

Camptothecin (2 mM)

Ready-to-use

Instruction Manual

Catalog Number	PK-CA577-1039-1
Description	Reversibly inhibits nuclear topoisomerase I by binding to and stabilizing the topoisomerase-DNA covalent complex. Inhibits Tat-mediated transcription of HIV-1. Induces apoptosis in Jurkat, osteosarcoma and hepatoma cells.
Quantity	1 ml (2 mM)
Sequence / Molecular Weight / Molecular Formula	348,4 Da; C ₂₀ H ₁₆ N ₂ O ₄
Purity	>98% by TLC
Appearance / Formulation / Solubility	2 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 6 months at -20°C 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	NA
Caution	Potential mutagen/reproductive hazard!

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