

# PRODUCT INFORMATION



## Ac-YVAD-CMK

Item No. 10014

**CAS Registry No.:** 178603-78-6  
**Formal Name:** N-acetyl-L-tyrosyl-L-valyl-N-[(1S)-1-(carboxymethyl)-3-chloro-2-oxo-propyl]-L-alaninamide

**Synonym:** N-Ac-Tyr-Val-Ala-Asp-CMK

**MF:** C<sub>24</sub>H<sub>33</sub>ClN<sub>4</sub>O<sub>8</sub>

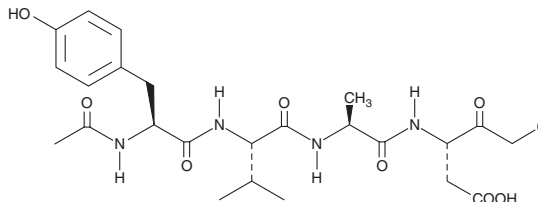
**FW:** 541.0

**Purity:** ≥98%

**Supplied as:** A crystalline 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-YVAD-CMK is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-YVAD-CMK in the solvent of choice, which should be purged with an inert gas. Ac-YVAD-CMK is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Ac-YVAD-CMK in these solvents is approximately 20 and 10 mg/ml, respectively.

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-CMK can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Ac-YVAD-CMK in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Ac-YVAD-CMK is a selective, irreversible inhibitor of interleukin-1 $\beta$  converting enzyme (ICE; Caspase 1). Ac-YVAD-CMK is neuroprotective in a rat model of cerebral ischemia.<sup>1</sup>

### Reference

1. Rabuffetti, M., Sciorati, C., Tarozzo, G., *et al.* Inhibition of caspase-1-like activity by Ac-Tyr-Val-Ala-Asp-chloromethyl ketone induces long-lasting neuroprotection in cerebral ischemia through apoptosis reduction and decrease of proinflammatory cytokines. *J. Neurosci.* **20**, 4398-4404 (2000).

#### 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, 09/28/2022

#### 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