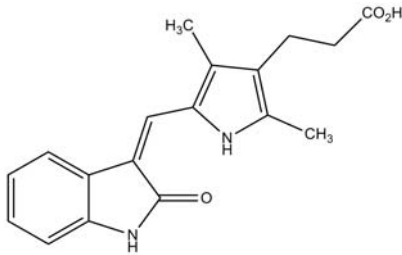


## 5-[1,2-Dihydro-2-oxo-3H-indol-3-ylidene)methyl]-2,4-dimethyl-1H-pyrrole-3-propanoic acid

## Instruction Manual

Catalog Number	PK-CA577-1931-25
Description	A potent, cell-permeable and ATP-competitive inhibitor of PDGF, VEGF and FGF receptor tyrosine kinases (RTKs) with IC50 values of 0.06, 2.43, 3.04 and > 100 $\mu$ M at PDGFR $\beta$ , VEGFR2, FGFR1 and EGFR respectively. Acts as a potent antiangiogenic and antitumor agent.
Quantity	25 mg
Sequence / Molecular Weight / Molecular Formula	310.35 Da; C <sub>18</sub> H <sub>18</sub> N <sub>2</sub> O <sub>3</sub>
Chemical Structure	
Purity	$\geq$ 98%
Appearance / Formulation / Solubility	Orange-red Solid; Soluble in DMSO (100 mM)
Storage & Stability	Store at -20°C. Protect from moisture.
Applications	see Description
References	Laird, D.A., et al. (2000). Cancer Res. 60, 4152-4260. Sun, L., et al. (1999). J. Med. Chem. 42, 5120-5130.
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.