

PRODUCT INFORMATION



Z-DEVD-FMK

Item No. 14414

CAS Registry No.: 210344-95-9

Formal Name: N-[(phenylmethoxy)carbonyl]-L- α -aspartyl-L- α -glutamyl-N-[(1S)-3-fluoro-1-(2-methoxy-2-oxoethyl)-2-oxopropyl]-L-valinamide, 1,2-dimethyl ester

MF: $C_{30}H_{41}FN_4O_{12}$

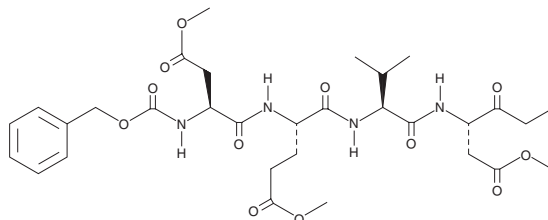
FW: 668.7

Purity: $\geq 95\%$

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-DEVD-FMK is supplied as a crystalline solid. A stock solution may be made by dissolving the Z-DEVD-FMK in the solvent of choice, which should be purged with an inert gas. Z-DEVD-FMK is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

Caspase-3 is a cysteinyl aspartate-specific protease which plays a central role in apoptosis.¹ Z-DEVD-FMK is a cell-permeable irreversible inhibitor of caspase-3 ($\text{IC}_{50} = 130 \text{ nM}$).^{2,3} It potently inhibits apoptosis.⁴

References

1. Granville, D.J., Carthy, C.M., Jiang, H., *et al.* Rapid cytochrome c release, activation of caspases 3, 6, 7, and 8 followed by Bap31 cleavage in HeLa cells treated with photodynamic therapy. *FEBS Lett.* **437**, 5-10 (1998).
2. Brockstedt, E., Rickers, A., Kostka, S., *et al.* Identification of apoptosis-associated proteins in a human Burkitt lymphoma cell line. Cleavage of heterogeneous nuclear ribonucleoprotein A1 by caspase 3. *J. Biol. Chem.* **273**(43), 28057-28064 (1998).
3. Lee, J. and Bogoy, M. Development of near-infrared fluorophore (NIRF)-labeled activity-based probes for *in vivo* imaging of legumain. *ACS Chem. Biol.* **5**(2), 233-243 (2010).
4. Kugawa, F., Ueno, A., and Aoki, M. Apoptosis of NG108-15 cells induced by buprenorphine hydrochloride occurs via the caspase-3 pathway. *Biol. Pharm. Bull.* **23**(8), 930-935 (2000).

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, 09/29/2022

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