

PRODUCT INFORMATION



Ac-FLTD-CMK (trifluoroacetate salt)

Item No. 34469

Formal Name: (4S,7S,10S,13S)-4-benzyl-13-(2-chloroacetyl)-10-((R)-1-hydroxyethyl)-7-isobutyl-2,5,8,11-tetraoxo-3,6,9,12-tetraazapentadecan-15-oic acid, trifluoroacetate salt

Synonym: Ac-Phe-Leu-Thr-Asp-CMK

MF: C₂₆H₃₇ClN₄O₈ • XCF₃COOH

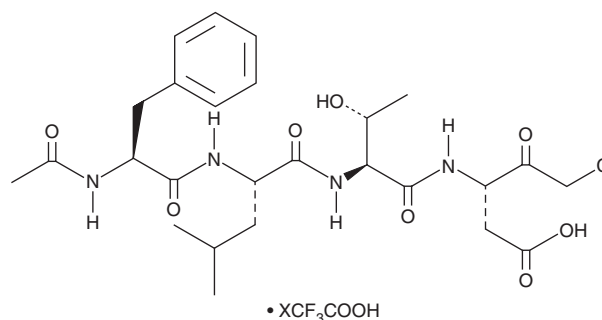
FW: 569.0

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

Ac-FLTD-CMK (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-FLTD-CMK (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-FLTD-CMK (trifluoroacetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

Ac-FLTD-CMK is an inhibitor of caspase-1, -4, -5, and -11.¹ It inhibits caspase-1, -4, and -5 (IC₅₀s = 0.0467, 1.49, and 0.329 μM, respectively, for the recombinant human enzymes) and cleavage of gasdermin D (GSDMD), an effector involved in pyroptosis and NETosis, by mouse caspase-11 when used at a concentration of 10 μM. Ac-FLTD-CMK (10 μM) inhibits LPS- and ATP- or LPS- and nigericin-induced pyroptosis and IL-1β release in mouse bone marrow-derived macrophages (BMDMs).

Reference

1. Yang, J., Liu, Z., Wang, C., *et al.* Mechanism of gasdermin D recognition by inflammatory caspases and their inhibition by a gasdermin D-derived peptide inhibitor. *Proc. Natl. Acad. Sci. USA* **115**(26), 6792-6797 (2018).

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, 12/22/2022

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