

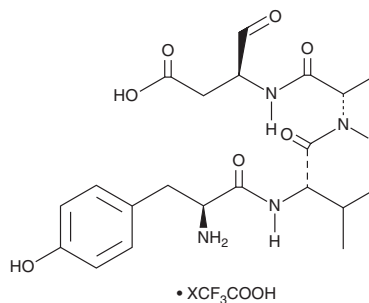
PRODUCT INFORMATION



YVAD-CHO (trifluoroacetate salt)

Item No. 27420

Formal Name: L-tyrosyl-L-valyl-N-[(1S)-2-carboxy-1-formylethyl]-L-alaninamide, trifluoroacetate salt
Synonym: H-Tyr-Val-Ala-Asp-CHO
MF: C₂₁H₃₀N₄O₇ • XCF₃COOH
FW: 450.5
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

YVAD-CHO (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the YVAD-CHO (trifluoroacetate salt) in water. The solubility of YVAD-CHO (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

YVAD-CHO is an inhibitor of caspase-1/interleukin-1 β converting enzyme (ICE) and a non-acetylated form of Ac-YVAD-CHO (Item No. 10016).¹ It prevents ICE-like and CPP32-like proteinase activity and inhibits apoptosis induced by ceramide or irinotecan (CPT-11; Item No. 14180) in 4B1 cells when used at a concentration of 600 μ M. YVAD-CHO (100 μ M) also prevents chromogranin A-induced cell death of primary rat microglia but not primary rat cerebellar granule neurons.²

References

1. Suzuki, A., Iwasaki, M., Kato, M., *et al.* Sequential operation of ceramide synthesis and ICE cascade in CPT-11-initiated apoptotic death signaling. *Exp. Cell Res.* **233(1)**, 41-47 (1997).
2. Kingham, P.J., Cuzner, M.L., and Pocock, J.M. Apoptotic pathways mobilized in microglia and neurones as a consequence of chromogranin A-induced microglial activation. *J. Neurochem.* **73(2)**, 538-547 (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, 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