

Instruction Manual

Catalog Number	PK-CA577-K121-5			
Description	The set contains 5 ready-to-use chemical reagents that induce apoptosis through different mechanisms. Actinomycin D, an antineoplastic antibiotic, inhibits RNA synthesis. Camptothecin, an inhibitor of nuclear topoisomerase, induces apoptosis in many types of cells. Cycloheximide, an active antibiotic against many yeast and fungi, inhibits protein synthesis. Dexamethasone, an active and highly stable glucocorticoid, probably induces apoptosis by binding and activating the intracellular glucocorticoid receptor. Etoposide, a derivative of podophyllotoxin, inhibits topoisomerase activity.			
Quantity	1 kit			
Kit Components	Component	Concentration	Volume	Cap color
	Actinomycin D	10 mM	50 µl	Violet
	Camptothecin	5.7 mM	1 ml	Red
	Cycloheximide	100 mM	1 ml	Green
	Dexamethasone	10 mM	1 ml	Blue
	Etoposide	100 mM	100 µl	Yellow
Applications / Assay Protocol	We recommend using 1000X dilutions for inducing apoptosis in cell cultures. However, the optimal doses may vary for different cells and culture conditions. See references: 1. Onishi, Y., et al. (1993) Biochem. Biophys. Acta 1175:147-154. 2. Fearnhead, H.O., et al. (1994) Biochem. Pharmacol. 48:1073-1079.			
Storage & Stability	Store at –20°C.			
Related Products	Please visit our website www.promocell.com to find a variety of products for apoptosis research (e.g. assays, inducers & inhibitors, antibodies, ELISAs as well as cytokines & growth factors).			

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