

## Instruction Manual

Catalog Number	PK-CA577-1571-5
Description	Selectively inhibits PDGF receptor autophosphorylation (IC <sub>50</sub> = 0.3-0.5 μM) and selectively inhibits PDGF dependent DNA synthesis. Potent inhibitor of human and porcine smooth muscle cell (SMC) proliferation in vitro and of rat SMC proliferation in vivo.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	234,3 kDa; C <sub>16</sub> H <sub>14</sub> N <sub>2</sub>
Purity	>99%
Appearance / Formulation / Solubility	Off-white solid. Soluble in DMSO.
Storage & Stability	Store at -20°C. Do not expose to sunlight or fluorescent lights for extended periods.
Applications	NA
References	NA
Caution	Do not ingest, swallow, or inhale. Do not get in eyes or on skin. Care should be taken to prevent contact through all routes of exposure.

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