

PRODUCT INFORMATION

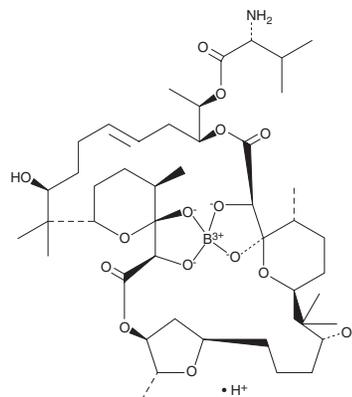


Boromycin

Item No. 28245

CAS Registry No.: 34524-20-4
Formal Name: (T-4)-[(1R)-1-[(1S,2R,5S,6R,8R,12R,14S,17R,18S,19R,22S,24Z,28S,30S,33R)-1,2,18,19-tetra(hydroxy-κO)-12,28-dihydroxy-6,13,13,17,29,29,33-heptamethyl-3,20-dioxo-4,7,21,34,35-pentaoxatetracyclo[28.3.1.1^{5,8}.1^{14,18}]hexatriacont-24-en-22-yl]ethyl D-valinato(4-)]-borate(1-), monohydrogen

Synonym: NSC 121380
MF: C₄₅H₇₃BNO₁₅ • H
FW: 879.9
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Boromycin is supplied as a solid. A stock solution may be made by dissolving the boromycin in the solvent of choice. Boromycin is soluble in organic solvents such as ethanol, methanol, and DMSO, which should be purged with an inert gas.

Description

Boromycin is a boron-containing macrolide antibiotic that has been found in *Streptomyces*.¹ Boromycin inhibits growth of *B. subtilis* (MIC = 0.05 μg/ml) and induces efflux of potassium ions from *B. subtilis* without affecting Na⁺/K⁺-ATPase activity.² It decreases the synthesis of protein, RNA, and DNA in *B. subtilis* when used at a concentration of 0.05 μg/ml. It inhibits the growth of *B. halodurans* (MIC = 10 ng/ml) and inhibits the futasoline pathway of menaquinone synthesis in *B. halodurans*.³ Boromycin (3.4 nM) reverses bleomycin-induced cell cycle arrest at the G₂ phase in Jurkat cells.⁴ It inhibits replication of the HIV-1 strains LAV-1 and RF and the HIV-2 strain LAV-2 in MT-4 cells (IC₅₀s = 0.008, 0.11, and 0.007 μM, respectively).¹ It also inhibits replication of a clinical isolate of HIV-1, strain KK-1, in MT-4 cells and peripheral blood mononuclear cells (PBMCs; IC₅₀s = 0.14 and <0.1 μM, respectively).

References

1. Kohno, J., Kawahata, T., Otake, T., *et al.* Boromycin, an anti-HIV antibiotic. *Biosci. Biotech. Biochem.* **60**(6), 1036-1037 (1996).
2. Pache, W. and Zähler, H. Metabolic products of microorganisms. 77. Studies on the mechanism of action of boromycin. *Arch. Mikrobiol.* **67**(2), 156-165 (1969).
3. Shimizu, Y., Ogasawara, Y., Matsumoto, A., *et al.* Aplasmomycin and boromycin are specific inhibitors of the futasoline pathway. *J. Antibiot. (Tokyo)* **71**(11), 968-970 (2018).
4. Arai, M., Koizumi, Y., Sato, H., *et al.* Boromycin abrogates bleomycin-induced G2 checkpoint. *J. Antibiot. (Tokyo)* **57**(10), 662-668 (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, 01/03/2023

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