

# Necrox-5™ methanesulfonate

## Necrosis inhibitor

Cell permeable necrosis inhibitor with antioxidant activity. It localizes mostly in the mitochondria. Selectively blocks oxidative stress-induced necrotic cell death (0.1µM NecroX™-5 prevented ~50% cell death in H9C2 cells exposed to 400µM t-BuOOH for 2 hours LDH assay used). Does not protect against staurosporine or etoposide-induced apoptosis. Protects cells against cold shock, hypoxia and oxidative stress *in vitro*, as well as CCl<sub>4</sub>-induced acute liver injury and chronic liver fibrosis in rodent models.

Citations: 5

[View Online »](#)

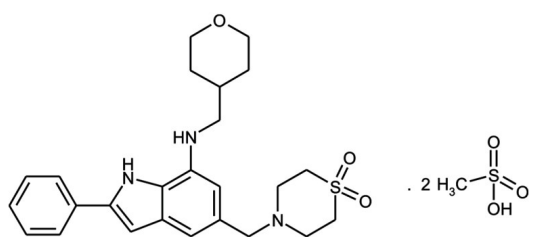
## Ordering Information

[Order Online »](#)

ALX-430-167-M001	1mg
ALX-430-167-M005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and keep under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	[5-(1,1-Dioxo-thiomorpholin-4-ylmethyl)-2-phenyl-1H-indol-7-yl]-(tetrahydro-pyran-4-ylmethyl)-amine
Appearance	Off-white to light pink solid.
Formula	$C_{25}H_{31}N_3O_3S \cdot 2 CH_4O_3S$
MW	645.8
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.



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-  
fr@enzolifesciences.com](mailto:info-fr@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)