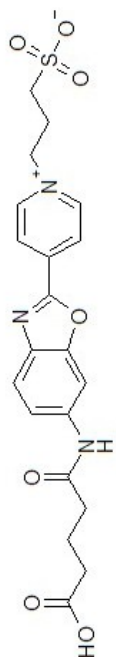
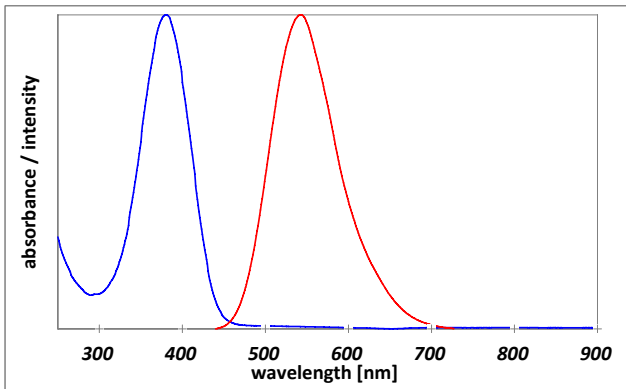


Instruction Manual

Catalog Number	PK-PF375LSS-6-01	PromoFluor-375LSS, biotin	1 mg
	PK-PF375LSS-7-01	PromoFluor-375LSS, phalloidin	10 nmole/300 units
	PK-PF375LSS-8-100	PromoFluor-375LSS-aadUTP	100 µl (1 mM)

Structure & Spectra	Chemical Structure		Absorption and Emission Spectra	
				
Absorption / Emission max.	376 nm / 543 nm (in ethanol)			
Molar Absorbance	16.000 M ⁻¹ cm ⁻¹			
Molecular Data	Modification	Molecular Weight	Molecular Formula	
	Phalloidin	1217.35 g·mol ⁻¹	C ₅₅ H ₆₈ N ₁₂ O ₁₆ S	
	dUTP	976.41 g·mol ⁻¹	C ₃₂ H ₃₅ N ₆ O ₂₀ P ₃ S * 4Li	
	Biotin	803.96 g·mol ⁻¹	C ₃₆ H ₄₉ N ₇ O ₁₀ S ₂	
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.