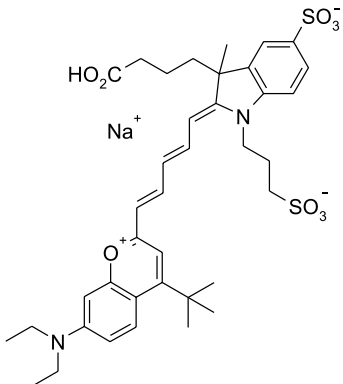
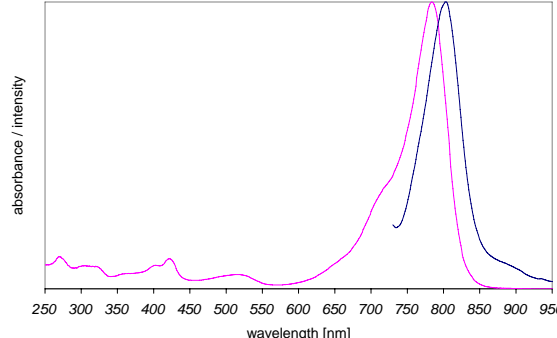


Instruction Manual

| | | | |
|----------------|----------------|---------------------------------|------------|
| Catalog Number | PK-PF780-0-01 | PromoFluor-780, carboxylic acid | 1 mg |
| | PK-PF780-0-05 | PromoFluor-780, carboxylic acid | 5 mg |
| | PK-PF780-1-01 | PromoFluor-780, NHS ester | 1 mg |
| | PK-PF780-1-01A | PromoFluor-780, NHS ester | 5 x 0.2 mg |
| | PK-PF780-1-05 | PromoFluor-780, NHS ester | 5 mg |
| | PK-PF780-2-01 | PromoFluor-780, amino-modified | 1 mg |
| | PK-PF780-3-01 | PromoFluor-780, maleimide | 1 mg |
| | PK-PF780-3-05 | PromoFluor-780, maleimide | 5 mg |

| | | |
|---------------------|--|---|
| Structure & Spectra | Chemical Structure | Absorption and Emission Spectra |
| |  |  |

| | |
|----------------------------|------------------------------|
| Absorption / Emission max. | 783 nm / 800 nm (in ethanol) |
|----------------------------|------------------------------|

| | |
|------------------|--|
| Molar Absorbance | 170.000 M ⁻¹ cm ⁻¹ |
|------------------|--|

| Molecular Data | Modification | Molecular Weight | Molecular Formula |
|----------------|------------------|----------------------------|--|
| | Carboxylic acid | 762.92 g·mol ⁻¹ | C ₃₈ H ₄₇ N ₂ O ₉ S ₂ Na |
| | NHS-ester | 860.00 g·mol ⁻¹ | C ₄₂ H ₅₀ N ₃ O ₁₁ S ₂ Na |
| | Maleimide | 885.05 g·mol ⁻¹ | C ₄₄ H ₅₃ N ₄ O ₁₀ S ₂ Na |
| | Amino-derivative | 783.03 g·mol ⁻¹ | C ₄₀ H ₅₄ N ₄ O ₈ S ₂ |

| | |
|-------------|-------|
| Formulation | Solid |
|-------------|-------|

| | |
|------------|--|
| Solubility | Soluble in water, methanol, DMF and 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 www.promokine.info . |
|--------------|--|

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