

# Necrox-5™ hydrochloride

## 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: 7

[View Online »](#)

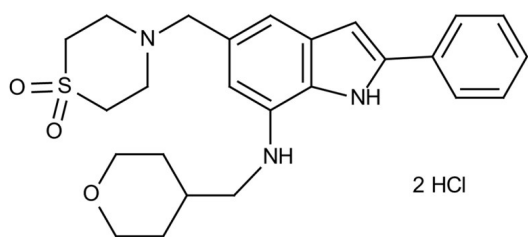
## Ordering Information

[Order Online »](#)

ALX-430-247-M001	1mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. 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	Yellow solid.
Application Notes	<p>Can be used for <i>in vitro</i> and <i>in vivo</i> applications.</p> <p>For <i>in vitro</i> experiments, dissolve Necrox-5™ in DMSO to 20-50mM stock solution and dilute further in cell culture media. Treat cells with 30µM to 0.1µM final concentration.</p> <p>For <i>in vivo</i> experiments, dissolve the product in DMSO at a concentration of 10mg/ml stock. Do not use PBS and Water!</p> <p>Optimal conditions must be determined individually for each application.</p>
Formula	$C_{25}H_{31}N_3O_3S \cdot 2 HCl$
Identity	Determined by LC-MS.
MW	453.6 . 72.9
Purity	≥96% (HPLC, TLC)
Solubility	Soluble in DMSO (>10mg/ml).



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)