

PRODUCT INFORMATION



Suc-YVAD-AMC (acetate)

Item No. 27138

Formal Name: (3S,6S,9S,12S)-12-(4-hydroxybenzyl)-9-isopropyl-6-methyl-3-((4-methyl-2-oxo-2H-chromen-7-yl)carbamoyl)-5,8,11,14-tetraoxo-4,7,10,13-tetraazaheptadecanedioic acid, acetate

Synonyms: Suc-Tyr-Val-Ala-Asp-AMC, Suc-Tyr-Val-Ala-Asp-7-amino-4-methylcoumarin, Suc-YVAD-NH-Mec

MF: $C_{35}H_{41}N_5O_{12} \cdot XC_2H_4O_2$

FW: 723.7

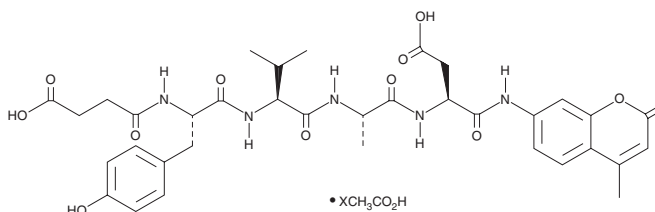
Purity: $\geq 95\%$

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

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

Description

Suc-YVAD-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. Chakravarthy, B.R., Walker, T., Rasquinha, I., *et al.* Activation of DNA-dependent protein kinase may play a role in apoptosis of human neuroblastoma cells. *J. Neurochem.* **72**(3), 933-942 (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, 11/04/2022

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