

PRODUCT INFORMATION

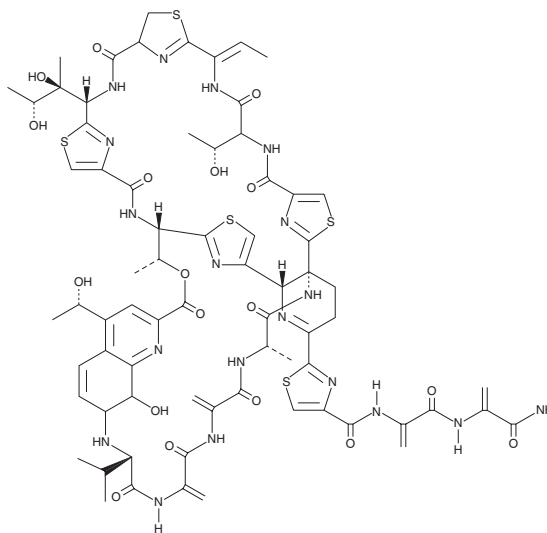


Siomycin A

Item No. 15378

CAS Registry No.: 12656-09-6
Formal Name: N-[(7R,8S)-2-carboxy-7,8-dihydro-8-hydroxy-4-[(1S)-1-hydroxyethyl]-7-quinoliny]-L-valyl-2,3-didehydroalanyl-2,3-didehydroalanyl-L-alanyl-2-[(4aR,11S,14Z,18S,21S,28S,32aS)-4a-amino-21-[(1S,2R)-1,2-dihydroxy-1-methylpropyl]-14-ethylidene-3,4,4a,9,10,11,12,13,14,18,19,20,21,27,28,32a-hexadecahydro-11,28-bis[(1R)-1-hydroxyethyl]-9,12,19,26-tetraoxo-17H,26H-8,5:18,15:25,22:32,29-tetranitrilo-5H,15H-pyrido[3,2-m][1,11,17,24,4,7,20,27]tetrathiatetraazacyclotriacontin-2-yl]-4-thiazolecarbonyl-2,3-didehydroalanyl-2,3-didehydroalaninamide, (1→5²⁸)-lactone

Synonym: NSC 285116
MF: C₇₁H₈₁N₁₉O₁₈S₅
FW: 1,648.9
Purity: ≥85%
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

Siomycin A is supplied as a solid. A stock solution may be made by dissolving the siomycin A in the solvent of choice, which should be purged with an inert gas. Siomycin A is soluble in DMSO and dimethyl formamide.

Description

The mammalian transcription factor forkhead box M1 (FoxM1) is induced during G₁ phase, with expression continuing through S phase and mitosis. Siomycin A is a peptide thiazole antibiotic, first isolated from *S. sioyaensis*, that interacts with the 23S rRNA of bacteria.¹ It inhibits FoxM1 (IC₅₀ = ~5 μM), preventing the expression of FoxM1-regulated genes, which includes FoxM1 itself.^{2,3} Through this mechanism, siomycin A prevents proliferation and induces apoptosis in certain types of stem cells as well as cancer cells.⁴⁻⁶

References

1. Tori, K., Tokura, K., Yoshimura, Y., et al. *J. Antibiot. (Tokyo)* **32(10)**, 1072-1077 (1979).
2. Radhakrishnan, S.K., Bhat, U.G., Hughes, D.E., et al. *Cancer Res.* **66(19)**, 9731-9735 (2006).
3. Gartel, A.L. *Front. Oncol.* **3**, 150 (2013).
4. Nakano, I., Joshi, K., Visnyei, K., et al. *Neuro. Oncol.* **13(6)**, 622-634 (2011).
5. Joshi, K., Banasavadi-Siddegowda, Y., Mo, X., et al. *Stem Cells* **31(6)**, 1051-1063 (2013).
6. Bhat, U.G., Halasi, M., and Gartel, A.L. *PLoS One* **4(5)**, e5592 (2009).

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, 11/14/2022

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