

Calpeptin

Calpain and cathepsin inhibitor

Membrane permeable inhibitor of calpain I and II, cathepsin L and cathepsin K ($IC_{50}=0.11$ nM). Promotes neurite elongation in differentiating PC12 cells. Prevents Ca^{2+} -ionophore induced degradation of actin binding protein and P235 in platelets. Inhibits myosin light chain phosphorylation in platelets stimulated by collagen, ionomycin (Prod. No. ALX-450-006) or thrombin. Inhibits LPS induced iNOS transcription and NF- κ B activation, possibly via proteasome inhibition.

Citations: 16

[View Online »](#)

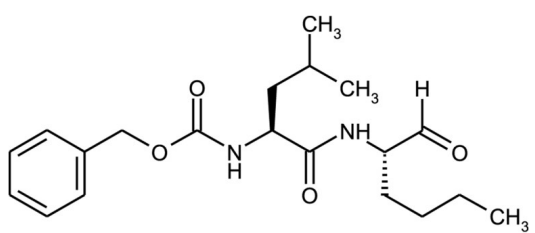
Ordering Information

[Order Online »](#)

| | |
|----------------|------|
| BML-PI101-0010 | 10mg |
| BML-PI101-0050 | 50mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|--------------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Stable for up to 1 year when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C. |
| Handling | Protect from light and moisture. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|---------------------------------------|---|
| Alternative Name | Z-Leu-norleucinal, N-Benzylloxycarbonyl-L-leucyl-norleucinal |
| Appearance | White to off-white powder. |
| CAS | 117591-20-5 |
| Couple Target | Calpain, Cathepsin, Proteasome |
| Couple Type | Inhibitor |
| Formula | $C_{20}H_{30}N_2O_4$ |
| MW | 362.5 |
| Purity | ≥98% (HPLC) |
| Solubility | Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml or dimethyl formamide (5mg/ml). |
| Technical Info / Product Notes | Concentrations for intact cells: 30μM (platelets), 10μM (myoblast cells) Note: Product is not sterile. |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com