

PRODUCT INFORMATION



Z-VAD(OMe)-FMK

Item No. 14463

CAS Registry No.: 187389-52-2
Formal Name: N-[(phenylmethoxy)carbonyl]-

Synonym: Z-Val-Ala-Asp-(OMe)-
Fluoromethyl Ketone

MF: C₂₂H₃₀FN₃O₇

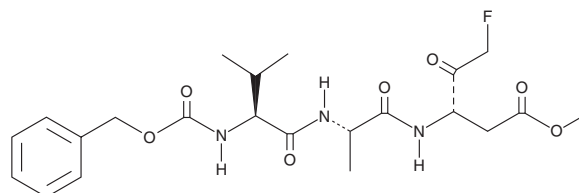
FW: 467.5

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Z-VAD(OMe)-FMK is supplied as a crystalline solid. A stock solution may be made by dissolving the Z-VAD(OMe)-FMK in the solvent of choice. Z-VAD(OMe)-FMK is soluble in DMSO at a concentration of approximately 5 mg/ml.

Z-VAD(OMe)-FMK is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Z-VAD(OMe)-FMK is a cell-permeable, competitive, and irreversible inhibitor of all caspases.^{1,2} Through this action, it inhibits cleavage of poly(ADP-ribose) polymerase, preventing apoptosis when used at 10-50 μM.^{1,3} It also blocks caspase-mediated apoptosis *in vivo*.⁴ Z-VAD(OMe)-FMK effectively prevents caspase action in inflammasomes.⁵

References

1. Xiang, J., Chao, D.T., and Korsmeyer, S.J. BAX-induced cell death may not require interleukin 1β-converting enzyme-like proteases. *Proc. Natl. Acad. Sci. USA* **93**, 14559-14563 (1996).
2. Garcia-Calvo, M., Peterson, E.P., Leiting, B., et al. Inhibition of human caspases by peptide-based and macromolecular inhibitors. *J. Biol. Chem.* **273**, 32608-32613 (1998).
3. Slee, E.A., Zhu, H., Chow, S.C., et al. Benzylloxycarbonyl-Val-Ala-Asp (OMe) fluoromethylketone (Z-VAD-FMK) inhibits apoptosis by blocking the processing of CPP32. *Biochem. J.* **315**, 21-24 (1996).
4. Künstle, G., Leist, M., Uhlig, S., et al. ICE-protease inhibitors block murine liver injury and apoptosis caused by CD95 or by TNF-α. *Immunol. Lett.* **55(1)**, 5-10 (1997).
5. Meraz, I.M., Melendez, B., Gu, J., et al. Activation of the inflammasome and enhanced migration of microparticle-stimulated dendritic cells to the draining lymph node. *Mol. Pharm.* **9**, 2049-2062 (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/05/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM