

# Z-AAD-CMK

## Caspase inhibitor

Irreversible, cell-permeable inhibitor of granzyme B ( $ID_{50}=0.3\text{ }\mu\text{M}$ ). Does not inhibit granzyme A. Inhibition of granzyme B with Z-AAD-CMK prevented caspase-3 activation in T cells.

Citations: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-P165-0001	1mg
---------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Granzyme B inhibitor
<b>Appearance</b>	White to off-white solid.
<b>Formula</b>	$C_{19}H_{24}ClN_3O_7$
<b>Formulation</b>	Lyophilized.
<b>MW</b>	441.9
<b>Purity</b>	≥95% (HPLC)
<b>Sequence</b>	Benzyloxycarbonyl-Ala-Ala-Asp-chloromethylketone
<b>Solubility</b>	Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml) or water (20mg/ml).
<b>Source</b>	Synthetic.



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)