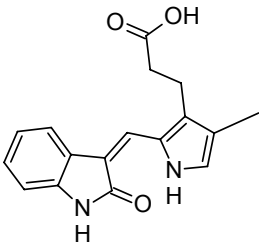


3-[3-(2-Carboxyethyl)-4-methylpyrrol-2-methylidanyl]-2-indolinone

Instruction Manual

Catalog Number	PK-CA577-1645-05
Description	A potent and selective inhibitor of fibroblast growth factor receptor (FGFR) (IC ₅₀ = 0.03 μM for FGFR1) and vascular endothelial growth factor receptors (VEGFR) (IC ₅₀ = 0.02 μM for VEGFR2). In addition, integrin β 4-induced differentiation of neural stem cells was attenuated by SU5402.
Quantity	500 μg
Sequence / Molecular Weight / Molecular Formula	296.32 Da; C ₁₇ H ₁₆ N ₂ O ₃
Chemical Structure	
Purity	≥ 97% as determined by HPLC
Appearance / Formulation / Solubility	Yellow Solid; Soluble in DMSO (100 mM)
Storage & Stability	Store at -20°C.
Applications	see Description
References	NA
Caution	Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

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