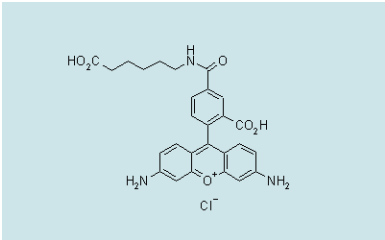
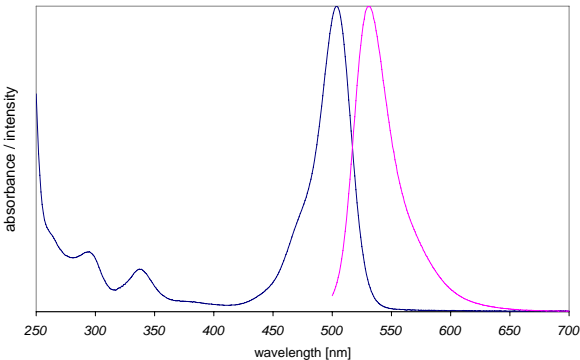


Instruction Manual

Catalog Number	PK-PF505-6-01	PromoFluor-505, biotin	1 mg
	PK-PF505-7-01	PromoFluor-505, phalloidin	10 nmole/300 units
	PK-PF505-8-100	PromoFluor-505-aadUTP	100 µl (1 mM)

Structure & Spectra	Chemical Structure		Absorption and Emission Spectra
			
Absorption / Emission max.	505 nm / 530 nm (in ethanol)		
Molar Absorbance	80.000 M ⁻¹ cm ⁻¹		
Molecular Data	Modification	Molecular Weight	Molecular Formula
	Phalloidin	1257.40 g·mol ⁻¹	C ₆₂ H ₇₂ N ₁₂ O ₁₅ S
	dUTP	1016.46 g·mol ⁻¹	C ₃₉ H ₃₉ N ₆ O ₁₉ P ₃ *4Li
	Biotin	644.10 g·mol ⁻¹	C ₄₃ H ₅₃ N ₇ O ₉ S
Formulation	<ul style="list-style-type: none"> - PromoFluor-Biotin Conjugate: lyophilized - PromoFluor-Phalloidin Conjugate: lyophilized (10 nmole, 300 units) - PromoFluor-dUTP Conjugate: 1 mM in aqueous solution (100 nmole) 		
Solubility	Phalloidin conjugates can be solubilized in methanol, biotin conjugates in methanol or DMF.		
Storage & Stability	Store at -20°C in the dark and protect solid and lyophilized conjugates from humidity.		

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