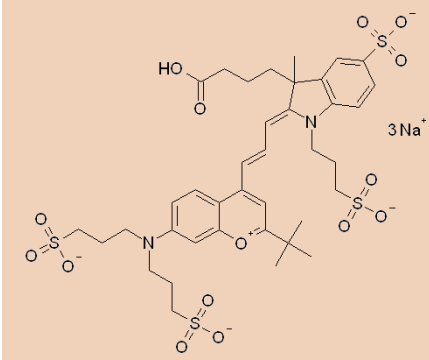
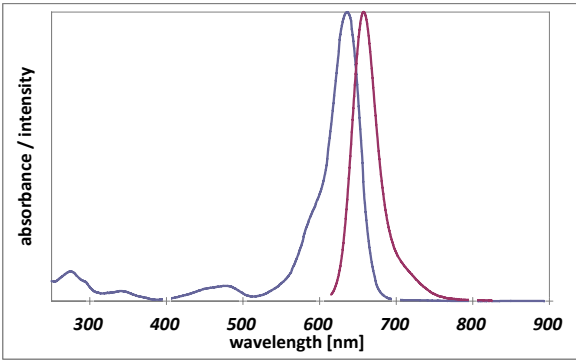


## Instruction Manual

Catalog Number	PK-PF633P-0-01	PromoFluor-633P, carboxylic acid	1 mg
	PK-PF633P-0-05	PromoFluor-633P, carboxylic acid	5 mg
	PK-PF633P-1-01	PromoFluor-633P, NHS ester	1 mg
	PK-PF633P-1-01A	PromoFluor-633P, NHS ester	5 x 0.2 mg
	PK-PF633P-1-05	PromoFluor-633P, NHS ester	5 mg
	PK-PF633P-2-01	PromoFluor-633P, amino-modified	1 mg
	PK-PF633P-3-01	PromoFluor-633P, maleimide	1 mg
	PK-PF633P-3-05	PromoFluor-633P, maleimide	5 mg
Structure & Spectra	<b>Chemical Structure</b>		<b>Absorption and Emission Spectra</b>
			
Absorption / Emission max.	635 nm / 658 nm (in ethanol); 622 nm / 647 nm (in PBS)		
Molar Absorbance	200.000 M <sup>-1</sup> cm <sup>-1</sup>		
Molecular Data	<b>Modification</b>	<b>Molecular Weight</b>	<b>Molecular Formula</b>
	Carboxylic acid	969.03 g·mol <sup>-1</sup>	C <sub>38</sub> H <sub>47</sub> N <sub>2</sub> O <sub>15</sub> S <sub>4</sub> * 3 Na
	NHS-ester	1066.1 g·mol <sup>-1</sup>	C <sub>42</sub> H <sub>50</sub> N <sub>3</sub> O <sub>17</sub> S <sub>4</sub> * 3 Na
	Maleimide	1091.16 g·mol <sup>-1</sup>	C <sub>44</sub> H <sub>53</sub> N <sub>4</sub> O <sub>16</sub> S <sub>4</sub> * 3 Na
	Amino-derivative	989.12 g·mol <sup>-1</sup>	C <sub>40</sub> H <sub>54</sub> N <sub>4</sub> O <sub>14</sub> S <sub>4</sub> * 2 Na
Formulation	Solid		
Solubility	Soluble in water, 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 <a href="http://www.promkine.info">www.promkine.info</a> .		

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