

## Instruction Manual

Catalog Number	PK-CA577-1043-100
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	100 µl
Sequence / Molecular Weight / Molecular Formula	588,6 Da; C <sub>29</sub> H <sub>32</sub> O <sub>13</sub>
Purity	>97% by HPLC
Appearance / Formulation / Solubility	100 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 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. 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.