

Instruction Manual

| | |
|---|---|
| Catalog Number | PK-CA577-1048-010 |
| Description | Potent inhibitor of protein kinase C and most other kinases, including tyrosine kinases. Inhibits topoisomerase II directly by blocking transfer of phosphodiester bonds from DNA to active site tyrosine. Induces apoptosis in CHO cells at 1 μ M. |
| Quantity | 100 μ g |
| Sequence / Molecular Weight / Molecular Formula | 466.5 Da; C ₂₅ H ₂₆ N ₄ O ₃ |
| Purity | >99% by TLC |
| Appearance / Formulation / Solubility | Soluble in DMSO (25 mg/ml). |
| Storage & Stability | Store powder desiccated at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months. |
| Applications | see Description |
| References | <ol style="list-style-type: none"> Seynaeve, C.M., et al. (1994) Mol. Pharmacology 45:1207. Wang, X., et al. (1996) EMBO J. 15:1012. Lassota, P., et al. (1996) J. Biol. Chem. 271:26418. |
| Caution | NA |

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