

PRODUCT INFORMATION

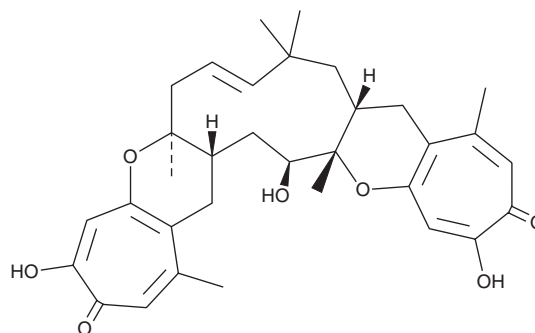


Eupenifeldin

Item No. 25114

CAS Registry No.: 151803-45-1
Formal Name: (6aS,8E,11aR,18aS,19S,20aS)-6a,7,10,11,11a,18a,19,20,20a,21-decahydro-4,16,19-trihydroxy-1,6a,10,10,13,18a-hexamethyl-3H-dicyclohepta[e,e']cycloundeca[1,2-b:5,6-b']dipyran-3,15(12H)-dione

MF: C₃₃H₄₀O₇
FW: 548.7
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

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

Description

Eupenifeldin is a pentacyclic bispolone fungal metabolite originally isolated from *E. brefeldianum* ATCC 74184.¹ It is cytotoxic against HCT116 and HCTVM46 colon carcinoma cells *in vitro* (IC₅₀s = 0.005 and 0.002 µg/ml, respectively). *In vivo*, eupenifeldin (0.3-2 mg/kg per day) increases median survival time in a mouse model of P388 leukemia.

Reference

1. Mayerl, F., Gao, Q., Huang, S., *et al.* Eupenifeldin, a novel cytotoxic bispolone from *Eupenicillium brefeldianum*. *J. Antibiot. (Tokyo)* **46(7)**, 1082-1088 (1993).

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, 12/22/2022

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