

PRODUCT INFORMATION



Caspase-9 Inhibitor III (trifluoroacetate salt)

Item No. 24306

Formal Name: N-acetyl-L-leucyl-L- α -glutamyl-N-[(1S)-1-(carboxymethyl)-3-chloro-2-oxopropyl]-L-histidinamide, trifluoroacetate salt

Synonym: Ac-LEHD-CMK

MF: C₂₄H₃₅ClN₆O₉ • XCF₃COOH

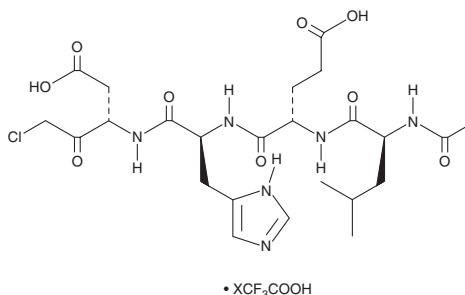
FW: 587.0

Purity: ≥95%

Supplied as: A 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

Caspase-9 inhibitor III (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the caspase-9 inhibitor III (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Caspase-9 inhibitor III is an inhibitor of caspase-9 with cardioprotective effects.¹ It reduces infarct size, lactate dehydrogenase (LDH) and creatine kinase (CK) release, and DNA smearing induced by ischemia and reperfusion in isolated rat hearts. Caspase-9 inhibitor III also inhibits Tnl proteolysis induced by matrix metalloproteinase-2 (MMP-2) when used at a concentration of 30 μ M.²

References

1. Sodhi, R.K., Singh, M., Singh, N., *et al.* Protective effects of caspase-9 and poly(ADP-ribose) polymerase inhibitors on ischemia-reperfusion-induced myocardial injury. *Arch. Pharm. Res.* **32(7)**, 1037-1043 (2009).
2. Castro, M.M., Fuah, J., Ali, M., *et al.* Inhibitory effects of caspase inhibitors on the activity of matrix metalloproteinase-2. *Biochem. Pharmacol.* **86(4)**, 469-475 (2013).

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, 12/01/2022

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