

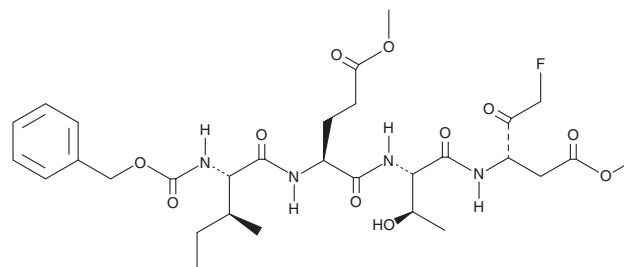
# PRODUCT INFORMATION



## Z-IETD-FMK

Item No. 42519

**CAS Registry No.:** 210344-98-2  
**Formal Name:** N-[(phenylmethoxy)carbonyl]-L-isoleucyl-L- $\alpha$ -glutamyl-L-[(1S)-3-fluoro-1-(2-methoxy-2-oxoethyl)-2-oxopropyl]-L-threoninamide, methyl ester  
**Synonyms:** Cbz-Ile-Glu(OMe)-Thr-Asp(OMe)-CH<sub>2</sub>F  
**Peptide Sequence:** ZIE(OMe)TD(OMe)-CH<sub>2</sub>F  
**MF:** C<sub>30</sub>H<sub>43</sub>FN<sub>4</sub>O<sub>11</sub>  
**FW:** 654.7  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Z-IETD-FMK is supplied as a solid. A stock solution may be made by dissolving the Z-IETD-FMK in the solvent of choice, which should be purged with an inert gas. Z-IETD-FMK is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

### Description

Z-IETD-FMK is a peptide inhibitor of caspase-8.<sup>1</sup> It inhibits TNF- $\alpha$ -induced apoptosis in isolated human neutrophils when used at a concentration of 3 or 10  $\mu$ M. Z-IETD-FMK (100  $\mu$ M) decreases human rhinovirus (HRV) replication and HRV-induced apoptosis in HeLa cervical cancer cells.<sup>2</sup> Intrathecal administration of Z-IETD-FMK (0.5  $\mu$ g/animal) reduces tumor growth and increases tumor macrophage and T cell infiltration in a mouse model of lung cancer induced by N-methyl-N-nitrosourea (NMU).<sup>3</sup>

### References

1. Cowburn, A.S., White, J.F., Deighton, J., *et al.* z-VAD-fmk augmentation of TNF $\alpha$ -stimulated neutrophil apoptosis is compound specific and does not involve the generation of reactive oxygen species. *Blood* **105**(7), 2970-2972 (2005).
2. Deszcz, L., Seipelt, J., Vassilieva, E., *et al.* Antiviral activity of caspase inhibitors: Effect on picornaviral 2A proteinase. *FEBS Lett.* **560**(1-3), 51-55 (2004).
3. Terlizzi, M., Di Crescenzo, V.G., Perillo, G., *et al.* Pharmacological inhibition of caspase-8 limits lung tumour outgrowth. *Br. J. Pharmacol.* **172**(15), 3917-3928 (2015).

#### 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, 6/26/2025

#### 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