

# PRODUCT INFORMATION



## Biotin-DEVD-CHO (trifluoroacetate salt)

Item No. 27133

**Formal Name:** N-[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]-L-α-aspartyl-L-α-glutamyl-N-[(1S)-2-carboxy-1-formylethyl]-L-valinamide, trifluoroacetate salt

**Synonym:** Biotin-Asp-Glu-Val-Asp-CHO

**MF:** C<sub>28</sub>H<sub>42</sub>N<sub>6</sub>O<sub>12</sub>S • XCF<sub>3</sub>COOH

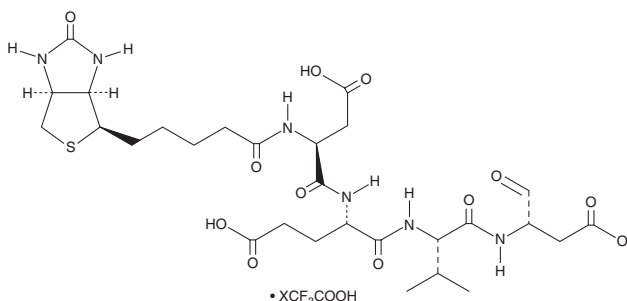
**FW:** 686.7

**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

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

### Description

Biotin-DEVD-CHO is a biotin-conjugated form of the caspase-3 and -7 inhibitor Ac-DEVD-CHO (Item No. 10017). It has been used for affinity purification of active caspase-3, -6, -7, and -8.<sup>1-3</sup> Biotin-DEVD-CHO has also been used for detection of active caspase-3 in tissue samples and permeabilized cells using affinity labeling with streptavidin conjugates.<sup>4,5</sup>

### References

1. Kumar, S. and Dorstyn, L. Analysing caspase activation and caspase activity in apoptotic cells. *Apoptosis: Methods and Protocols*. Erhardt, P. and Toth, A., eds., 2nd edition, *Humana Press* (2009).
2. Nicholson, D.W., Ali, A., Thornberry, N.A., et al. Identification and inhibition of the ICE/CED-3 protease necessary for mammalian apoptosis. *Nature* **376**, 37-43 (1995).
3. Sohn, D., Schulze-Osthoff, K., and Jänicke, R.U. Caspase-8 can be activated by interchain proteolysis without receptor-triggered dimerization during drug-induced apoptosis. *J. Biol. Chem.* **280**(7), 5267-5273 (2005).
4. Suria, H., Chau, L.A., Negrou, E., et al. Cytoskeletal disruption induces T cell apoptosis by a caspase-3 mediated mechanism. *Life Sci.* **65**(25), 2697-26707 (1999).
5. Huesmann, G.R. and Clayton, D.F. Dynamic role of postsynaptic caspase-3 and BIRC4 in zebra finch song-response habituation. *Neuron* **52**(6), 1061-1072 (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, 10/26/2022

#### 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