

PRODUCT INFORMATION



Xanthoquinodin A1

Item No. 28792

CAS Registry No.: 151063-27-3
Formal Name: (4R,4aS,7aS,15R)-rel-2,3,4,7,8,13,15,17-octahydro-1,4,12,14,16-pentahydroxy-10-methyl-8,13,17-trioxo-4aH-7a,15-ethenonaphtho[2',3':4,5]cyclohepta[1,2-b]xanthene-4a-carboxylic acid, methyl ester

Synonym: Xanthoquinidone C

MF: C₃₁H₂₄O₁₁

FW: 572.5

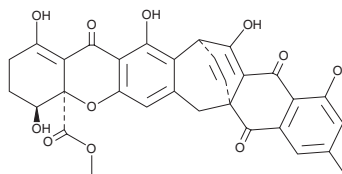
Purity: ≥70%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Chaetomium* sp.



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

Laboratory Procedures

Xanthoquinodin A1 is supplied as a solid. A stock solution may be made by dissolving the xanthoquinodin A1 in the solvent of choice, which should be purged with an inert gas. Xanthoquinodin A1 is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

Xanthoquinodin A1 is a fungal metabolite that has been found in *Humicola* and has diverse biological activities.^{1,2} It inhibits *E. tenella* schizont formation in BHK-21 cells with a minimum effective concentration (MEC) value of 0.02 µg/ml.¹ Xanthoquinodin A1 is active against *B. subtilis*, *M. luteus*, *S. aureus*, *A. laidlawii*, and *B. fragilis* in a disc assay when used at a concentration of 1 mg/ml. It is also active against *B. cereus* (MIC = 0.44 µM).² Xanthoquinodin A1 is cytotoxic to KB, MCF-7, and NCI H187 cancer cells.

References

1. Tabata, N., Suzumura, Y., Tomoda, H., et al. Xanthoquinodins, new anticoccidial agents produced by *Humicola* sp. Production, isolation and physico-chemical and biological properties. *J. Antibiot. (Tokyo)* **46**(5), 749-755 (1993).
2. Tantapakul, C., Promgool, T., Kanokmedhakul, K., et al. Bioactive xanthoquinodins and epipolythiodioxopiperazines from *Chaetomium globosum* 7s-1, an endophytic fungus isolated from *Rhapis cochinchinensis* (Lour.) Mart. *Nat. Prod. Res.* **34**(4), 494-502 (2020).

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, 08/16/2023

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