

PRODUCT INFORMATION



Ac-DEVD-AMC

Item No. 14986

CAS Registry No.: 169332-61-0

Formal Name: N-acetyl-L- α -aspartyl-L- α -glutamyl-L-valyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L- α -asparagine

Synonyms: N-Acetyl-Asp-Glu-Val-Asp-7-amido-4-Methylcoumarin, Caspase-3 Substrate (Fluorogenic)

MF: C₃₀H₃₇N₅O₁₃

FW: 675.6

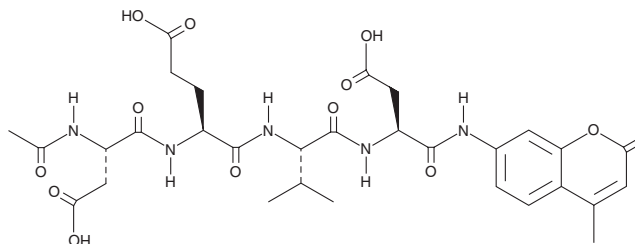
Purity: $\geq 95\%$

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

Supplied as: A solution in DMSO

Storage: -20°C

Stability: ≥ 1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ac-DEVD-AMC is supplied as a solution in DMSO. The solubility of Ac-DEVD-AMC in DMSO is approximately 10 mg/ml.

Description

Ac-DEVD-AMC is a fluorogenic substrate that can be cleaved specifically by caspase-3.¹ 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.

Reference

1. Kravchenko, D.V., Kysil, V.V., Ilyn, A.P., *et al.* 1,3-Dioxo-4-methyl-2,3-dihydro-1H-pyrrolo[3,4-c]quinolines as potent caspase-3 inhibitors. *Bioorg. Med. Chem. Lett.* **15**, 1841-1845 (2005).

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, 12/20/2017

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