

Instruction Manual

Catalog Number	PK-CA577-1551
Description	A potent synthetic anti-estrogenic agent. A reversible inhibitor of protein kinase C (IC ₅₀ = 10 mM). Induces apoptosis in human malignant glioma cell lines and inhibits prostate cancer cell growth by induction of p21 protein.
Quantity	1 g
Sequence / Molecular Weight / Molecular Formula	371.5 x 192.1 Da; C ₂₅ H ₂₉ NO x C ₆ H ₈ O ₇
Purity	>99%
Appearance / Formulation / Solubility	White, crystalline solid. Soluble in methanol, slightly soluble in water.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	see Description
References	O'Brian, C.A., et al. (1985) Cancer Res. 45, 2462. Couldwell, W.T. et al. (1994) FEBS Lett. 345, 43.
Caution	Toxic/Irritant.

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