

# PRODUCT INFORMATION



## Z-AEVD-FMK

Item No. 14987

CAS Registry No.: 1135688-47-9

Formal Name: methyl-14S-(2-fluoroacetyl)-11S-isopropyl-8S-(3-methoxy-3-oxopropyl)-5S-methyl-3,6,9,12-tetraoxo-1-phenyl-2-oxa-4,7,10,13-tetraazahexadecan-16-oate

Synonym: Z-Ala-Glu-Val-Asp-Fluoromethyl Ketone

MF:  $C_{28}H_{39}FN_4O_{10}$

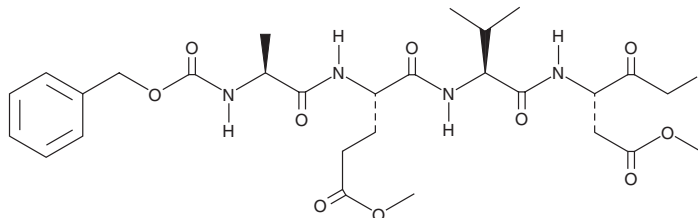
FW: 610.6

Purity:  $\geq 95\%$

Supplied as: A crystalline solid

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

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Z-AEVD-FMK is supplied as a crystalline solid. A stock solution may be made by dissolving the Z-AEVD-FMK in the solvent of choice. Z-AEVD-FMK is soluble in DMSO at a concentration of approximately 10 mg/ml.

### Description

Z-AEVD-FMK is an irreversible inhibitor of caspase-10 and related caspases.<sup>1</sup> At 10  $\mu\text{M}$ , it can prevent the initiation of Fas signaling by caspase-10 in Jurkat T lymphoma cells, preventing Bid cleavage into its active form, caspase cascade activation, and apoptosis.<sup>2</sup>

### References

1. Park, S.-J., Wu, C.-H., Gordon, J.D., *et al.* Taxol induces caspase-10-dependent apoptosis. *J. Biol. Chem.* **279**(49), 51057-51067 (2004).
2. Milhas, D., Cuvillier, O., Therville, N., *et al.* Caspase-10 triggers Bid cleavage and caspase cascade activation in FasL-induced apoptosis. *J. Biol. Chem.* **280**(20), 19836-19842 (2005).

#### 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, 11/15/2022

#### 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