

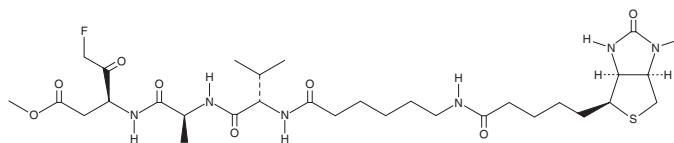
PRODUCT INFORMATION



Biotin-VAD-FMK

Item No. 24305

CAS Registry No.: 1135688-15-1
Formal Name: methyl (3S,6S,9S)-3-(2-fluoroacetyl)-9-isopropyl-6-methyl-5,8,11,18-tetraoxo-22-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)-4,7,10,17-tetraazadocosanoate
MF: C₃₀H₄₉FN₆O₈S
FW: 672.8
Purity: ≥95%
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

Biotin-VAD-FMK is supplied as a solid. A stock solution may be made by dissolving the biotin-VAD-FMK in the solvent of choice. Biotin-VAD-FMK is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 20 mM.

Description

Biotin-VAD-FMK is a biotin-conjugated form of the cell-permeable caspase inhibitor Z-VAD(OMe)-FMK (Item No. 14463).¹ It is an affinity probe that allows for detection and immobilization of caspases using the biotin ligand. Biotin-VAD-FMK inhibits caspase-3-like activity in a dose-dependent manner in hepatocytes stimulated with TNF- α and ActD. It also binds to caspase-3 and caspase-8 in hepatocyte lysates, an effect that can be blocked by the caspase inhibitor Ac-DEVD-CHO (Item No. 10017).

Reference

1. Li, J., Bombeck, C.A., Yang, S., *et al.* Nitric oxide suppresses apoptosis via interrupting caspase activation and mitochondrial dysfunction in cultured hepatocytes. *J. Biol. Chem.* **274**(24), 17325-17333 (1999).

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, 05/07/2024

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