Ac-Leu-Leu-NIe-CHO

Cysteine protease inhibitor

Cell-permeable, peptide aldehyde inhibitor of calpain I (K_i =190nM), calpain II (K_i =150nM), cathepsin L (K_i =0.5nM) 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 IkB α and IkB 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.

Last modified: May 29, 2024