

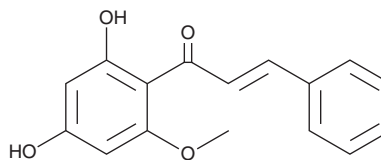
PRODUCT INFORMATION



Cardamonin

Item No. 18310

CAS Registry No.: 19309-14-9
Formal Name: (2E)-1-(2,4-dihydroxy-6-methoxyphenyl)-3-phenyl-2-propen-1-one
Synonyms: Alpinetin chalcone, Cardamomin
MF: C₁₆H₁₄O₄
FW: 270.3
Purity: ≥98%
UV/Vis.: λ_{max}: 343 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

Cardamonin is supplied as a crystalline solid. A stock solution may be made by dissolving the cardamonin in the solvent of choice, which should be purged with an inert gas. Cardamonin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cardamonin in ethanol is approximately 1 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Cardamonin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cardamonin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cardamonin has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cardamonin, isolated from the fruits of *Alpinia* species, is a chalconoid with anti-inflammatory and anti-tumor activity.^{1,2} It has been shown to suppress nitric oxide and prostaglandin E₂ synthesis, to suppress cyclooxygenase-2 expression, and to inhibit NF-κB signaling.^{1,2} It can also target the Bcl-2 protein, inducing apoptosis in cancer cells.¹

References

1. Yadav, V.R., Prasad, S., Sung, B., *et al.* The role of chalcones in suppression of NF-κB-mediated inflammation and cancer. *Int. Immunopharmacol.* **11(3)**, 295-309 (2011).
2. Sung, B., Prasad, S., Yadav, V.R., *et al.* Cancer cell signaling pathways targeted by spice-derived nutraceuticals. *Nutr. Cancer* **64(2)**, 173-197 (2012).

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, 12/13/2022

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