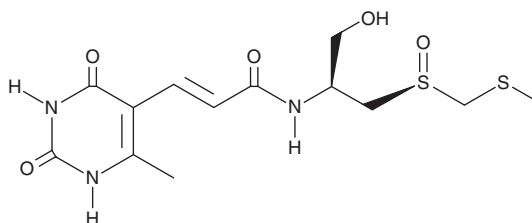


PRODUCT INFORMATION



Sparsomycin Item No. 29934

CAS Registry No.: 1404-64-4
Formal Name: (2E)-N-[(1S)-1-(hydroxymethyl)-2-[(R)-[(methylthio)methyl]sulfinyl]ethyl]-3-(1,2,3,4-tetrahydro-6-methyl-2,4-dioxo-5-pyrimidinyl)-2-propenamide
Synonyms: NSC 059729, NSC 59729, (+)-Sparsomycin, U-19183
MF: C₁₃H₁₉N₃O₅S₂
FW: 361.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

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

Description

Sparsomycin is a bacterial metabolite and a nucleoside analog of uracil that has been found in *S. sparsogenes* and has diverse biological activities.¹⁻³ It is active against KB carcinoma cells, Gram-positive and Gram-negative bacteria, and fungi.⁴ Sparsomycin is an inhibitor of peptidyl transferase that interferes with tRNA binding to the A-site of the peptidyl transfer center and increases the binding of peptidyl-tRNA to the P-site.² It inhibits protein synthesis in bacteria, archaea, and eukaryotes.^{2,3} Sparsomycin reduces tumor growth in a P388 mouse leukemia model and in a Walker 256 carcinosarcoma rat model.⁵

References

- Ottenheijm, H.C., van den Broek, L.A., Ballesta, J.P., *et al.* Chemical and biological aspects of sparsomycin, an antibiotic from *Streptomyces*. *Prog. Med. Chem.* **23**, 219-268 (1986).
- Wilson, D.N. The A-Z of bacterial translation inhibitors. *Crit. Rev. Biochem. Mol. Biol.* **44(6)**, 393-433 (2009).
- Lazaro, E., Van den Broek, L.A., San Felix, A., *et al.* Chemical, biochemical and genetic endeavours characterizing the interaction of sparsomycin with the ribosome. *Biochimie* **73(7-8)**, 1137-1143 (1991).
- Owen, S.P., Dietz, A., and Camiener, G.W. Sparsomycin, a new antitumor antibiotic. I. Discovery and biological properties. *Antimicrob. Agents Chemother.* **1962**, 772-779 (1963).
- Zylicz, Z., Wagener, D.J., van Rennes, H., *et al.* *In vivo* antitumor activity of sparsomycin and its analogues in eight murine tumor models. *Invest. New Drugs* **6(4)**, 285-292 (1988).

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, 03/05/2020

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