

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-850-235-KI01	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 *ready-to-use* reagents:

- Actinomycin D, 50µl (10mM), violet cap
- Camptothecin, 1ml (2mM), red cap
- Cycloheximide, 1ml (100mM), green cap
- Dexamethasone, 1ml (10mM), blue cap
- Etoposide, 100µl (100mM), yellow cap

Formulation Liquid. In DMSO

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



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)