



## **PRODUCT INFORMATION & MANUAL**

# **Apoptosis Inducer Compounds Kit**

***NBP2-54825***

For research use only.

Not for diagnostic or therapeutic procedures.

[www.novusbio.com](http://www.novusbio.com) - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - [technical@novusbio.com](mailto:technical@novusbio.com)

Novus kits are guaranteed for 6 months from date of receipt

# Ready-to-use Apoptosis Inducer Set

**CATALOG #:** NBP2-54825

**LOT #:**

**STORAGE CONDITIONS:** Store at –20°C.

**SHELF LIFE:** 6 months under proper storage conditions

## 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.

## PACKAGE CONTENTS:

Concentration	Description	Volume	Cap Color	Part Number
10 mM	Actinomycin D	50 µl	Violet	1036-50
2 mM	Camptothecin	1 ml	Red	1039-1
100 mM	Cycloheximide	1 ml	Green	1041-1
10 mM	Dexamethasone	1 ml	Blue	1042-1
100 mM	Etoposide	100 µl	Yellow	1043-100

**Note:** All components were formulated in pure DMSO.

## RECOMMENDED USAGE:

We recommend using 1000X dilutions for inducing apoptosis in cell cultures. However, the optimal doses may vary for different cells and culture conditions.

**FOR RESEARCH USE ONLY! Not to be used in humans.**