



## Apoptosis Inducers & Set

### INTRODUCTION

Apoptosis is commonly defined as programmed mechanism of cell death and is a highly regulated process that can be initiated by a variety of internal and external stimuli (inducers). G-Biosciences's Apoptosis Inducers are ready-to-use solutions for inducing apoptosis in various cell types. They are supplied individually as well as in a set of 6 inducers. These inducers include: Actinomycin D, Camptothecin, Cycloheximide, Dexamethasone, Doxorubicin and Etoposide and can be used in 1000X dilution. These agents induce apoptosis by different mechanisms. Depending on the need, individual as well as set of inducers may be ordered.

### KIT COMPONENTS:

Catalog #	Name of Inducer	Size
API-01	Actinomycin D (10mM)	1 x 50µl
API-02	Camptothecin (2mM)	1 x 1ml
API-03	Cycloheximide (100mM)	1 x 1ml
API-04	Dexamethasone (10mM)	1 x 1ml
API-05	Doxorubicin (2mM)	1 x 50µl
API-06	Etoposide (100mM)	1 x 50µl
API-16S	Apoptosis Inducer Set	Set of all 6 inducers

### STORAGE CONDITION

It is shipped in Blue Ice. Upon arrival, store at -20°C upon arrival and is stable for 6-8 months if stored properly.

### PROTOCOL

#### Preparation before Use:

Take out the vial(s) from refrigerator and thaw to RT. Vortex to mix the contents.

These apoptotic agents induce apoptosis by different mechanisms, e.g. ActinomycinD inhibits RNA synthesis; Camptothecin inhibits enzyme nuclear topoisomerase; Cycloheximide inhibits protein synthesis; Dexamethasone activates intracellular glucocorticoid receptor; Doxorubicin inhibits nucleic acid synthesis presumably by specific intercalation into the DNA double helix; and Etoposide inhibits topoisomerase activity. Depending on the condition and cell type, these inducers can be used in 1000X dilution, i.e. 1µl to 1ml cell culture (researchers may also decide their own concentrations).

### RELATED PRODUCTS:

#### 1. CasPASE™ -Apoptosis Assay Kit (Cat # 786-200 to 786-205)

The assay kit provides a simple method for assaying various caspase activity. For details please visit our web site at [www.GBiosciences.com](http://www.GBiosciences.com).

#### 2. CasPASE Substrates and Inhibitors

Ready to use caspase substrates and inhibitors are provided individually as well as in set. For details please visit our web site at [www.GBiosciences.com](http://www.GBiosciences.com)

For other related products, visit our web site [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Rev 11.03.08-SA/MM/IA

