

PRODUCT INFORMATION



Ac-DNLD-AMC

Item No. 24566

CAS Registry No.: 958001-92-8

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

Synonyms: Ac-Asp-Asn-Leu-Asp-MCA, N-Acetyl-Asp-Asn-Leu-Asp-7-amido-Methylcoumarin, Caspase-3 Substrate

MF: $C_{30}H_{38}N_6O_{12}$

FW: 674.7

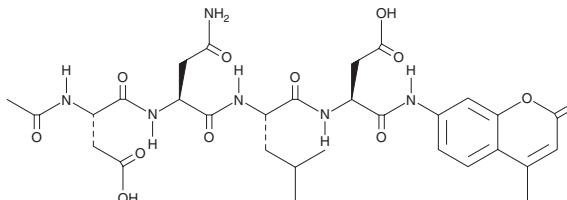
Purity: $\geq 95\%$

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

Supplied as: A lyophilized powder

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Ac-DNLD-AMC is supplied as a lyophilized powder. A stock solution may be made by dissolving the Ac-DNLD-AMC in the solvent of choice, which should be purged with an inert gas. Ac-DNLD-AMC is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Ac-DNLD-AMC is a fluorogenic caspase-3 substrate.¹ Upon enzymatic cleavage by caspase-3, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify caspase-3 activity. AMC displays excitation/emission maxima of 340-360 and 440-460 nm, respectively.

Reference

1. Yoshimori, A., Sakai, J., Sunaga, S., *et al.* Structural and functional definition of the specificity of a novel caspase-3 inhibitor, Ac-DNLD-CHO. *BMC Pharmacol.* 7, 8 (2007).

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, 12/17/2022

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