

Instruction Manual

Catalog Number	PK-CA577-1039-50MG
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	50 mg
Sequence / Molecular Weight / Molecular Formula	348.4 Da; C ₂₀ H ₁₆ N ₂ O ₄
Purity	>98% by TLC
Appearance / Formulation / Solubility	Pale-yellow solid. Soluble in DMSO (10 mg/ml).
Storage & Stability	Store powder desiccated at -20°C for up to one year. Store solutions at -20°C for up to 3 months.
Applications	see Description
References	1. Nicholson, D.W., et al. (1995) Nature 376:37-40. 2. Suzuki, A. and M. Kato (1996) Exp. Cell. Res. 227:154-163.
Caution	Potential mutagen/reproductive hazard!

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