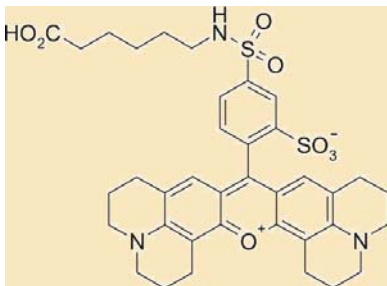
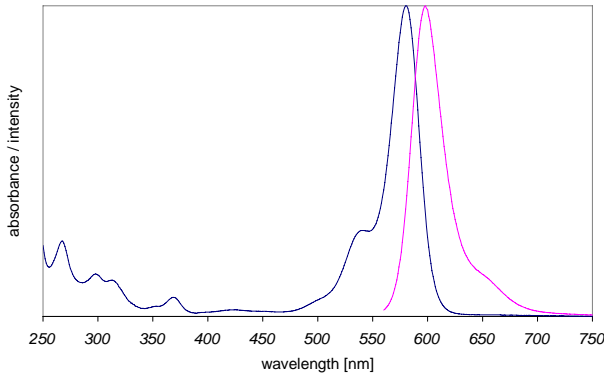


Instruction Manual

Catalog Number	PK-PF568-0-01	PromoFluor-568, carboxylic acid	1 mg
	PK-PF568-0-05	PromoFluor-568, carboxylic acid	5 mg
	PK-PF568-1-01	PromoFluor-568, NHS ester	1 mg
	PK-PF568-1-01A	PromoFluor-568, NHS ester	5 x 0.2 mg
	PK-PF568-1-05	PromoFluor-568, NHS ester	5 mg
	PK-PF568-2-01	PromoFluor-568, amino-modified	1 mg
	PK-PF568-3-01	PromoFluor-568, maleimide	1 mg
	PK-PF568-3-05	PromoFluor-568, maleimide	5 mg

Structure & Spectra	Chemical Structure	Absorption and Emission Spectra
		

Absorption / Emission max.	580 nm / 598 nm (in ethanol)		
Molar Absorbance	120.000 M ⁻¹ cm ⁻¹		
Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	719.88 g·mol ⁻¹	C ₃₇ H ₄₁ N ₃ O ₈ S ₂
	NHS-ester	816.96 g·mol ⁻¹	C ₄₁ H ₄₄ N ₄ O ₁₀ S ₂
	Maleimide	842.01 g·mol ⁻¹	C ₄₃ H ₄₇ N ₅ O ₉ S ₂
	Amino-derivative	798.42 g·mol ⁻¹	C ₃₉ H ₄₈ N ₅ O ₇ S ₂ Cl
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.promokine.info .		

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