

Instruction Manual

Catalog Number	PK-CA577-1543-25
Description	Potent inhibitor of protein phosphatases 1 and 2A in numerous cell types. Does not affect activity of acid phosphatase, alkaline phosphatase and tyrosine phosphatase. Non-phorbol type tumor promoter. Induces apoptosis in human breast carcinoma cells (MB-231 and MCF-7) and in myeloid cells, but inhibits glucocorticoid-induced apoptosis in T-cell hybridomas. Has shown contractile effect on smooth muscle and heart muscle.
Quantity	25 µg
Sequence / Molecular Weight / Molecular Formula	805 Da; C ₄₄ H ₅₅ O ₁₃
Purity	>98%
Appearance / Formulation / Solubility	White crystalline solid. Soluble in DMSO or ethanol. Use freshly.
Storage & Stability	Store at -20°C in a dry environment. Protect from light.
Applications	Can be used for cell viability and cytotoxicity assays as well as for drug screening, NADH/NADPH and SOD detection etc.
References	NA
Caution	Highly irritant tumor promoter.

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