MDL-28170

Calpain and cathepsin inhibitor

A potent and selective cell permeable calpain inhibitor. Inhibits oxidative damage-induced apoptosis in PC12 cells and capsaicin-induced apoptosis in dorsal root ganglion neurons. It protects against anoxia in rat optic nerves and reduces neuronal damage in an animal model of global cerebral ischemia(50 mg/kg). In isolated hippocampal pyramidal neurons, it inhibits Ca^{2+} -induced suppression of neurite outgrowth (complete inhibition at 20 µM). Calpain I K_i = 8 nM; Cathepsin B Ki = 24 nM in isolated enzyme assays. IC50 = 0.3 µM in intact cell assay.

Citations: 6

View Online »

Ordering Information

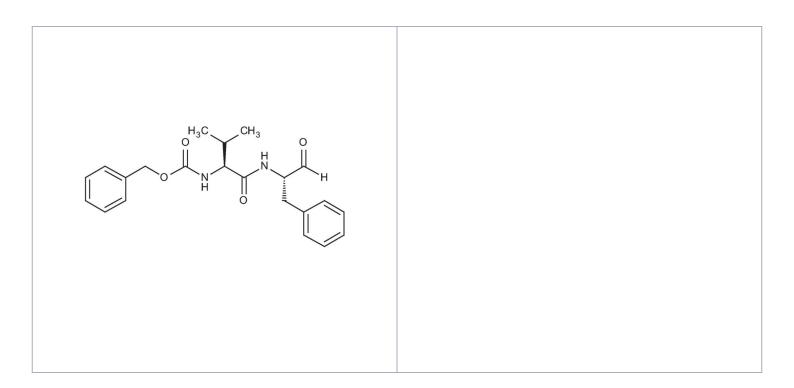
Order Online »

BML-PI130-0050

50mg

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-Benzyloxycarbonylvalylphenylalaninal

Appearance White solid.

CAS 88191-84-8

Formula $C_{22}H_{26}N_2O_4$

MW 382.5

Purity ≥98% (HPLC)

Sequence Z-Val-Phe-CHO

Solubility Soluble in DMSO (200mM).

Last modified: May 29, 2024