

E-64

Protease inhibitor

Irreversible inhibitor of cysteine proteinases like papain, calpain, cathepsin B, L and H. Acts by forming a thioether bond with the thiol of the active cysteine. Does not inhibit serine proteinases. Inhibits degradation of autophagic cargo in autophagolysosomes.

Citations: 7

[View Online »](#)

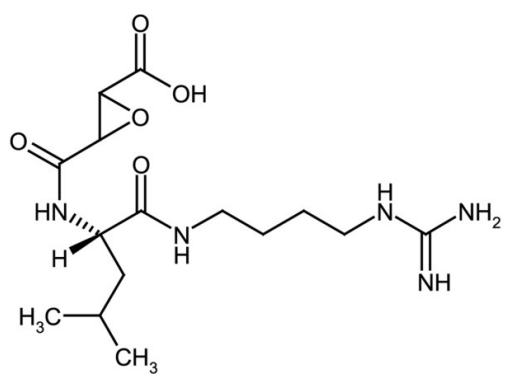
Ordering Information

[Order Online »](#)

ALX-260-007-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



260-007
E-64

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-[N-(L-3-trans-Carboxyoxirane-2-carbonyl)-L-leucyl]- agmatine, L-trans-Epoxy succinyl-leucylamide-(4-guanido)- butane
Appearance	White to off-white solid.
CAS	66701-25-5
Couple Target	Calpain, Cathepsin, Papain
Couple Type	Inhibitor
Formula	$C_{15}H_{27}N_5O_5 \cdot 0.5 H_2O$
MW	357.4
Purity	≥98%
Solubility	Soluble in water/ethanol 1:1 mixture (15 mg/ml).
Source	Synthetic.

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