

PRODUCT INFORMATION



Ac-AAVALLPAVLLALLAP-DEVD-CHO (trifluoroacetate salt)

Item No. 27427

Formal Name:	N-acetyl-L-alanyl-L-alanyl-L-valyl-L-alanyl-L-leucyl-L-leucyl-L-prolyl-L-alanyl-L-valyl-L-leucyl-L-leucyl-L-alanyl-L-leucyl-L-leucyl-L-alanyl-L-prolyl-L- α -aspartyl-L- α -glutamyl-N-[(1S)-2-carboxy-1-formylethyl]-L-valinamide, trifluoroacetate salt	Ac-Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-Leu-Ala-Pro-Asp-Glu-Val-Asp-CHO
Synonyms:	Ac-Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-Leu-Ala-Pro-Asp-Glu-Val-Asp-CHO, Ac-AAVALLPAVLLALLAP-DEVD-aldehyde, Ac-AAVALLPAVLLALLAPDEVD-CHO, Caspase-3 Inhibitor I, DEVD-CHO-CPP 32	Leu-Ala-Leu-Leu-Ala-Pro-Asp-Glu-Val-Asp-CHO
MF:	$C_{94}H_{158}N_{20}O_{27} \cdot XCF_3COOH$	• XCF_3COOH
FW:	2,000.4	
Purity:	$\geq 95\%$	
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

Ac-AAVALLPAVLLALLAP-DEVD-CHO (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-AAVALLPAVLLALLAP-DEVD-CHO (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Ac-AAVALLPAVLLALLAP-DEVD-CHO (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Ac-AAVALLPAVLLALLAP-DEVD-CHO is a composite of Ac-DEVD-CHO (Item No. 10017), a peptide inhibitor of caspase-3 and -7, and a cell-permeable hydrophobic sequence derived from K-FGF.¹⁻³ It is an inhibitor of caspase-3 that reduces caspase-3 activity and apoptosis induced by prostaglandin E_2 (PGE_2 ; Item No. 14010) in rat cortical neurons when used at concentrations ranging from 1 to 2.5 μM .⁴

References

1. Garcia-Calvo, M., Peterson, E.P., Leiting, B., *et al.* Inhibition of human caspases by peptide-based and macromolecular inhibitors. *J. Biol. Chem.* **273**(49), 32608-32613 (1998).
2. Talanian, R.V., Quinlan, C., Trautz, S., *et al.* Substrate specificities of caspase family proteases. *J. Biol. Chem.* **272**(15), 9677-9682 (1997).
3. Lin, Y.Z., Tao, S., Veach, R.A., *et al.* Inhibition of nuclear translocation of transcription factor NF- κB by a synthetic peptide containing a cell membrane-permeable motif and nuclear localization sequence. *J. Biol. Chem.* **270**(24), 14255-14258 (1995).
4. Takadera, T., Yumoto, H., Tozuka, Y., *et al.* Prostaglandin E_2 induces caspase-dependent apoptosis in rat cortical cells. *Neurosci. Lett.* **317**(2), 61-64 (2002).

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, 11/30/2022

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