

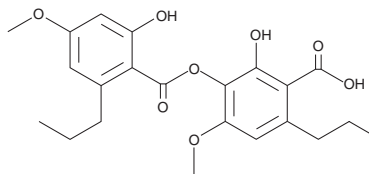
PRODUCT INFORMATION



Sekikaic Acid

Item No. 13993

CAS Registry No.: 607-11-4
Formal Name: 2-hydroxy-3-[(2-hydroxy-4-methoxy-6-propylbenzoyl)oxy]-4-methoxy-6-propylbenzoic acid
MF: C₂₂H₂₆O₈
FW: 418.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Ramalina* sp.



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

Laboratory Procedures

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

Description

Sekikaic acid is a phenolic lichen metabolite that has been found in *H. obscurata* and has diverse biological activities.¹⁻³ It scavenges superoxide and DPPH (Item No. 14805) radicals in cell-free assays when used at a concentration of 0.5 μM.¹ Sekikaic acid is active against the rg recombinant strain of respiratory syncytial virus (RSV; IC₅₀ = 5.69 μg/ml).² It inhibits the protein-protein interaction between mixed lineage leukemia 1 (MLL1) and the GACKIX domain of CREB-binding protein (CBP; IC₅₀ = 34 μM).³

References

1. Thadhani, V.M., Choudhary, M.I., Ali, S., *et al.* Antioxidant activity of some lichen metabolites. *Nat. Prod. Res.* **25(19)**, 1827-1837 (2011).
2. Lai, D., Odimegwu, D.C., Esimone, C., *et al.* Phenolic compounds with *in vitro* activity against respiratory syncytial virus from the Nigerian lichen *Ramalina farinacea*. *Planta Med.* **79(15)**, 1440-1446 (2013).
3. Majmudar, C.Y., Højfeldt, J.W., Arevang, C.J., *et al.* Sekikaic acid and lobaric acid target a dynamic interface of the coactivator CBP/p300. *Angew. Chem. Int. Ed. Engl.* **51(45)**, 11258-11262 (2012).

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, 12/13/2022

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