

## Instruction Manual

Catalog Number	PK-CA577-1533-100
Description	A reported tyrosine kinase inhibitor.
Quantity	100 mg
Sequence / Molecular Weight / Molecular Formula	270.2 Da; C <sub>15</sub> H <sub>10</sub> O <sub>5</sub>
Purity	>99%
Appearance / Formulation / Solubility	Off-white powder. Soluble in DMSO.
Storage & Stability	Store desiccated at -20°C. Stable for at least 1 year under proper storage conditions.
Applications	NA
References	1. Shertzner, H.O., et al. (1999) Chem. Biol. Interact. 123:31 2. Piskula, M.K., et al. (1999) FEBS Lett. 447:287.
Caution	NA

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