

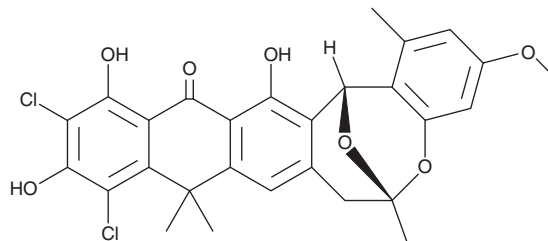
PRODUCT INFORMATION



Bischloroanthrabenzoxocinone

Item No. 19439

CAS Registry No.: 866022-28-8
Formal Name: (6R,16R)-10,12-dichloro-6,7,9,16-tetrahydro-11,13,15-trihydroxy-3-methoxy-1,6,9,9-tetramethyl-6,16-epoxy-14H-anthra[2,3-d][1]benzoxocin-14-one
Synonyms: BABX, (-)-Bischloroanthrabenzoxocinone
MF: C₂₈H₂₄Cl₂O₇
FW: 543.4
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

Bischloroanthrabenzoxocinone (BABX) is supplied as a residue. A stock solution may be made by dissolving the BABX in the solvent of choice, which should be purged with an inert gas. BABX is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Bacterial type II fatty acid synthesis (FAS-II) is mediated by a series of enzymes, each of which may be targeted by potential antibiotics.¹ BABX is an inhibitor of FAS-II, blocking fatty acid synthesis in *S. aureus* and *E. coli* with IC₅₀ values of 11.4 and 35.3 µg/ml, respectively.² It inhibits the growth of *S. aureus* and permeable *E. coli* strains with minimum inhibitory concentrations ranging from 0.2-0.4 µg/ml.^{2,3} BABX also displays binding to liver X receptors (LXRs), inhibiting agonist binding in an LXRA-scintillation proximity assay (IC₅₀ = 10 µM).³

References

1. Yao, J. and Rock, C.O. How bacterial pathogens eat host lipids: Implications for the development of fatty acid synthesis therapeutics. *J. Biol. Chem.* **290**(10), 5940-5946 (2015).
2. Kodali, S., Galgoci, A., Young, K., et al. Determination of selectivity and efficacy of fatty acid synthesis inhibitors. *J. Biol. Chem.* **280**(2), 1669-1677 (2005).
3. Herath, K.B., Jayasuriya, H., Guan, Z., et al. Anthrabenzoxocinones from *Streptomyces* sp. as liver X receptor ligands and antibacterial agents. *J. Nat. Prod.* **68**(9), 1437-1440 (2005).

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, 11/04/2022

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