

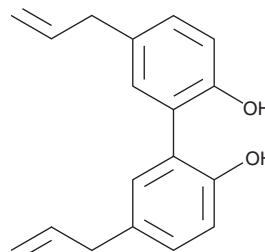
PRODUCT INFORMATION



Magnolol

Item No. 14233

CAS Registry No.: 528-43-8
Formal Name: 5,5'-di-2-propen-1-yl-[1,1'-biphenyl]-2,2'-diol
Synonym: NSC 293099
MF: C₁₈H₁₈O₂
FW: 266.3
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 292 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Magnolol is supplied as a crystalline solid. A stock solution may be made by dissolving the magnolol in the solvent of choice, which should be purged with an inert gas. Magnolol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of magnolol in DMSO is approximately 16 mg/ml and approximately 20 mg/ml in ethanol and DMF.

Magnolol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, magnolol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Magnolol has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Magnolol is a bioactive compound isolated from the bark of *M. officinalis* that has been used in Asian traditional medicine for the treatment of anxiety, sleep disorders, and allergic diseases. Magnolol can activate cannabinoid (CB) receptors, behaving as a partial agonist with selectivity for the peripheral CB₂ subtype (EC₅₀ = 3.28 μM; K_i = 1.44 μM) versus central CB₁ (EC₅₀ = 18.3 μM; K_i = 3.15 μM).¹

Reference

1. Rempel, V., Fuchs, A., Hinz, S., *et al.* Magnolia extract, magnolol, and metabolites: Activation of cannabinoid CB₂ receptors and blockade of the related GPR55. *ACS Med. Chem. Lett.* **4**, 41-45 (2013).

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, 11/23/2022

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