

Ac-Leu-Leu-Nle-CHO

Cysteine protease inhibitor

Cell-permeable, peptide aldehyde inhibitor of calpain I ($K_i=190\text{nM}$), calpain II ($K_i=150\text{nM}$), cathepsin L ($K_i=0.5\text{nM}$) and other neutral cysteine proteases. Inhibits cell cycle progression at G1/S and metaphase/anaphase in CHO cells by inhibiting cyclin B degradation. Also stimulates HMG-CoA synthase transcription by inhibiting degradation of active SREBP-1 (sterol regulatory element-binding protein 1). Protects against neuronal damage caused by hypoxia and ischemia. Inhibits apoptosis in thymocytes and metamyelocytes. Also prevents nitric oxide production by activated macrophages by interfering with the transcription of inducible nitric oxide synthase (iNOS; NOS II). Inhibits proteolytic degradation of I κ B α and I κ B β in RAW macrophages induced with LPS. It also prolong association of MHC class I molecules with the transporters associated with antigen processing.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P120-0005	5mg
BML-P120-0025	25mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Store dry and protect from light.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ALLN, Calpain inhibitor I, Ac-LL-norleucinal
Appearance	White powder.
CAS	110044-82-1
Formula	$C_{20}H_{37}N_3O_4$
MW	383.5
Purity	≥95% (HPLC)
Sequence	N-Acetyl-Leu-Leu-Nle-CHO
Solubility	Soluble in ethanol (5 mg/ml) or DMSO (10 mg/ml).
Source	Synthetic.



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)