

Z-AAF-CMK

Negative control for AAF-CMK

Negative control for serine protease inhibitor AAF-CMK which inhibits tripeptidyl peptidase II (TPPII), a giant protease which may substitute for some proteasome functions. Blockage of the N-terminus by the Benzyloxycarbonyl or “Z” group abolishes the activity of the AAF-CMK inhibitor.

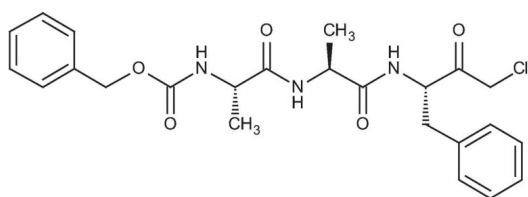
Ordering Information

[Order Online »](#)

BML-PI124-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



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 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Z-Ala-Ala-Phe-CMK
Appearance	Pale yellow solid.
MW	474.0
Purity	≥98% (TLC)
Sequence	Benzyloxycarbonyl – Ala-Ala-Phe-chloromethylketone
Solubility	Soluble in DMSO(2 mg/ml) or ethanol (2 mg/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-
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)