

# Apoptosis Activator 2

## Apoptosis inducer

Induces apoptosis by promoting the oligomerization of Apaf-1 into the mature, functional apoptosome, triggering caspase-9 activation resulting in activated caspase-3. Shows strong selectivity for cancer cells against normal cell lines. Acts in a similar way like apoptosis inducer PETCM (Prod. No. ALX-420-031).

Citations: 2

[View Online »](#)

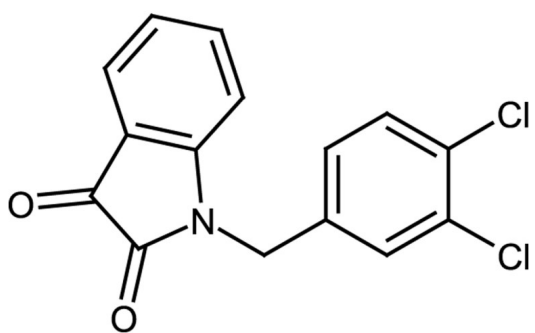
## Ordering Information

[Order Online »](#)

ALX-420-035-M005	5mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 1-[(3,4-Dichlorophenyl)methyl]-1H-indole-2,3-dione

**Appearance** Orange to red solid.

**CAS** 79183-19-0

**Formula**  $C_{15}H_9NO_2Cl_2$

**MW** 306.1

**MeltingPoint** 174-175°C

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (5mg/ml), methanol (5mg/ml) or 100% ethanol.

**Last modified: May 29, 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)