

# N-(1-Naphthalenylsulfonyl)-Ile-Trp-aldehyde

## Cathepsin-L inhibitor

Potent, selective and reversible inhibitor of human cathepsin L ( $IC_{50}$  =1.9nM). Inhibits release of  $Ca^{2+}$  and hydroxyproline from bone *in vitro*.

Citations: 7

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-260-133-M001	1mg
------------------	-----

## 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** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N-(1Naphthalenylsulfonyl)- IW-CHO

**Appearance** White to off-white powder.

**CAS** 161709-56-4

**Formula**  $C_{27}H_{29}N_3O_4S$

**MW** 491.6

**Purity** ≥97% (HPLC)

**Solubility** Soluble in DMSO or ethyl acetate (50mg/ml).

**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)