

# Necrostatin-7

## Necroptosis inhibitor

Necroptosis inhibitor. The biological activity of necrostatin-7 (Nec-7) is different from that of necrostatin-1 to -5 (Nec-1 to -5), as Nec-7 does not inhibit RIP1 kinase. Nec-7 may target an additional regulatory molecule in the pathway.

Citations: 2

[View Online »](#)

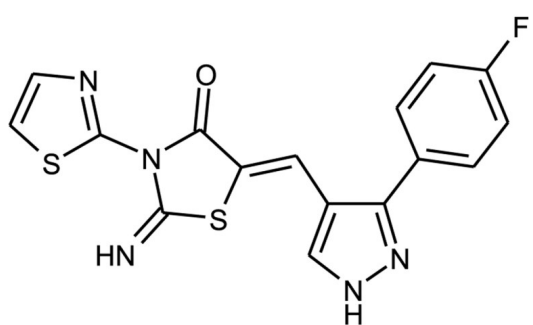
## Ordering Information

[Order Online »](#)

ALX-430-170-M005	5mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep cool and dry. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	Yellow solid.
CAS	351062-08-3
Formula	$C_{16}H_{10}N_5OS_2F$
MW	371.4
Purity	≥98% (HPLC, TLC)
Solubility	Soluble in DMSO or 100% ethanol.



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)