

Ready-to-use Apoptosis Inducer Set

CATALOG #: K121-5

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
10 mM	Actinomycin D	50 µl	Violet
2 mM	Camptothecin	1 ml	Red
100 mM	Cycloheximide	1 ml	Green
10 mM	Dexamethasone	1 ml	Blue
100 mM	Etoposide	100 µl	Yellow

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.

REFERENCES:

- Onishi, Y., *et al.* (1993) *Biochem. Biophys. Acta* **1175**:147-154.
- Fearnhead, H.O., *et al.* (1994) *Biochem. Pharmacol.* **48**:1073-1079.

FOR RESEARCH USE ONLY! Not to be used in human.

RELATED PRODUCTS:

Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents
- Apoptosis Inducers and Set
- Apoptosis siRNA Vectors

Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Fractionation kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- High Throughput Apoptosis/Cell Viability Assay Kits
- LDH-Cytotoxicity Assay Kit
- Bioluminescence Cytotoxicity Assay Kit
- Live/Dead Cell Staining Kit

Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Fluorometric & Colorimetric Assays & Drug Discovery Kits
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits

Adipocyte & Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP & PLTP Activity Assay & Drug Discovery Kits
- Total Cholesterol Quantification Kit

Molecular Biology & Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Minutes DNA Ligation Kit
- 20 Minutes Gel Staining/Destaining Kit
- β-Galactosidase Staining Kit & Luciferase Reporter Assay Kit

Growth Factors and Cytokines

- Adiponectin/Resistin/Leptin and their Antibodies
- Recombinant Protein A and Protein G
- Recombinant Complement C5a
- Recombinant Cytokines and Growth Factors

Monoclonal and Polyclonal Antibodies