

PRODUCT INFORMATION



Z-VAD-AMC (acetate)

Item No. 27142

Formal Name: (5S,8S,11S)-5-isopropyl-8-methyl-11-((4-methyl-2-oxo-2H-chromen-7-yl)carbamoyl)-3,6,9-trioxo-1-phenyl-2-oxa-4,7,10-triazatridecan-13-oic acid, monoacetate

Synonyms: Z-Val-Ala-Asp-AMC,
Z-Val-Ala-Asp-7-amino-4-methylcoumarin

MF: C₃₀H₃₄N₄O₉ • XCH₃COOH

FW: 594.6

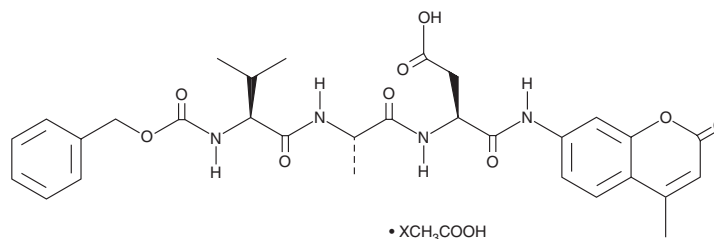
Purity: ≥95%

EX./Em. Max: 340-360/440-460 nm

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

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

Description

Z-VAD-AMC is a fluorogenic substrate for caspase-1.¹ Upon enzymatic cleavage by caspase-1, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify caspase-1 activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Küntzle, G., Leist, M., Uhlig, S., *et al.* ICE-protease inhibitors block murine liver injury and apoptosis caused by CD95 or by TNF- α . *Immunol. Lett.* **55**(1), 5-10 (1997).

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, 12/19/2022

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