

Leupeptin (synthetic)

Cathepsin B inhibitor

Leupeptin is a competitive and reversible inhibitor of serine and cysteine proteinases like calpain, porcine kallikrein, trypsin, plasmin, papain and cathepsin B. Does not inhibit chymotrypsin and thrombin. Also known to inhibit activation-induced programmed cell death and restoration of defective immune responses of HIV⁺ donors.

Citations: 19

[View Online »](#)

Ordering Information

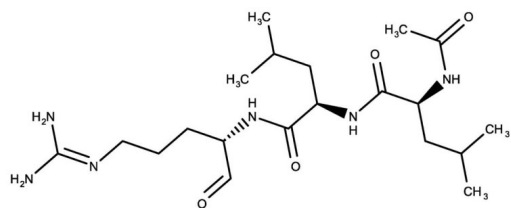
[Order Online »](#)

ALX-260-009-M025	25mg
ALX-260-009-M100	100mg

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Acetyl-L-leucyl-L-leucyl-L-argininal . hemisulfate
Appearance	White to off-white powder.
CAS	103476-89-7
Formula	$C_{20}H_{38}N_6O_4 \cdot 0.5 H_2SO_4$
MI	14: 5456
MW	426.6 . 49.0
Purity	≥97% (HPLC)
Solubility	Soluble in water, methanol or 100% ethanol.
Source	Synthetic.
Technical Info / Product Notes	Leupeptin gives multiple peaks on HPLC due to equilibria among three forms in solution. Purity determined using three main peaks. Majority of contaminating peptide is racemized leupeptin.

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-
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)