

PRODUCT INFORMATION



(2S)-6-Prenylnaringenin

Item No. 35138

CAS Registry No.: 68236-13-5
Formal Name: (2S)-2,3-dihydro-5,7-dihydroxy-2-(4-hydroxyphenyl)-6-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one

Synonym: (2S)-6-PN

MF: C₂₀H₂₀O₅

FW: 340.4

Purity: ≥98%

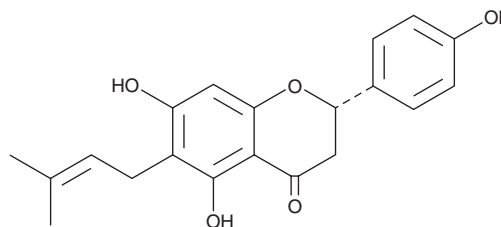
UV/Vis.: λ_{max}: 215, 295 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Humulus lupulus*



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

Laboratory Procedures

(2S)-6-Prenylnaringenin is supplied as a solid. A stock solution may be made by dissolving the (2S)-6-prenylnaringenin in the solvent of choice, which should be purged with an inert gas. (2S)-6-Prenylnaringenin is soluble in DMSO.

Description

(2S)-6-Prenylnaringenin is a prenylated 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 in HEK293 cells and high-voltage-activated calcium currents in differentiated NG 108-15 cells in whole-cell patch-clamp assays (IC₅₀s = 0.928 and 2.005 μM, respectively).¹ (2S)-6-Prenylnaringenin reduces the activity of aldose reductase 2 (ALR2; IC₅₀ = 6.2 μM for the recombinant human enzyme).² It reduces mechanical allodynia in a mouse model of sodium hydrosulfide-induced somatic pain when administered at a dose of 10 pmol/paw.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. Shim, S.H., Kim, Y., Lee, J.Y., *et al.* Aldose reductase inhibitory activity of the compounds from the seed of *Psoralea corylifolia*. *J. Korean Soc. Appl. Biol. Chem.* **52(5)**, 568-572 (2009).

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, 11/07/2022

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