

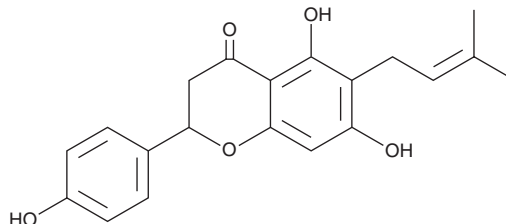
PRODUCT INFORMATION



(±)-6-Prenylnaringenin

Item No. 34499

CAS Registry No.: 68682-01-9
Formal Name: 2,3-dihydro-5,7-dihydroxy-2-(4-hydroxyphenyl)-6-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one
Synonym: (±)-6-PN
MF: C₂₀H₂₀O₅
FW: 340.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Synthetic



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

Laboratory Procedures

(±)-6-Prenylnaringenin ((±)-6-PN) is supplied as a solid. A stock solution may be made by dissolving the (±)-6-PN in the solvent of choice, which should be purged with an inert gas. (±)-6-PN is soluble in methanol.

Description

(±)-6-PN is a flavonoid that has been found in the hop plant, *H. lupulus*, and has diverse biological activities.^{1,2} It inhibits the T-type voltage-gated calcium (Ca_v) channel Ca_v3.2 and high-voltage-activated calcium currents in whole-cell patch-clamp assays using HEK293 cells (IC₅₀s = 0.912 and 4.765 μM, respectively).¹ It inhibits proliferation of PC3 and DU145 prostate cancer cells (IC₅₀s = 18.4 and 29.1 μM, respectively).² (±)-6-PN (20 and 30 mg/kg) reduces mechanical allodynia in a mouse model of neuropathic pain induced by partial sciatic nerve ligation.¹

References

1. Sekiguchi, F., Fujita, T., Deguchi, T., *et al.* Blockade of T-type calcium channels by 6-prenylnaringenin, a hop component, alleviates neuropathic and visceral pain in mice. *Neuropharmacology* **138**, 232-244 (2018).
2. Delmulle, L., Bellahcène, A., Dhooze, W., *et al.* Anti-proliferative properties of prenylated flavonoids from hops (*Humulus lupulus* L.) in human prostate cancer cell lines. *Phytomedicine* **13(9-10)**, 732-734 (2006).

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, 10/14/2021

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