

PRODUCT INFORMATION



Streptimidone

Item No. 29984

CAS Registry No.: 738-72-7
Formal Name: 4-[(2R,5S,6E)-2-hydroxy-5,7-dimethyl-4-oxo-6,8-nonadien-1-yl]-2,6-piperidinedione

Synonyms: Ao58A, NSC 66645

MF: C₁₆H₂₃NO₄

FW: 293.4

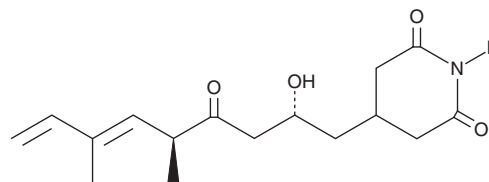
Purity: ≥95%

Supplied as: A residue

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

Streptimidone is supplied as a residue. A stock solution may be made by dissolving the streptimidone in the solvent of choice, which should be purged with an inert gas. Streptimidone is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Streptimidone is a bacterial metabolite that has been found in *Streptomyces* and has diverse biological activities.¹⁻³ It inhibits protein synthesis in a cell-free assay when used at a concentration of 50 µg/ml.² It reduces tumor growth in H.S. No. 1 human sarcoma and H.Ep. No. 3 human epidermoid carcinoma xenograft models when administered at doses of 12.5 and 25 mg/kg per day.³

References

1. Kondo, H., Oritani, T., and Kiyota, H. Synthesis and antifungal activity of the four stereoisomers of streptimidone, a glutarimide antibiotic from *Streptomyces rimosus* forma *paromomycinus*. *Eur. J. Org. Chem.* (20), 3459-3462 (2000).
2. Bennett, L.L., Ward, V.L., and Brockman, R.W. Inhibition of protein synthesis *in vitro* by cycloheximide and related glutarimide antibiotics. *Biochim. Biophys. Acta.* **103**(3), 478-485 (1965).
3. Teller, M.N. Antibiotics in experimental cancer chemotherapy. *Trans. N. Y. Acad. Sc.* **24**, 158-166 (1961).

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, 08/25/2023

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