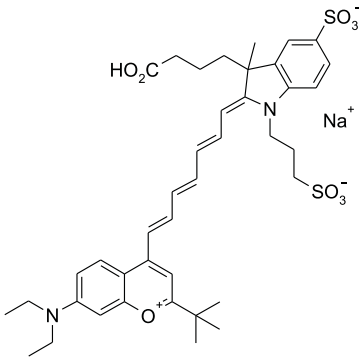


Instruction Manual

| | | | |
|---|--|---|--|
| Catalog Number | PK-PF840-0-01 | PromoFluor-840, carboxylic acid | 1 mg |
| | PK-PF840-0-05 | PromoFluor-840, carboxylic acid | 5 mg |
| | PK-PF840-1-01 | PromoFluor-840, NHS ester | 1 mg |
| | PK-PF840-1-01A | PromoFluor-840, NHS ester | 5 x 0.2 mg |
| | PK-PF840-1-05 | PromoFluor-840, NHS ester | 5 mg |
| | PK-PF840-3-01 | PromoFluor-840, maleimide | 1 mg |
| | PK-PF840-3-05 | PromoFluor-840, maleimide | 5 mg |
| | Structure & Spectra | Chemical Structure  | |
| Absorption / Emission max. 844 nm / 884 nm (in ethanol) | | | |
| Molar Absorbance | 220.000 M ⁻¹ cm ⁻¹ | | |
| Molecular Data | Modification | Molecular Weight | Molecular Formula |
| | Carboxylic acid | 788.96 g·mol ⁻¹ | C ₄₀ H ₄₉ N ₂ O ₉ S ₂ Na |
| | NHS-ester | 886.04 g·mol ⁻¹ | C ₄₄ H ₅₂ N ₃ O ₁₁ S ₂ Na |
| | Maleimide | 911.09 g·mol ⁻¹ | C ₄₆ H ₅₅ N ₄ O ₁₀ S ₂ Na |
| | Amino-derivative | 809.06 g·mol ⁻¹ | C ₄₂ H ₅₆ N ₄ O ₈ S ₂ |
| 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 www.promokine.info . | | |

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