

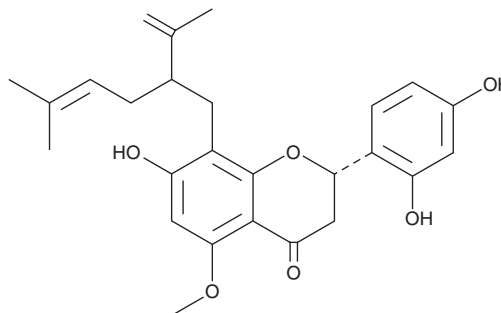
PRODUCT INFORMATION



Kurarinone

Item No. 38966

CAS Registry No.: 1049662-21-6
Formal Name: (2S)-2-(2,4-dihydroxyphenyl)-2,3-dihydro-7-hydroxy-5-methoxy-8-[5-methyl-2-(1-methylethenyl)-4-hexen-1-yl]-4H-1-benzopyran-4-one
MF: C₂₆H₃₀O₆
FW: 438.5
Purity: ≥98%
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

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

Description

Kurarinone is a lavandulylated flavanone that has been found in *S. flavescens* and has diverse biological activities.¹⁻⁴ It inhibits COX-1 and 5-lipoxygenase (5-LO; IC₅₀s = 0.6-1 and 22 μM, respectively).¹ Kurarinone inhibits LPS-induced nitric oxide (NO) production in RAW 264.7 cells (IC₅₀ = 26.6 μM).² It reduces potassium chloride- or norepinephrine-induced contractions in rat thoracic aortic helical strips when used at concentrations of 10 and 100 μM.³ *In vivo*, kurarinone (20 and 40 mg/kg) decreases disease severity and retinal neutrophil infiltration in a mouse model of experimental autoimmune uveitis induced by IRBP (1-20) (Item No. 36727).⁴

References

1. Chi, Y.S., Jong, H.G., Son, K.H., *et al.* Effects of naturally occurring prenylated flavonoids on enzymes metabolizing arachidonic acid: Cyclooxygenases and lipoxygenases. *Biochem. Pharmacol.* **62**(9), 1185-1191 (2001).
2. Kwon, J., Basnet, S., Lee, J.W., *et al.* Chemical constituents isolated from the Mongolian medicinal plant *Sophora alopecuroides* L. and their inhibitory effects on LPS-induced nitric oxide production in RAW 264.7 macrophages. *Bioorg. Med. Chem. Lett.* **25**(16), 3314-3318 (2015).
3. Yamahara, J., Kobayashi, G., Iwamoto, M., *et al.* Vasodilatory active principles of *Sophora flavescens* root. *J. Ethnopharmacol.* **29**(1), 79-85 (1990).
4. Gu, C., Liu, Y., Lv, J., *et al.* Kurarinone regulates Th17/Treg balance and ameliorates autoimmune uveitis via Rac1 inhibition. *J. Adv. Res.* **69**, 381-398 (2025).

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, 06/23/2025

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