

Instruction Manual

Catalog Number	PK-CA577-1048-1
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	1 mg
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"> 1. Seynaeve, C.M., et al. (1994) Mol. Pharmacology 45:1207. 2. Wang, X., et al. (1996) EMBO J. 15:1012. 3. Lassota, P., et al. (1996) J. Biol. Chem. 271:26418.
Caution	NA

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