

PRODUCT INFORMATION



Z-VEID-FMK

Item No. 44015

CAS Registry No.: 210344-96-0
Formal Name: methyl (5S,8S,11S,14S)-11-((S)-sec-butyl)-14-(2-fluoroacetyl)-5-isopropyl-8-(3-methoxy-3-oxopropyl)-3,6,9,12-tetraoxo-1-phenyl-2-oxa-4,7,10,13-tetraazahexadecan-16-oate

Synonym: Cbz-Val-Glu(OMe)-Ile-Asp(OMe)-CH₂F

Peptide Sequence: ZVXIJ-CH₂F (X = E(OMe), J = D(OMe))

MF: C₃₁H₄₅FN₄O₁₀

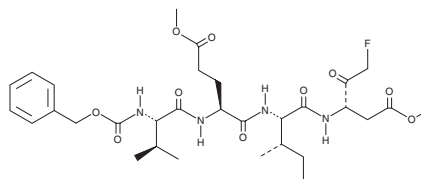
FW: 652.7

Purity: ≥98%

Supplied as: A 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-VEID-FMK is supplied as a solid. A stock solution may be made by dissolving the Z-VEID-FMK in the solvent of choice, which should be purged with an inert gas. Z-VEID-FMK is soluble in DMSO.

Description

Z-VEID-FMK is a peptide inhibitor of caspase-3 and -6 (apparent IC₅₀s = 45.1 and 128.6 nM, respectively).¹ It decreases lamin cleavage induced by caspase-6 in a cell-free assay (IC₅₀ = 45 μM). Z-VEID-FMK (160 μM) decreases apoptosis in primary human erythroblasts cultured in erythropoietin-deficient media. It increases survival of primary rat retinal whole-mount cultures when used at a concentration of 10 μM.² Intraocular administration of Z-VEID-FMK decreases apoptosis in retinal ganglion cells and increases axonal regeneration in a rat model of optic nerve crush.

References

1. Mintzer, R., Ramaswamy, S., Shah, K., *et al.* A whole cell assay to measure caspase-6 activity by detecting cleavage of lamin A/C. *PLoS One* **7(1)**, e30376 (2012).
2. Monnier, P.P., D'Onofrio, P.M., Magharious, M., *et al.* Involvement of caspase-6 and caspase-8 in neuronal apoptosis and the regenerative failure of injured retinal ganglion cells. *J. Neurosci.* **31(29)**, 10494-10505 (2011).

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, 08/15/2025

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