

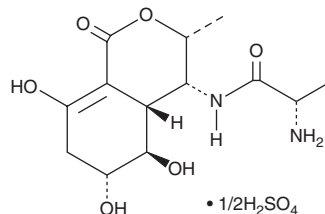
PRODUCT INFORMATION



Actinobolin (hemisulfate)

Item No. 42359

CAS Registry No.: 18802-17-0
Formal Name: 2S-amino-N-[(3R,4R,4aR,5R,6R)-3,4,4a,5,6,7-hexahydro-5,6,8-trihydroxy-3-methyl-1-oxo-1H-2-benzopyran-4-yl]-propanamide, hemisulfate
MF: $C_{13}H_{20}N_2O_6 \cdot 1/2H_2SO_4$
FW: 349.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
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

Actinobolin (hemisulfate) is supplied as a solid. A stock solution may be made by dissolving the actinobolin (hemisulfate) in the solvent of choice, which should be purged with an inert gas. Actinobolin (hemisulfate) is soluble in methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Actinobolin is a microbial metabolite that has been found in *Pseudomonas* and has antibiotic and anticancer activities.¹ It is active against *E. coli* (MIC = 170 μM). Actinobolin decreases protein synthesis in *E. coli* (IC₅₀ = 30 μM). It inhibits the proliferation of adenocarcinoma 755 cells (IC₅₀ = 70 μM).²

References

1. Hori, M., Suzukake, K., Ishikawa, C., *et al.* Biochemical studies on bactobolin in relation to actinobolin. *J. Antibiot. (Tokyo)* **34**(4), 465-468 (1981).
2. Smithers, D., Bennett, L.L., Jr., and Struck, R.F. Inhibition of protein synthesis in mammalian cells by actinobolin. *Mol. Pharmacol.* **5**(5), 433-445 (1969).

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, 07/0/2025

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