

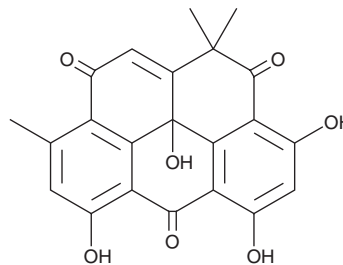
PRODUCT INFORMATION



Resistoflavine

Item No. 24178

CAS Registry No.: 29706-96-5
Formal Name: (-)-3,5,7,11b-tetrahydroxy-1,1,9-trimethyl-2H-benzo[cd]pyrene-2,6,10(1H,11bH)-trione
MF: C₂₂H₁₆O₇
FW: 392.4
Purity: ≥98%
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

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

Description

Resistoflavine is a metabolite of the marine actinomycete *S. chibaensis*.¹ It slows the growth of and is cytotoxic to HMO2 and HepG2 cells with mean lethal concentrations (LC₅₀s) of 0.013 and 0.016 µg/ml, respectively.

Reference

1. Gorajana, A., Venkatesan, M., Vinjamuri, S., et al. Resistoflavine, cytotoxic compound from a marine actinomycete, *Streptomyces chibaensis* AUBN₁/7. *Microb. Res.* **162**(4), 322-327 (2007).

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, 11/29/2022

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