

E-64-D

Calpain and cathepsin inhibitor

E-64-D is a cell-permeable ethyl ester of E-64-C (Prod. No. BML-PI106). The ester is hydrolyzed by intracellular esterases to produce E-64-C that inhibits calpain and cathepsins B, F, H, K, and L. E-64-D inhibits degradation of autophagic cargo inside autophagolysosomes, especially when used in combination with pepstatin (Prod. No. ALX-260-085). E-64-D combined with camostat mesylate also inhibits entry of SARS-CoV-2 in cells *in vitro*.

Citations: 41

[View Online »](#)

Ordering Information

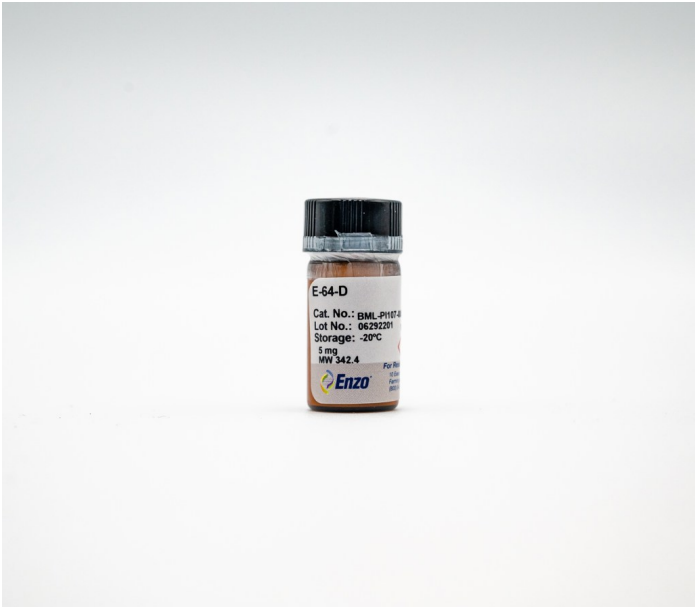
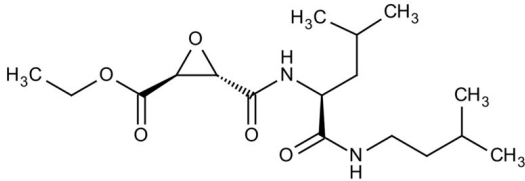
[Order Online »](#)

| | |
|----------------|-----|
| BML-PI107-0001 | 1mg |
| BML-PI107-0005 | 5mg |

Manuals, SDS & CofA

[View Online »](#)

- Potent inhibitor of calpain and cathepsins
- Used in autophagy research
- highly cited



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 2 months. |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|---|
| Alternative Name | 2S,3S-trans-(Ethoxycarbonyloxirane-2-carbonyl)-L-leucine-(3-methylbutyl) amide, Aloxistatin |
| Appearance | White solid. |
| CAS | 88321-09-9 |
| Couple Target | Calpain, Cathepsin, Protease |
| Couple Type | Inhibitor |
| Formula | $C_{17}H_{30}N_2O_5$ |
| Identity | Identity determined by MS. |
| MW | 342.4 |
| Purity | ≥98% (HPLC) |
| Solubility | Soluble in DMSO (20mg/ml) or 100% ethanol (>10mg/ml). |
| Source | Synthetic. |

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)