

## Instruction Manual

Catalog Number	PK-CA577-1036-50
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	50 µl
Sequence / Molecular Weight / Molecular Formula	1255,5 Da; C <sub>62</sub> H <sub>86</sub> N <sub>12</sub> O <sub>16</sub>
Purity	>98% by HPLC
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Stable for up to 3 months under proper storage conditions.
Applications	We recommend using 1000X dilutions for inducing apoptosis in Jurkat cells (e.g., Add 1 µl to 1 ml of culture medium). However, the optimal doses may vary for different cells and culture conditions.
References	1. Martin, S.J., et al. (1995) J. Exp. Med. 182:1545-1556.
Caution	NA

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