

Instruction Manual

Catalog Number	PK-CA577-1043-250MG
Description	A topoisomerase II inhibitor. A derivative of podophyllotoxin that has major activity against a number of tumors, including germ cell neoplasms, small cell lung cancer, and malignant lymphoma. Induces apoptosis in human T cells, mouse thymocytes, and HL-60 human leukemia cells (1).
Quantity	250 mg
Sequence / Molecular Weight / Molecular Formula	588.6 Da; C ₂₉ H ₃₂ O ₁₃
Purity	>98% by HPLC
Appearance / Formulation / Solubility	White crystalline powder. Soluble in DMSO.
Storage & Stability	Store powder desiccated at -20°C for up to one year. Store solutions at -20°C for up to 3 months.
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. Onishi, Y., et al. (1993) Biochem. Biophys. Acta 1175:147-154.
Caution	NA

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