1 mg
	PK-PF594-1-01A	PromoFluor-594, NHS ester	5 x 0.2 mg
	PK-PF594-1-05	PromoFluor-594, NHS ester	5 mg
	PK-PF594-2-01	PromoFluor-594, amino-modified	1 mg
	PK-PF594-3-01	PromoFluor-594, maleimide	1 mg
	PK-PF594-3-05	PromoFluor-594, maleimide	5 mg

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

Absorption / Emission max.	587 nm / 614 nm (in PBS); 594 nm / 615 nm (in ethanol)
----------------------------	--

Molar Absorbance	92.000 M ⁻¹ cm ⁻¹
------------------	---

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	937.06 g·mol ⁻¹	C ₃₉ H ₄₄ N ₄ O ₁₅ S ₄
	NHS-ester	1078.10 g·mol ⁻¹	C ₄₃ H ₄₅ N ₅ O ₁₇ S ₄ * 2Na
	Maleimide	1059.19 g·mol ⁻¹	C ₄₅ H ₅₀ N ₆ O ₁₆ S ₄
	Amino-derivative	979.14 g·mol ⁻¹	C ₄₁ H ₅₀ N ₆ O ₁₄ S ₄

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

Solubility	Soluble in methanol, 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.promkine.info .
--------------	--

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