

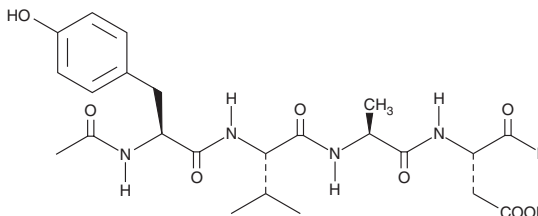
PRODUCT INFORMATION



Ac-YVAD-CHO

Item No. 10016

CAS Registry No.: 143313-51-3
Formal Name: N-acetyl-L-tyrosyl-L-valyl-N-
[(1S)-2-carboxyl-1-formylethyl]-L-
alaninamide
Synonym: N-Ac-Tyr-Val-Ala-Asp-CHO
MF: C₂₃H₃₂N₄O₈
FW: 492.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that Ac-YVAD-CHO be stored as supplied at -20°C. It should be stable for at least two years.

Ac-YVAD-CHO is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-YVAD-CHO in an organic solvent purged with an inert gas. Ac-YVAD-CHO is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Ac-YVAD-CHO in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Ac-YVAD-CHO can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of Ac-YVAD-CHO in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-YVAD-CHO is a selective inhibitor of interleukin-1 β converting enzyme (ICE; Caspase 1).¹ Caspase inhibitors interfere with the initiation of apoptosis and have numerous possible clinical applications.²

References

1. Thornberry, N.A., Peterson, E.P., Zhao, J.J., *et al.* Inactivation of interleukin-1 β converting enzyme by peptide (Acyloxy)methyl ketones. *Biochemistry* **33**, 3934-3940 (1994).
2. Nicholson, D.W., Ali, A., Thornberry, N.A., *et al.* Identification and inhibition of the ICE/CED-3 protease necessary for mammalian apoptosis. *Nature* **376**, 37-43 (1995).

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, 07/28/2015

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