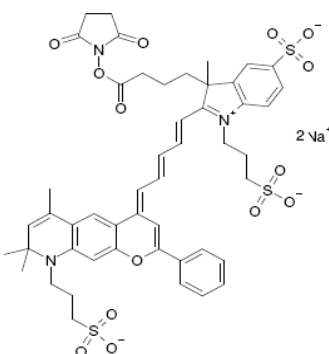
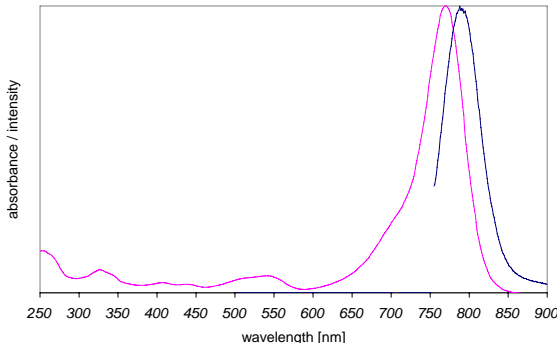


## Instruction Manual

Catalog Number	PK-PF770-0-01	PromoFluor-770, carboxylic acid	1 mg
	PK-PF770-0-05	PromoFluor-770, carboxylic acid	5 mg
	PK-PF770-1-01	PromoFluor-770, NHS ester	1 mg
	PK-PF770-1-01A	PromoFluor-770, NHS ester	5 x 0.2 mg
	PK-PF770-1-05	PromoFluor-770, NHS ester	5 mg
	PK-PF770-2-01	PromoFluor-770, amino-modified	1 mg
	PK-PF770-3-01	PromoFluor-770, maleimide	1 mg
	PK-PF770-3-05	PromoFluor-770, maleimide	5 mg
	Structure & Spectra	<b>Chemical Structure</b>	
			
Absorption / Emission max.	769 nm / 796 nm (in ethanol)		
Molar Absorbance	220.000 M <sup>-1</sup> cm <sup>-1</sup>		
Molecular Data	<b>Modification</b>	<b>Molecular Weight</b>	<b>Molecular Formula</b>
	Carboxylic acid	951.06 g·mol <sup>-1</sup>	C <sub>45</sub> H <sub>48</sub> N <sub>2</sub> O <sub>12</sub> S <sub>3</sub> Na <sub>2</sub>
	NHS-ester	1048.14 g·mol <sup>-1</sup>	C <sub>49</sub> H <sub>51</sub> N <sub>3</sub> O <sub>14</sub> S <sub>3</sub> Na <sub>2</sub>
	Maleimide	1073.19 g·mol <sup>-1</sup>	C <sub>51</sub> H <sub>54</sub> N <sub>4</sub> O <sub>13</sub> S <sub>3</sub> Na <sub>2</sub>
	Amino-derivative	971.16 g·mol <sup>-1</sup>	C <sub>47</sub> H <sub>55</sub> N <sub>4</sub> O <sub>11</sub> S <sub>3</sub> Na
Formulation	Solid		
Solubility	Soluble in water or methanol (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 <a href="http://www.promokine.info">www.promokine.info</a> .		

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