

# PRODUCT INFORMATION

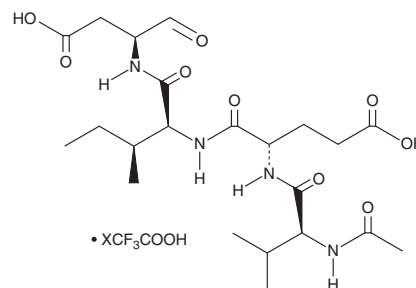


## Ac-VEID-CHO (trifluoroacetate salt)

Item No. 27437

**Formal Name:** (4S,7S,10S,13S)-10-((S)-sec-butyl)-7-(2-carboxyethyl)-13-formyl-4-isopropyl-2,5,8,11-tetraoxo-3,6,9,12-tetraazapentadecan-15-oic acid, trifluoroacetate salt

**Synonym:** Ac-Val-Glu-Ile-Asp-CHO  
**MF:** C<sub>22</sub>H<sub>36</sub>N<sub>4</sub>O<sub>9</sub> • XCF<sub>3</sub>COOH  
**FW:** 500.5  
**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

Ac-VEID-CHO (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Ac-VEID-CHO (trifluoroacetate salt) in water. The solubility of Ac-VEID-CHO (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Ac-VEID-CHO is an inhibitor of caspase-6 (IC<sub>50</sub> = 16.2 nM).<sup>1</sup> It is selective for caspase-6 over caspase-7 (IC<sub>50</sub> = 162.1 nM) but can also inhibit caspase-3 (IC<sub>50</sub> = 13.6 nM). Ac-VEID-CHO (100 μM) inhibits ricin-induced caspase-6-, but not caspase-3-, like activity in U937 cells.<sup>2</sup> It also inhibits caspase-6-like activity in Norway spruce cell extracts when used at a concentration of 2 μM, as well as in staurosporine-treated SKNAS neuroblastoma cells with an IC<sub>50</sub> value of 0.49 μM.<sup>1,3</sup>

### References

1. Mintzer, R., Ramaswamy, S., Shah, K., *et al.* A whole cell assay to measure caspase-6 activity by detecting cleavage of lamin A/C. *PLoS One* **7(1)**, e30376 (2012).
2. Komatsu, N., Oda, T., and Muramatsu, T. Involvement of both caspase-like proteases and serine proteases in apoptotic cell death induced by ricin, modeccin, diphtheria toxin, and *Pseudomonas* toxin. *J. Biochem.* **124(5)**, 1038-1044 (1998).
3. Bozhkov, P.V., Filonova, L.H., Suarez, M.F., *et al.* VEIDase is a principal caspase-like activity involved in plant programmed cell death and essential for embryonic pattern formation. *Cell Death Differ.* **11(2)**, 175-182 (2004).

#### 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, 12/08/2022

#### 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