

PRODUCT INFORMATION

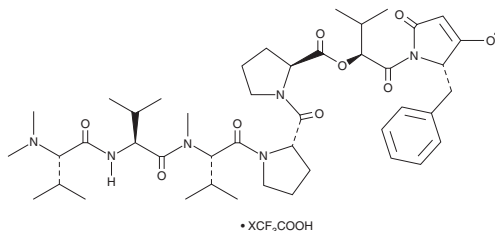


Dolastatin 15 (trifluoroacetate salt)

Item No. 30144

Formal Name: N,N-dimethyl-L-valyl-L-valyl-N-methyl-L-valyl-L-prolyl-L-proline, (1S)-1-[[[(2S)-2,5-dihydro-3-methoxy-5-oxo-2-(phenylmethyl)-1H-pyrrol-1-yl]carbonyl]-2-methylpropyl ester, trifluoroacetate salt

Synonym: NSC 617668
MF: C₄₅H₆₈N₆O₉ • XCF₃COOH
FW: 837.1
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Dolastatin 15 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the dolastatin 15 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Dolastatin 15 (trifluoroacetate salt) is soluble in the organic solvent methanol.

Description

Dolastatin 15 is a peptide that has been found in *D. auricularia* and has anticancer activity.¹⁻³ It induces cell cycle arrest at the G₂/M phase and inhibits proliferation of JVM-2 and EHEB human B cell leukemia cells when used at a concentration of 1 nM.¹ Dolastatin 15 also inhibits proliferation of RPMI-8826, U266, and IM-9 human multiple myeloma cells (IC₅₀s = 0.5-1 nM), as well as a panel of colon and ovarian cancer cell lines (IC₅₀s = 0.031-1.8 nM).^{2,3} It reduces tumor growth in a CH1 mouse xenograft model when administered at a dose of 5.25 mg/kg per day.³

References

1. Hu, Z.-B., Gignac, S.M., Quentmeier, H., *et al.* Effects of dolastatins on human b-lymphocytic leukemia cell lines. *Leuk. Res.* **17(4)**, 333-339 (1993).
2. Sato, M., Sagawa, M., Nakazato, T., *et al.* A natural peptide, dolastatin 15, induces G₂/M cell cycle arrest and apoptosis of human multiple myeloma cells. *Int. J. Oncol.* **30(6)**, 1453-1459 (2007).
3. Aherne, G.W., Hardcastle, A., Valenti, M., *et al.* Antitumour evaluation of dolastatins 10 and 15 and their measurement in plasma by radioimmunoassay. *Cancer Chemother. Pharmacol.* **38(3)**, 225-232 (1996).

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, 02/01/2023

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