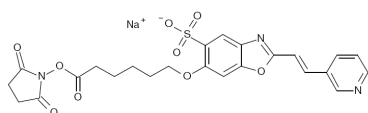


## Instruction Manual

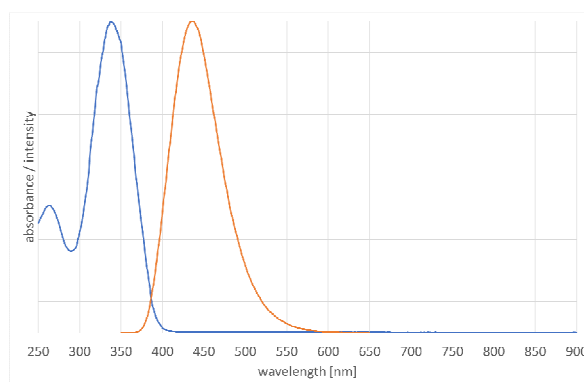
Catalog Number	PK-PF350P-0-01	PromoFluor-350P, carboxylic acid	1 mg
	PK-PF350P-0-05	PromoFluor-350P, carboxylic acid	5 mg
	PK-PF350P-1-01	PromoFluor-350P, NHS-ester	1 mg
	PK-PF350P-1-01A	PromoFluor-350P, NHS-ester	5 x 0.2 mg
	PK-PF350P-1-05	PromoFluor-350P, NHS-ester	5 mg
	PK-PF350P-2-01	PromoFluor-350P, amino-modified	1 mg
	PK-PF350P-3-01	PromoFluor-350P, maleimide	1 mg
	PK-PF350P-3-05	PromoFluor-350P, maleimide	5 mg

## Structure & Spectra

### Chemical Structure



### Absorption and Emission Spectra



**Absorption / Emission max.** 338 nm / 436 nm (in PBS)

**Molar Absorbance** 33.000 M<sup>-1</sup>cm<sup>-1</sup>

### Molecular Data

Modification	Molecular Weight	Molecular Formula
Carboxylic acid	454.43 g·mol <sup>-1</sup>	C <sub>20</sub> H <sub>19</sub> N <sub>2</sub> O <sub>7</sub> S * Na
NHS-ester	551.50 g·mol <sup>-1</sup>	C <sub>24</sub> H <sub>22</sub> N <sub>3</sub> O <sub>9</sub> S * Na
Maleimide	576.55 g·mol <sup>-1</sup>	C <sub>26</sub> H <sub>25</sub> N <sub>4</sub> O <sub>8</sub> S * Na
Amino-derivative	474.53 g·mol <sup>-1</sup>	C <sub>22</sub> H <sub>26</sub> N <sub>4</sub> O <sub>6</sub> S

**Formulation** Solid

**Solubility** Soluble in water or aqueous buffers such as PBS.

**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 [www.promokine.info](http://www.promokine.info).

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