Z-AAD-CMK

Caspase inhibitor

Irreversible, cell-permeable inhibitor of granzyme B (ID $_{50}$ =0.3 μ 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

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.

Handling 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 Granzyme B inhibitor

Appearance White to off-white solid.

Formula $C_{19}H_{24}CIN_3O_7$

Formulation Lyophilized.

MW 441.9

Purity ≥95% (HPLC)

Sequence Benzyloxycarbonyl-Ala-Ala-Asp-chloromethylketone

Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml) or

water (20mg/ml).

Source Synthetic.

Last modified: May 29, 2024



 ENZO LIFE SCIENCES,
 EL

 INC.
 EI

 Phone: 800.942.0430
 (E

 info Pl

 usa@enzolifesciences.com
 inf

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com