

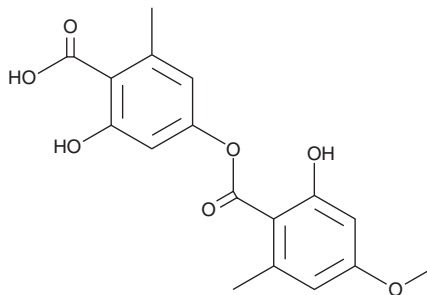
PRODUCT INFORMATION



Evernic Acid

Item No. 17430

CAS Registry No.: 537-09-7
Formal Name: 2-hydroxy-4-[(2-hydroxy-4-methoxy-6-methylbenzoyl)oxy]-6-methylbenzoic acid
Synonym: NSC 81164
MF: C₁₇H₁₆O₇
FW: 332.3
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 269, 306 nm
Supplied as: A crystalline 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

Evernic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the evernic acid in the solvent of choice, which should be purged with an inert gas. Evernic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of evernic acid in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Evernic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, evernic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Evernic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Evernic acid is a secondary metabolite produced by some species of lichen that can bind to allosteric sites on the protein surface of FAS-II enzymes and produce antibacterial and antiplasmodial effects.¹ It inhibits two key plasmodial FAS-II enzymes *PfFabZ* and *PfFabI* (IC₅₀ = 10.7 and 36.1 μM, respectively) but shows low efficacy against the malaria parasite *P. berghei* (IC₅₀ = 77.3 μM).¹

Reference

1. Lauinger, I.L., Vivas, L., Perozzo, R., *et al.* Potential of lichen secondary metabolites against *Plasmodium* liver stage parasites with FAS-II as the potential target. *J. Nat. Prod.* **76**(6), 1064-1070 (2013).

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, 10/13/2022

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