

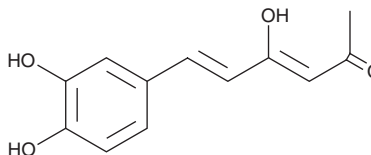
PRODUCT INFORMATION



Hispolon

Item No. 38415

CAS Registry No.: 173933-40-9
Formal Name: 6-(3,4-dihydroxyphenyl)-4-hydroxy-3Z,5E-hexadien-2-one
MF: $C_{12}H_{12}O_4$
FW: 220.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Hispolon is supplied as a solid. A stock solution may be made by dissolving the hispolon in the solvent of choice, which should be purged with an inert gas. Hispolon is soluble in methanol and acetone.

Description

Hispolon is a phenol originally isolated from *I. hispidus* that has diverse biological activities.¹⁻³ It scavenges superoxide and hydroxyl radicals in cell-free assays, as well as reduces LPS-induced production of reactive oxygen species (ROS) in RAW 264.7 cells when used at concentrations of 5 and 10 μM .¹ Hispolon (60-80 μM) inhibits cell proliferation and induces apoptosis in C6 rat glioma and DBTRG-05MG human glioblastoma cells.² Hispolon (5 and 10 mg/kg) reduces tumor volume in a DBTRG-05MG mouse xenograft model. It also reduces acetic acid-induced writhing in mice when administered at a dose of 20 mg/kg.³

References

1. Lee, E.K., Koh, E.M., Kim, Y.N., *et al.* Immunomodulatory effect of hispolon on LPS-induced RAW264.7 cells and mitogen/alloanitgen-stimulated spleen lymphocytes of mice. *Pharmaceutics* **14**(7), 1423 (2022).
2. Liao, K.-F., Chiu, T.-L., Chang, S.-F., *et al.* Hispolon induces apoptosis, suppresses migration and invasion of glioblastoma cells and inhibits gbm xenograft tumor growth in vivo. *Molecules* **26**(15), 4497 (2021).
3. Chang, H.-Y., Sheu, M.-J., Yang, C.-H., *et al.* Analgesic effects and the mechanisms of anti-inflammation of hispolon in mice. *Evid. Based Complement. Alternat. Med.* **2011**, 478246 (2009).

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, 04/03/2024

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