

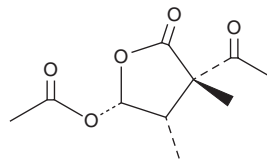
PRODUCT INFORMATION



Acetomycin

Item No. 31410

CAS Registry No.: 510-18-9
Formal Name: (3S,4S,5R)-3-acetyl-5-(acetyloxy) dihydro-3,4-dimethyl-2(3H)-furanone
Synonyms: (-)-Acetomycin, NSC 350598
MF: C₁₀H₁₄O₅
FW: 214.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces ramulosus*



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

Laboratory Procedures

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

Description

Acetomycin is a γ -lactone microbial metabolite originally isolated from *S. ramulosus* that has anticancer activity.¹ It inhibits proliferation of HCT-8 human colon and L1210 mouse leukemia cancer cells (IC₅₀s = 1.5 and 2.2 μ g/ml, respectively).

Reference

1. Mamber, S.W., Mitulski, J.D., Hamelchle, K.L., *et al.* Biological effects of acetomycin. I. Activity against tumor cells *in vitro* and *in vivo*. *J. Antibiot.* (Tokyo) **49(1)**, 73-76 (1987).

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, 12/08/2022

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