

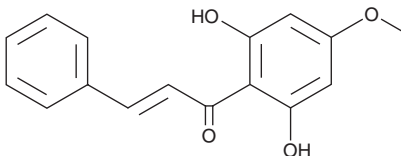
PRODUCT INFORMATION



Pinostrobin Chalcone

Item No. 43023

CAS Registry No.: 18956-15-5
Formal Name: (2E)-1-(2,6-dihydroxy-4-methoxyphenyl)-3-phenyl-2-propen-1-one
Synonyms: 2',6'-Dihydroxy 4'-methoxychalcone, DMC
MF: C₁₆H₁₄O₄
FW: 270.3
Purity: ≥98%
Supplied as: A 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

Pinostrobin chalcone is supplied as a solid. A stock solution may be made by dissolving the pinostrobin chalcone in the solvent of choice, which should be purged with an inert gas. Pinostrobin chalcone is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Pinostrobin chalcone is a chalcone that has been found in *H. cymosum* and has diverse biological activities.¹⁻⁴ It reduces ferric to ferrous iron in a ferric reducing antioxidant power (FRAP) assay (IC₅₀ = 334 μM).¹ Pinostrobin chalcone is an inhibitor of α-glucosidase (IC₅₀ = 12.94 μM). It decreases the proliferation of *L. amazonensis* promastigotes when used at a concentration of 1 μg/ml.² Pinostrobin chalcone is cytotoxic to KB squamous cell carcinoma, MCF-7 breast, Ca Ski cervical, and HCT116 colorectal cancer cells (IC₅₀s = 6.2, 7.3, 7.7, and 54.1 μM, respectively).³ It inhibits LPS-induced increases in TNF-α and IL-1β in J774A.1 macrophages (IC₅₀s = 69 and 43 μM, respectively).⁴

References

1. Jadalla, B.M.I.S., Moser, J.J., Sharma, R., *et al.* In vitro alpha-glucosidase and alpha-amylase inhibitory activities and antioxidant capacity of *Helichrysum cymosum* and *Helichrysum pandurifolium* schrank constituents. *Separations* **9**(8), 190 (2022).
2. Torres-Santos, E.C., Moreira, D.L., Kaplan, M.A., *et al.* Selective effect of 2',6'-dihydroxy-4'-methoxychalcone isolated from *Piper aduncum* on *Leishmania amazonensis*. *Antimicrob Agents Chemother.* **43**(5), 1234-1241 (1999).
3. Malek, S.N., Phang, C.W., Ibrahim, H., *et al.* Phytochemical and cytotoxic investigations of *Alpinia mutica* rhizomes. *Molecules* **16**(1), 583-589 (2011).
4. Patel, N.K. and Bhutani, K.K. Pinostrobin and Cajanus lactone isolated from *Cajanus cajan* (L.) leaves inhibits TNF-α and IL-1 production: *In vitro* and *in vivo* experimentation. *Phytomedicine* **21**(7), 946-953 (2014).

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, 02/10/2025

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