

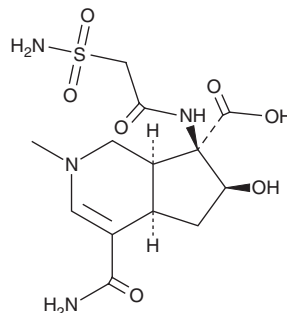
PRODUCT INFORMATION



Altemicidin

Item No. 32897

CAS Registry No.: 125399-82-8
Formal Name: 4-(aminocarbonyl)-7-[[2-(aminosulfonyl)acetyl]amino]-2,4aR,5,6S,7R,7aS-hexahydro-6-hydroxy-2-methyl-1H-cyclopenta[c]pyridine-7-carboxylic acid
Synonym: (-)-Altemicidin
MF: C₁₃H₂₀N₄O₇S
FW: 376.4
Purity: ≥98%
UV/Vis.: λ_{max}: 294 nm
Supplied as: A powder
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

Altemicidin is supplied as a powder. A stock solution may be made by dissolving the altemicidin in the solvent of choice, which should be purged with an inert gas. Altemicidin is soluble in ethanol.

Description

Altemicidin is a monoterpene alkaloid originally isolated from *S. siyoaensis* with acaricidal and anticancer activities.¹ It is acaricidal to two-spotted spider mites (*T. urticae*) in a greenhouse pot test at concentrations of 10 and 100 ppm. Altemicidin inhibits the growth of murine L1210 lymphocytic leukemia and IMC carcinoma cells (IC₅₀s = 0.84 and 0.82 μg/ml, respectively). It is toxic to mice with an LD₅₀ value of 0.3 mg/kg.

Reference

1. Takahashi, A., Kurasawa, S., Ikeda, D., *et al.* Altemicidin, a new acaricidal and antitumor substance. I. Taxonomy, fermentation, isolation and physico-chemical and biological properties. *J. Antibiot. (Tokyo)* **42(11)**, 1556-1561 (1989).

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, 12/02/2022

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