

Instruction Manual

Catalog Number	PK-CA577-1527-5
Description	Anti-tumor antibiotic. Inhibitor of Reverse Transcriptase and RNA Polymerase. Immunosuppressive agent. Highly effective myotoxin that inhibits topoisomerase II. Binds to nucleic acids, presumably by specific intercalation into the DNA double helix, thereby inhibiting nucleic acid synthesis. Induces apoptosis.
Quantity	5 mg
Sequence / Molecular Weight / Molecular Formula	543.5 Da; C ₂₅ H ₂₉ NO ₁₁ .HCl
Purity	>98% by TLC
Appearance / Formulation / Solubility	Orange-Red Power (m.p. 216°C). Soluble in water.
Storage & Stability	Store at -20°C in a dry environment. Protect it from light. Stable for 1 year under proper storage conditions.
Applications	We recommend using 1000X dilutions for inducing apoptosis in Jurkat cells. However, the optimal doses may vary for different cells and culture conditions.
References	1. Skladanowski A. & Konopa J. (1993) Biochem. Pharmacol. 46: 375-383.
Caution	Very toxic, may cause cancer!

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