

Instruction Manual

Catalog Number	PK-CA577-1036-5MG
Description	An antineoplastic antibiotic. Inhibits DNA-primed RNA polymerase by complexing with DNA via deoxyguanosine residues. A DNA-intercalator with growth inhibitory activity against certain leukemias and sarcomas. At higher concentrations, DNA polymerase is inhibited. Inhibits cell growth and colony formation in synchronized HeLa cells. A potent inducer of apoptosis in many cell lines (1).
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	1255,5 Da; C ₆₂ H ₈₆ N ₁₂ O ₁₆
Purity	>98% by TLC
Appearance / Formulation / Solubility	Red-orange solid. Soluble in DMSO (20 mg/ml).
Storage & Stability	Store at -20°C. Stable for up to 1 year as solid and up to 3 months in solution at -20°C.
Applications	see Description
References	1. Verhaegen, S., et al. (1995) Biochem. Pharmacol. 50:1021-1029.
Caution	NA

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