

PRODUCT INFORMATION



Galbinic Acid

Item No. 29584

CAS Registry No.: 56691-88-4
Formal Name: 5-[(acetyloxy)methyl]-1,3-dihydro-1,4,10-trihydroxy-8-methyl-3,7-dioxo-7H-isobenzofuro[4,5-b][1,4]benzodioxepin-11-carboxaldehyde

MF: C₂₀H₁₄O₁₁

FW: 430.3

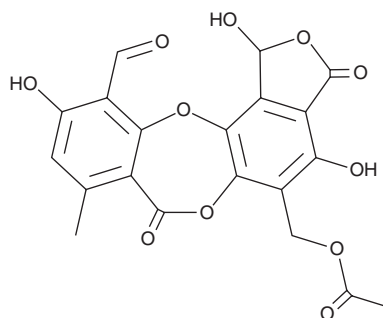
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Usnea undulata*



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

Laboratory Procedures

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

Description

Galbinic acid is a depsidone lichen metabolite that has been found in *U. undulata*.^{1,2} It is active against the Gram-positive bacteria *B. cereus*, *B. subtilis*, and *S. aureus*, but not *S. epidermidis* (MICs = 62.5, 62.5, 250, and >250 µg/ml, respectively), as well as the Gram-negative bacterium *E. coli*, but not *S. sonnei* (MICs = 125 and >250 µg/ml, respectively).³

References

1. Salgado, F., Albornoz, L., Cortéz, C., *et al.* Secondary metabolite profiling of species of the genus *Usnea* by UHPLC-ESI-OT-MS-MS. *Molecules* **23**(1), E54 (2017).
2. Elix, J.A. and Engkaninan, U. The structure of galbinic acid. A depsidone from the lichen *Usnea undulata*. *Aust. J. Chem.* **28**(8), 1793-1797 (1975).
3. Sultana, N. and Afolayan, A.J. A new depsidone and antibacterial activities of compounds from *Usnea undulata* Stirton. *J. Asian Nat. Prod. Res.* **13**(12), 1158-1164 (2011).

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, 08/15/2023

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