

Instruction Manual

Catalog Number	PK-CA577-1549-10
Description	Activator of p54 and MAP kinases. Synergistic with growth factors and phorbol esters to superinduce cFas and c-Jun, by acting as a potent signaling agonist. Induces apoptosis in the human monoblastoid cell line. Isolated from <i>Streptomyces griseolus</i> .
Quantity	10 mg
Sequence / Molecular Weight / Molecular Formula	[2-(p-Methoxybenzyl)-3,4-pyrrolidinediol-3-acetate]; 265,3 Da; C ₁₄ H ₁₉ NO ₄
Purity	>98%
Appearance / Formulation / Solubility	White crystalline solid. Soluble in DMSO
Storage & Stability	Store at -20°C. Store solution at -20°C for up to 3 months.
Applications	see Description
References	1. Kardalidou, E., et al. (1994) Mol. Cell. Biol. 14:1066. 2. Kochi, S.K. and Collier, R.J. (1993) Exp. Cell. Res. 208:296.
Caution	NA

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