

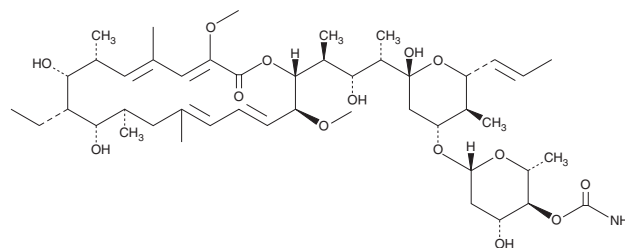
PRODUCT INFORMATION



Concanamycin A

Item No. 11050

CAS Registry No.: 80890-47-7
Formal Name: (3Z,5E,7R,8R,9S,10S,11R,13E,15E,17S,18R)-18-[[[(1S,2R,3S)-3-[(2R,4R,5S,6R)-4-[[4-O-(aminocarbonyl)-2,6-dideoxy-β-D-arabino-hexopyranosyl]oxy]tetrahydro-2-hydroxy-5-methyl-6-(1E)-1-propen-1-yl-2H-pyran-2-yl]-2-hydroxy-1-methylbutyl]-9-ethyl-8,10-dihydroxy-3,17-dimethoxy-5,7,11,13-tetramethyl-oxacyclooctadeca-3,5,13,15-tetraen-2-one



Synonyms: Antibiotic X 4357B, NSC 674620, X 4357B

MF: C₄₆H₇₅NO₁₄

FW: 866.1

Purity: ≥95%

Supplied as: A solution in acetonitrile

Storage: -20°C

Stability: ≥2 years

Item Origin: *Streptomyces* sp.

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

Laboratory Procedures

Concanamycin A is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as methanol and DMSO, purged with an inert gas, can be used.

Description

Concanamycin A is a plecomacrolide antibiotic produced by *Streptomyces* that blocks lysosomal acidification through selective inhibition of the vacuolar H⁺-ATPase (V-ATPase; EC_{50S} = 2.1-2.3 μM).¹⁻³ Concanamycin A blocks cell surface expression of viral glycoproteins without affecting their synthesis and, at 0.8 μM, prevents entry of influenza virus into cells.⁴

References

1. Sobota, J.A., Bäck, N., Eipper, B.A., *et al.* Inhibitors of the V₀ subunit of the vacuolar H⁺-ATPase prevent segregation of lysosomal- and secretory-pathway proteins. *J. Cell. Sci.* **122(19)**, 3542-3553 (2009).
2. Johnson, R.M., Allen, C., Melman, S.D., *et al.* Identification of inhibitors of V-ATPase pumps in yeast by HTS flow cytometry. *Anal. Biochem.* **398(2)**, 203-211 (2010).
3. Dröse, S., and Altendorf, K. Bafilomycins and concanamycins as inhibitors of V-ATPases and P-ATPases. *J. Exp. Biol.* **200(Pt 1)**, 1-8 (1997).
4. Guinea, R., and Carrasco, L. Requirement for vacuolar proton-ATPase activity during entry of influenza virus into cells. *J. Virol.* **69(4)**, 2306-2312 (1995).

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, 07/12/2019

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