

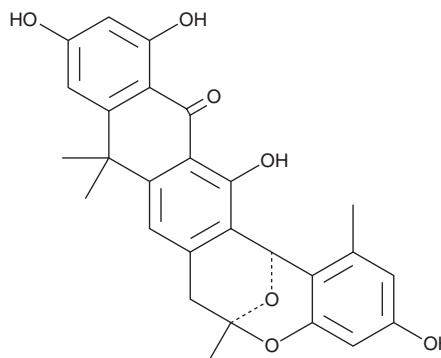
PRODUCT INFORMATION



BE-24566B

Item No. 29955

CAS Registry No.: 149466-04-6
Formal Name: (6S,16S)-6,7,9,16-tetrahydro-3,11,13,15-tetrahydroxy-1,6,9,9-tetramethyl-6,16-epoxy-14H-anthra[2,3-d][1]benzoxocin-14-one
Synonym: L-755,805
MF: C₂₇H₂₄O₇
FW: 460.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

BE-24566B is supplied as a solid. A stock solution may be made by dissolving the BE-24566B in the solvent of choice, which should be purged with an inert gas. BE-24566B is soluble in methanol, ethanol, dichloromethane, and DMSO.

Description

BE-24566B is a polyketide fungal metabolite originally isolated from *S. violaceusniger*.¹ It is active against *B. subtilis*, *B. cereus*, *S. aureus*, *M. luteus*, *E. faecalis*, and *S. thermophilus* (MICs = 1.56, 1.56, 1.56, 1.56, 3.13, and 3.13 µg/ml, respectively). BE-24566B is an endothelin (ET) receptor antagonist (IC₅₀s = 11 and 3.9 µM for ET_A and ET_B receptors, respectively).²

References

1. Kojiri, K., Nakajima, S., Fuse, A., *et al.* BE-24566B, a new antibiotic produced by *Streptomyces violaceusniger*. *J. Antibiot. (Tokyo)* **48(12)**, 1506-1508 (1995).
2. Lam, Y.K.T., Hensens, O., Helms, G., *et al.* L-755,805, a new polyketide endothelin binding inhibitor from an actinomycete. *Tetrahedron Lett.* **36(12)**, 2013-2016 (1995).

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, 05/18/2020

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