

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
ALX-260-133-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 -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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com