

PRODUCT INFORMATION



(-)-8-Prenylnaringenin

Item No. 37906

CAS Registry No.: 53846-50-7

Formal Name: (2S)-2,3-dihydro-5,7-dihydroxy-2-(4-hydroxyphenyl)-8-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one

Synonyms: Flavaprenin, (-)-8-PN, (S)-8-Prenylnaringenin, Sophoraflavanone B

MF: $C_{20}H_{20}O_5$

FW: 340.4

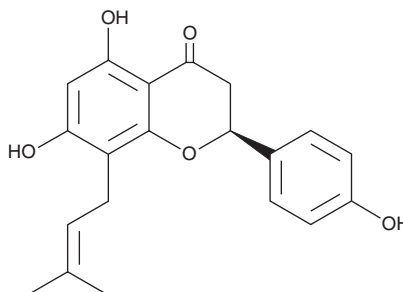
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Synthetic



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

Laboratory Procedures

(-)-8-Prenylnaringenin is supplied as a solid. A stock solution may be made by dissolving the (-)-8-prenylnaringenin in the solvent of choice, which should be purged with an inert gas. (-)-8-Prenylnaringenin is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

(-)-8-Prenylnaringenin is a prenylated flavonoid and phytoestrogen that has been found in *H. lupulus* and has estrogenic activity.¹ It binds to recombinant human estrogen receptor α (ER α) and ER β in cell-free assays and induces ER activation in a reporter assay using yeast expressing human ER in a concentration-dependent manner. (-)-8-Prenylnaringenin increases uterine vascular permeability in ovariectomized mice in a dose-dependent manner.

Reference

1. Milligan, S., Kalita, J., Pocock, V., *et al.* Oestrogenic activity of the hop phyto-oestrogen, 8-prenylnaringenin. *Reproduction* **123**(2), 235-242 (2002).

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, 01/16/2025

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