

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 K_i = 24 nM in isolated enzyme assays. IC_{50} = 0.3 μM in intact cell assay.

Citations: 6

[View Online »](#)

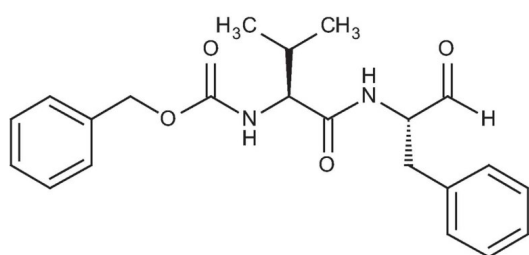
Ordering Information

[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-Benzylloxycarbonylvalylphenylalaninal

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).



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)