

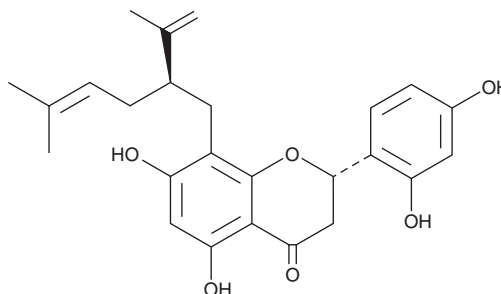
PRODUCT INFORMATION



Sophoraflavanone G

Item No. 34755

CAS Registry No.: 97938-30-2
Formal Name: (2S)-2-(2,4-dihydroxyphenyl)-2,3-dihydro-5,7-dihydroxy-8-[(2R)-5-methyl-2-(1-methylethenyl)-4-hexen-1-yl]-4H-1-benzopyran-4-one
Synonyms: Kushenol F, Norkurarinone, Vexibinol
MF: C₂₅H₂₆O₆
FW: 424.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Sophora flavescens*



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

Laboratory Procedures

Sophoraflavanone G is supplied as a solid. A stock solution may be made by dissolving the sophoraflavanone G in the solvent of choice, which should be purged with an inert gas. Sophoraflavanone G is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of sophoraflavanone G in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Sophoraflavanone G is a flavonoid that has been found in *S. flavescens* and has diverse biological activities.¹⁻⁴ It is active against 21 strains of methicillin-resistant *S. aureus* (MRSA; MICs = 3.13-6.25 µg/ml) and inhibits the activities of α-glucosidase and β-amylase (K_is = 5.6 and 30.6 µM, respectively).^{1,2} Sophoraflavanone G is cytotoxic to HL-60 and HepG2 cancer cells (IC₅₀s = 12.5 and 13.3 µM, respectively) and induces apoptosis in HL-60 cells when used at concentrations of 10 or 25 µM.³ It reduces LPS-induced production of prostaglandin E₂ (PGE₂; Item No. 14010) in RAW 264.7 cells when used at concentrations ranging from 1 to 50 µM.⁴ Sophoraflavanone G (250 mg/kg) inhibits carrageenan-induced paw edema in rats.

References

1. Sato, M., Tsuchiya, H., Takase, I., *et al.* Antibacterial activity of flavanone isolated from *Sophora exigua* against methicillin-resistant *Staphylococcus aureus* and its combination with antibiotics. *Phytother. Res.* **9(7)**, 509-512 (1995).
2. Kim, J.H., Ryu, Y.B., Kang, N.S., *et al.* Glycosidase inhibitory flavonoids from *Sophora flavescens*. *Biol. Pharm. Bull.* **29(2)**, 302-305 (2006).
3. Ko, W.G., Kang, T.H., Kim, N.Y., *et al.* Lavandulylfavonoids: A new class of *in vitro* apoptogenic agents from *Sophora flavescens*. *Toxicol. In Vitro* **14(5)**, 429-433 (2000).
4. Kim, D.W., Chi, Y.S., Son, K.H., *et al.* Effects of sophoraflavanone G, a prenylated flavonoid from *Sophora flavescens*, on cyclooxygenase-2 and *in vivo* inflammatory response. *Arch. Pharm. Res.* **25(3)**, 329-335 (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, 12/20/2022

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