

# PRODUCT INFORMATION



## Ac-AAVALLPAVLLALLAP-IETD-CHO (trifluoroacetate salt)

Item No. 27438

<b>Formal Name:</b>	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-isoleucyl-L- $\alpha$ -glutamyl-N-[(1S)-2-carboxy-1-formylethyl]-L-threoninamide, trifluoroacetate salt	Ac—Ala—Ala—Val—Ala—Leu—Leu—Pro—Ala—Val—Leu—Leu—Ala—Leu—Leu—Ala—Pro—Ile—Glu—Thr—Asp—CHO
<b>Synonyms:</b>	Ac-Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-Leu-Ala-Pro-Ile-Glu-Thr-Asp-CHO, Caspase-8 Inhibitor I	• XCF <sub>3</sub> COOH
<b>MF:</b>	C <sub>95</sub> H <sub>162</sub> N <sub>20</sub> O <sub>26</sub> • XCF <sub>3</sub> COOH	
<b>FW:</b>	2,000.4	
<b>Purity:</b>	≥95%	
<b>Supplied as:</b>	A solid	
<b>Storage:</b>	-20°C	
<b>Stability:</b>	≥4 years	

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

### Laboratory Procedures

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

### Description

Ac-AAVALLPAVLLALLAP-IETD-CHO is a composite of Ac-IETD-CHO (Item No. 27100), a peptide inhibitor of caspase-8, and a cell-permeable hydrophobic sequence derived from K-FGF.<sup>1,2</sup> It is an inhibitor of both caspase-8 and granzyme B.<sup>1,3,4</sup> It prevents death of SKOV3 ovarian carcinoma cells transfected with *Survivin* and granzyme B-induced apoptosis (SAGA) when used at a concentration of 20  $\mu$ M.<sup>4</sup>

### 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. Lin, Y.Z., Tao, S., Veach, R.A., *et al.* Inhibition of nuclear translocation of transcription factor NF- $\kappa$ B by a synthetic peptide containing a cell membrane-permeable motif and nuclear localization sequence. *J. Biol. Chem.* **270**(24), 14255-14258 (1995).
3. Li, N. and Tan, W. Intelligent nanomedicine integrating diagnosis and therapy. *The Regents of the University of California. US 8,257,918 B2* (2012).
4. Caldas, H., Jaynes, F.O., Boyer, M.W., *et al.* *Survivin* and Granzyme B-induced apoptosis, a novel anticancer therapy. *Mol. Cancer. Ther.* **5**(3), 693-703 (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, 12/22/2022

#### 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