

PRODUCT INFORMATION



Z-VA-DL-D(OMe)-FMK (trifluoroacetate salt)

Item No. 27421

Formal Name: N-[(phenylmethoxy)carbonyl]-L-valyl-N-[3-fluoro-1-(2-methoxy-2-oxoethyl)-2-oxopropyl]-L-alaninamide, trifluoroacetate salt

Synonyms: Caspase-1 Inhibitor V, Z-VAD-FMK, Z-VAD(OMe)-FMK, Z-Val-Ala-DL-Asp(OMe)-Fluoromethyl Ketone

MF: $C_{22}H_{30}FN_3O_7 \cdot XCF_3COOH$

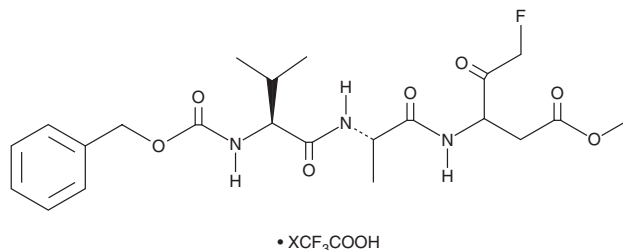
FW: 467.5

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Z-VA-DL-D(OMe)-FMK (Z-VAD-FMK) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Z-VAD-FMK (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Z-VAD-FMK (trifluoroacetate salt) is soluble in organic solvents such as DMSO and formic acid. The solubility of Z-VAD-FMK (trifluoroacetate salt) in these solvents is approximately 1 mg/ml.

Description

Z-VAD-FMK is a pan-caspase inhibitor.¹ It inhibits anti-Fas antibody-induced cleavage of caspase-3 and PARP in Jurkat T lymphocytes when used at a concentration of 0.5 μM . Z-VAD-FMK (50 μM) inhibits cell death induced by TNF- α and cycloheximide co-application in HeLa cells. It also inhibits radiation-induced increases in the number of TUNEL-positive neurons in the rat hypoglossal nucleus when administered intracerebroventricularly at a dose of 2 μg per animal.²

References

1. Yang, W., Guastella, J., Huang, J.-C., *et al.* MX1013, a dipeptide caspase inhibitor with potent *in vivo* antiapoptotic activity. *Br. J. Pharmacol.* **140**(2), 402-412 (2003).
2. Li, J., Wang, Y., Du, L., *et al.* Radiation-induced cytochrome c release and the neuroprotective effects of the pan-caspase inhibitor z-VAD-fmk in the hypoglossal nucleus. *Exp. Ther. Med.* **7**(2), 383-388 (2014).

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, 12/15/2022

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