

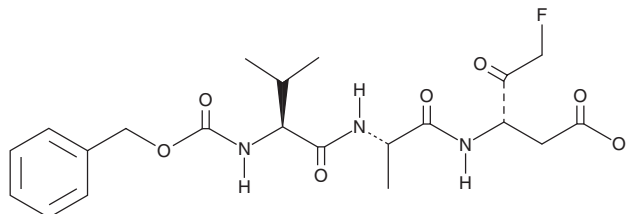
# PRODUCT INFORMATION



## Z-VAD(OH)-FMK

Item No. 14467

**CAS Registry No.:** 161401-82-7  
**Formal Name:** N-[(phenylmethoxy)carbonyl]-L-valyl-L-[[1S]-1-(carboxymethyl)-3-fluoro-2-oxopropyl]-L-alaninamide  
**Synonym:** Z-Val-Ala-Asp-(OH)-Fluoromethyl Ketone  
**MF:** C<sub>21</sub>H<sub>28</sub>FN<sub>3</sub>O<sub>7</sub>  
**FW:** 453.5  
**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(OH)-FMK is supplied as a crystalline solid. A stock solution may be made by dissolving the Z-VAD(OH)-FMK in the solvent of choice, which should be purged with an inert gas. Z-VAD(OH)-FMK is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

### Description

Z-VAD(OH)-FMK is an irreversible tripeptide inhibitor of all caspases.<sup>1-3</sup> This non-methylated form is useful in studies involving recombinant, isolated, or purified enzymes. Unlike the methylated form, Z-VAD(OMe)-FMK (Item No. 14463), this compound does not require pretreatment with esterases prior to *in vitro* use.

### References

1. Cain, K., Inayat-Hussain, S.H., Couet, C., *et al.* A cleavage-site-directed inhibitor of interleukin-1 $\beta$ -converting enzyme-like proteases inhibits apoptosis in primary cultures of rat hepatocytes. *Biochem. J.* **314**(Pt 1), 27-32 (1996).
2. Jacobson, M.D., Weil, M., and Raff, M.C. Role of Ced-3/ICE-family proteases in staurosporine-induced programmed cell death. *J. Cell Biol.* **133**(5), 1041-1051 (1996).
3. Garcia-Calvo, M., Peterson, E.P., Leiting, B., *et al.* Inhibition of human caspases by peptide-based and macromolecular inhibitors. *J. Biol. Chem.* **273**(49), 32608-32613 (1998).

#### 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, 10/10/2022

#### 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