

# Apoptosis inducer set

The set contains 5 *ready-to-use* 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.

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-850-2350	1Set
--------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Contents	Contains 5 <i>ready-to-use</i> reagents: <ul style="list-style-type: none"><li>– Actinomycin D, 50µl (10mM), violet cap</li><li>– Camptothecin, 1ml (2mM), red cap</li><li>– Cycloheximide, 1ml (100mM), green cap</li><li>– Dexamethasone, 1ml (10mM), blue cap</li><li>– Etoposide, 100µl (100mM), yellow cap</li></ul>
----------	---

Formulation	Liquid. In DMSO
-------------	-----------------

Recommendation Dilutions/Conditions	We recommend using 1,000x dilutions for inducing apoptosis in cell cultures. Optimal doses may vary for different cells and culture conditions and must be determined individually.
-------------------------------------	---

Last modified: June 13, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)