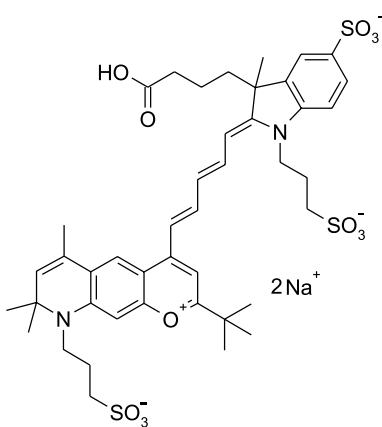
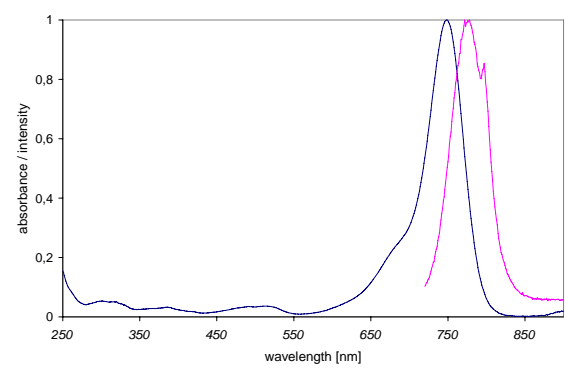


## Instruction Manual

Catalog Number	PK-PF750-0-01	PromoFluor-750, carboxylic acid	1 mg
	PK-PF750-0-05	PromoFluor-750, carboxylic acid	5 mg
	PK-PF750-1-01	PromoFluor-750, NHS ester	1 mg
	PK-PF750-1-01A	PromoFluor-750, NHS ester	5 x 0.2 mg
	PK-PF750-1-05	PromoFluor-750, NHS ester	5 mg
	PK-PF750-2-01	PromoFluor-750, amino-modified	1 mg
	PK-PF750-3-01	PromoFluor-750, maleimide	1 mg
	PK-PF750-3-05	PromoFluor-750, maleimide	5 mg

Structure & Spectra	<b>Chemical Structure</b>	<b>Absorption and Emission Spectra</b>
		

Absorption / Emission max.	748 nm / 772 nm (in ethanol)
----------------------------	------------------------------

Molar Absorbance	270.000 M <sup>-1</sup> cm <sup>-1</sup>
------------------	--

Molecular Data	Modification	Molecular Weight	Molecular Formula
	Carboxylic acid	931.07 g·mol <sup>-1</sup>	C <sub>43</sub> H <sub>52</sub> N <sub>2</sub> O <sub>12</sub> S <sub>3</sub> Na <sub>2</sub>
	NHS-ester	1028.15 g·mol <sup>-1</sup>	C <sub>47</sub> H <sub>55</sub> N <sub>3</sub> O <sub>14</sub> S <sub>3</sub> Na <sub>2</sub>
	Maleimide	1053.20 g·mol <sup>-1</sup>	C <sub>49</sub> H <sub>58</sub> N <sub>4</sub> O <sub>13</sub> S <sub>3</sub> Na <sub>2</sub>
	Amino-derivative	951.17 g·mol <sup>-1</sup>	C <sub>45</sub> H <sub>59</sub> N <sub>4</sub> O <sub>11</sub> S <sub>3</sub> Na

Formulation	Solid
-------------	-------

Solubility	Soluble in water and 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.promkine.info">www.promkine.info</a> .
--------------	--

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