

PRODUCT INFORMATION



Ac-LEVD-CHO (trifluoroacetate salt)

Item No. 27433

Formal Name: N-acetyl-L-leucyl-L- α -glutamyl-N-[(1S)-2-carboxy-1-formylethyl]-L-valinamide, trifluoroacetate salt

Synonyms: Ac-Leu-Glu-Val-Asp-CHO, Caspase-4 Inhibitor I

MF: C₂₂H₃₆N₄O₉ • XCF₃COOH

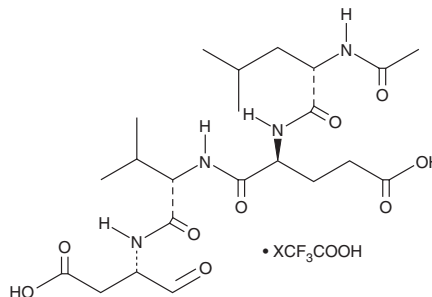
FW: 500.5

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

Description

Ac-LEVD-CHO is a caspase-4 inhibitor.¹ It inhibits IL-1 α expression and secretion and caspase-4 activation induced by the *T. denticola* periodontal pathogen surface protein Td92 in human gingival fibroblasts when used at a concentration of 30 μ M. Ac-LEVD-CHO (10 μ M) also reduces apoptosis induced by expression of the dominant negative adenoviral RNA-dependent protein kinase (Ad- Δ 6PKR) in A549 and PC3 cancer cells.²

References

1. Jun, H.-K., Jung, Y.-J., An, S.-J., *et al.* Caspase-4 activation by a bacterial surface protein is mediated by cathepsin G in human gingival fibroblasts. *Cell Death Differ.* **25(2)**, 380-391 (2018).
2. Pataer, A., Swisher, S.G., Roth, J.A., *et al.* Inhibition of RNA-dependent protein kinase (PKR) leads to cancer cell death and increases chemosensitivity. *Cancer Biol. Ther.* **8(3)**, 245-252 (2009).

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, 12/14/2022

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