

## Instruction Manual

|   |  |
|---|--|
| Catalog Number                                  | PK-CA577-1533-10   |
| Description                                     | A reported tyrosine kinase inhibitor.  |
| Quantity  | 10 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.<br>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.