

PRODUCT INFORMATION



Ac-VEID-AMC (ammonium acetate salt)

Item No. 14989

Formal Name: N-acetyl-L-valyl-L- α -glutamyl-L-isoleucyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L- α -asparagine, ammonium acetate salt

Synonyms: N-Acetyl-Val-Glu-Ile-Asp-7-amido-4-Methylcoumarin, Caspase-6 Substrate (Fluorogenic)

MF: C₃₂H₄₃N₅O₁₁ • XC₂H₇NO₂
FW: 673.7

Purity: ≥95%

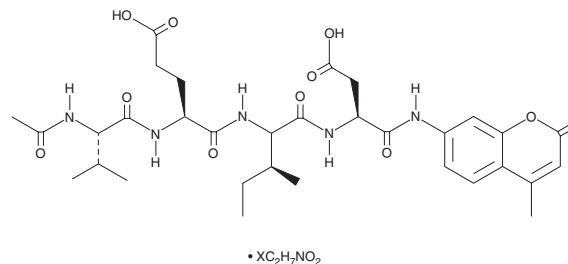
UV/Vis.: λ_{max} : 328 nm

Ex./Em. Max: 328 nm

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-VEID-AMC (ammonium acetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-VEID-AMC (ammonium acetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-VEID-AMC (ammonium acetate salt) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

Ac-VEID-AMC is a fluorogenic substrate based on the caspase-6 cleavage site in lamin A at amino acids VEID during apoptosis.¹ It has also been reported to be cleaved by related proteases, including caspase-8.² Caspase activity can be quantified by fluorescent detection of free AMC (also known as 7-amino-4-methylcoumarin), which is excited at 340-360 nm and emits at 440-460 nm.

References

1. Talanian, R.V., Quinlan, C., Trautz, S., et al. Substrate specificities of caspase family proteases. *J. Biol. Chem.* **272**(15), 9677-9682 (1997).
2. Chae, H.-J., Park, K.-M., Lee, G.-Y., et al. Je-Chun-Jun induced apoptosis of human cervical carcinoma HeLa cells. *Acta Pharmacol. Sin.* **25**(10), 1372-1379 (2004).

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, 11/07/2022

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