

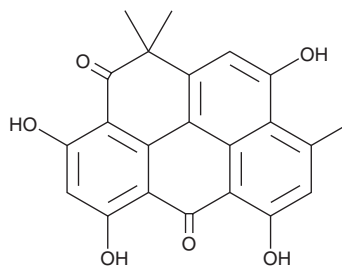
PRODUCT INFORMATION



Resistomycin

Item No. 20685

CAS Registry No.: 20004-62-0
Formal Name: 3,5,7,10-tetrahydroxy-1,1,9-trimethyl-2H-benzo[cf]pyrene-2,6(1H)-dione
Synonyms: Antibiotic A-3733A, Heliomycin
MF: C₂₂H₁₆O₆
FW: 376.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Resistomycin is a natural antibiotic produced by several actinobacteriae, including *Streptomyces*. Resistomycin, as well as culture broths rich in resistomycin, have antioxidant activity, modulate apoptosis, and can be cytotoxic against cancer cell lines.^{1,2} Resistomycin also has activity against HIV-1 protease.³

References

1. Shiono, Y., Shiono, N., Seo, S., *et al.* Effects of polyphenolic anthrone derivatives, resistomycin and hypercin, on apoptosis in human megakaryoblastic leukemia CMK-7 cell line. *Z. Naturforsch. C.* **57(9-10)**, 923-929 (2002).
2. Vijayabharathi, R., Bruheim, P., Andreassen, T., *et al.* Assessment of resistomycin, as an anticancer compound isolated and characterized from *Streptomyces aurantiacus* AAA5. *J. Microbiol.* **49(6)**, 920-926 (2011).
3. Roggo, B. E., Petersen, F., Delmendo, R., *et al.* 3-Alkanoyl-5-hydroxymethyl tetronic acid homologues and resistomycin: New inhibitors of HIV-1 protease. I. Fermentation, isolation and biological activity. *J. Antibiot. (Tokyo)* **47(2)**, 136-142 (1994).

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, 07/01/2025

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