

PRODUCT INFORMATION



Vanoxonin

Item No. 32931

CAS Registry No.: 86933-99-5

Formal Name: N-(2,3-dihydroxybenzoyl)-L-threonyl-N⁵-acetyl-N⁵-hydroxy-L-ornithine

Synonym: Antibiotic MG 245CF2A

MF: C₁₈H₂₅N₃O₉

FW: 427.4

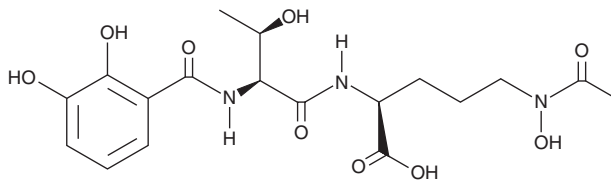
Purity: ≥80%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Saccharopolyspora hirsuta* MG245-CF2



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

Laboratory Procedures

Vanoxonin is supplied as a solid. A stock solution may be made by dissolving the vanoxonin in the solvent of choice, which should be purged with an inert gas. Vanoxonin is soluble in organic solvents such as methanol, DMSO, and acetic acid. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Vanoxonin is a bacterial metabolite that has been found in *S. hirsuta*.¹ It forms a complex with vanadium, and the complex inhibits thymidylate synthetase with an IC₅₀ value of 0.7 µg/ml. The vanoxonin-vanadium complex is cytotoxic to L1210 leukemia cells (MIC = 25 µg/ml) and increases survival in Ehrlich murine spontaneous adenocarcinoma and L1210 mouse tumor models.

Reference

1. Kanai, F., Sawa, T., Hamada, M., *et al.* Vanoxonin, a new inhibitor of thymidylate synthetase. *J. Antibiot. (Tokyo)* **36**(6), 656-660 (1983).

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, 11/10/2022

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