

PRODUCT INFORMATION



Ac-LEHD-AMC (trifluoroacetate salt)

Item No. 24569

Formal Name: N-acetyl-L-leucyl-L- α -glutamyl-L-histidyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L- α -asparagine, trifluoroacetate salt

Synonyms: Ac-Leu-Glu-His-Asp-AMC, Caspase-9 Substrate

MF: $C_{33}H_{41}N_7O_{11} \cdot XCF_3COOH$

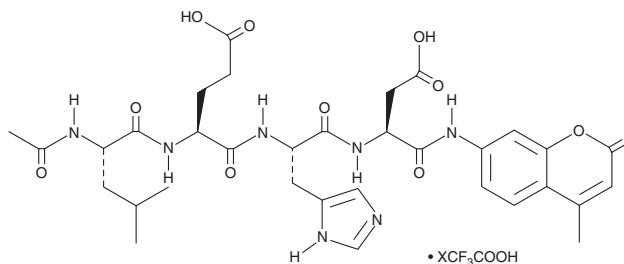
FW: 711.7

Purity: $\geq 95\%$

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-LEHD-AMC (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Ac-LEHD-AMC (trifluoroacetate salt) in water. The solubility of Ac-LEHD-AMC (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-LEHD-AMC is a fluorogenic substrate for caspase-9.¹ Upon cleavage by caspase-9, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify caspase-9 activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Prudnikov, I.M. and Smirnov, A.N. Short peptide tools for monitoring caspase and proteasome activities in embryonal and adult rat brain lysates: An approach for the differential identification of proteases. *J. Biochem.* **151**(3), 299-316 (2012).

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, 10/17/2022

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