

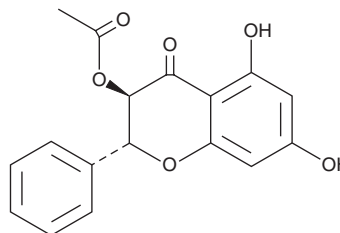
PRODUCT INFORMATION



3-O-Acetyl Pinobanksin

Item No. 43409

CAS Registry No.: 52117-69-8
Formal Name: 3-(acetyloxy)-2R,3R-dihydro-5,7-dihydroxy-2-phenyl-4H-1-benzopyran-4-one
Synonym: Pinobanksin 3-O-Acetate
MF: C₁₇H₁₄O₆
FW: 314.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant



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

Laboratory Procedures

3-O-Acetyl pinobanksin is supplied as a solid. A stock solution may be made by dissolving the 3-O-acetyl pinobanksin in the solvent of choice, which should be purged with an inert gas. 3-O-Acetyl pinobanksin is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

3-O-Acetyl pinobanksin is a flavonoid that has been found in propolis, a resin produced by honeybees, and has antioxidant and antiparasitic activities.^{1,2} It reduces the formation of advanced glycation end products (AGEs) in a cell-free assay (IC₅₀ = 0.06 mM).¹ 3-O-Acetyl pinobanksin is active against chloroquine-resistant *P. falciparum* (IC₅₀ = 15.62 μM).²

References

1. Boisard, S., Le Ray, A.M., Gatto, J., *et al.* Chemical composition, antioxidant and anti-AGEs activities of a French poplar type propolis. *J. Agric. Food Chem.* **62**(6), 1344-1351 (2014).
2. Ranaivoarisoa, R.H., Rasoarivelo Ralambonirina, S.T., Randriamialinoro, F., *et al.* Antiplasmodial activity of the extracts and flavonoids Isolated from *Helichrysum gymnocephalum* humbert (Asteraceae) from Madagascar. *African natural plant products, Volume III: Discoveries and innovations in chemistry, bioactivity, and applications.* American Chemical Society (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, 06/24/2025

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