

PRODUCT INFORMATION



Ac-LEHD-pNA (trifluoroacetate salt)

Item No. 24570

Formal Name: (2S,5S,8S,11S)-5-((1H-imidazol-5-yl)methyl)-8-(2-carboxyethyl)-11-isobutyl-2-(2-((4-nitrophenyl)amino)-2-oxoethyl)-4,7,10,13-tetraoxo-3,6,9,12-tetraazatetradecanoic acid, 2,2,2-trifluoroacetate

Synonyms: Ac-Leu-Glu-His-Asp-pNA, Caspase-9 Chromogenic Substrate I

MF: C₂₉H₃₈N₈O₁₁

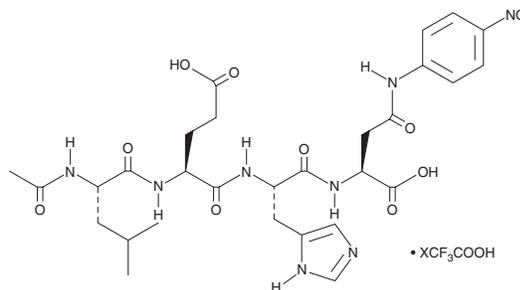
FW: 674.7

Purity: ≥95%

Supplied as: A lyophilized powder

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-LEHD-pNA (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Ac-LEHD-pNA (trifluoroacetate salt) in water. The solubility of Ac-LEHD-pNA (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-pNA is a substrate for caspase-9.¹ Caspase-9 preferentially binds to and cleaves the Leu-Glu-His-Asp (LEHD) peptide sequence to release p-nitroalanine, which can be quantified by colorimetric detection at 405 nm as a measure of enzyme activity.¹⁻³

References

1. Hsu, Y.L., Kuo, P.L., Lin, L.T., *et al.* Asiatic acid, a triterpene, induces apoptosis and cell cycle arrest through activation of extracellular signal-regulated kinase and p38 mitogen-activated protein kinase pathways in human breast cancer cells. *J. Pharmacol. Exp. Ther.* **313**(1), 333-344 (2005).
2. Poręba, M., Stróżyk, A., Salvesen, G.S., *et al.* Caspase substrates and inhibitors. *Cold Spring Harb. Perspect. Biol.* **5**(8), A008680 (2013).
3. Yu, M.-S., Suen, K.-C., Kwok, N.-S., *et al.* Beta-amyloid peptides induces neuronal apoptosis via a mechanism independent of unfolded protein responses. *Apoptosis* **11**(5), 687-700 (2006).

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, 10/28/2022

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