

# PRODUCT INFORMATION



## Boc-AEVD-CHO (trifluoroacetate salt)

Item No. 27148

**Formal Name:** N-[(1,1-dimethylethoxy)carbonyl]-L-alanyl-L- $\alpha$ -glutamyl-N-[(1S)-2-carboxy-1-formylethyl]-L-valinamide, trifluoroacetate salt

**Synonym:** Boc-Ala-Glu-Val-Asp-CHO

**MF:**  $C_{22}H_{36}N_4O_{10} \cdot XCF_3COOH$

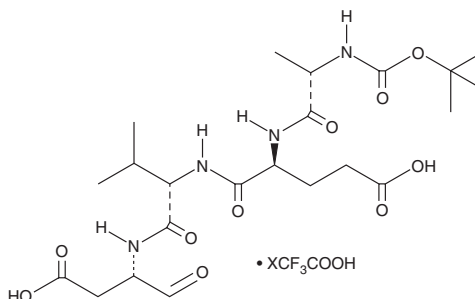
**FW:** 516.5

**Purity:**  $\geq 95\%$

**Supplied as:** A solid

**Storage:**  $-20^\circ C$

**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

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

### Description

Boc-AEVD-CHO is an inhibitor of caspase-8 ( $K_i = 1.6$  nM) that also inhibits caspase-1, -3, -6, and -9 ( $K_i$ s = <12, 42, 52, and 48 nM, respectively).<sup>1</sup> It is selective for these caspases over caspase-2, -4, -5, -7, and -10 ( $K_i$ s = >10,000, 375, 438, 425, and 320 nM, respectively). Boc-AEVD-CHO has been used to study the potential interference of caspase-8 activity in micronuclei detection.<sup>2</sup> It inhibits apoptosis induced by the aneugen nocodazole (Item No. 13857) in isolated human peripheral blood mononuclear cells (PBMCs) when used at a concentration of 300  $\mu M$ .

### 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. Decordier, I., Cundari, E., and Kirsch-Volders, M. Influence of caspase activity on micronuclei detection: A possible role for caspase-3 in micronucleation. *Mutagenesis* **20**(3), 173-179 (2005).

#### 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, 11/07/2022

#### 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