

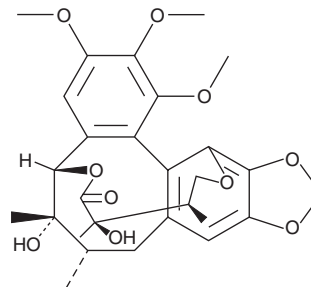
PRODUCT INFORMATION



Gomisin D

Item No. 30697

CAS Registry No.: 60546-10-3
Formal Name: (5S,6S,7S,13aR,16R,17R)-5,6,7,8-tetrahydro-6,17-dihydroxy-1,2,3-trimethoxy-6,7,16,17-tetramethyl-5,13-(epoxybutanoxy)benzo[3,4]cycloocta[1,2-f][1,3]benzodioxol-18-one
MF: C₂₈H₃₄O₁₀
FW: 530.6
Purity: ≥98%
UV/Vis.: λ_{max}: 217 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Schisandra chinensis*



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

Laboratory Procedures

Gomisin D is supplied as a crystalline solid. A stock solution may be made by dissolving the gomisin D in the solvent of choice, which should be purged with an inert gas. Gomisin D is soluble in DMSO at a concentration of approximately 10 mM.

Description

Gomisin D is a lignan that has been found in *S. chinensis* and has diverse biological activities, including enzyme inhibitory and antioxidant properties.¹⁻³ It inhibits acetylcholinesterase (AChE) with an IC₅₀ value of 7.84 μM.¹ Gomisin D scavenges ABTS (Item No. 27317) and 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals (IC₅₀s = 24.5 and 203 μM, respectively).³ It inhibits the UDP-glucuronosyltransferase (UGT) isoforms UGT1A1 and UGT1A3 by 89.9 and 83.1%, respectively, when used at a concentration of 100 μM.² It is selective for UGT1A1 and UGT1A3 over other UGT isoforms at 100 μM.

References

1. Hung, T.M., Na, M., Min, B.S., *et al.* Acetylcholinesterase inhibitory effect of lignans isolated from *Schisandra chinensis*. *Arch. Pharm. Res.* **30**(6), 685-690 (2007).
2. Song, J.-H., Cui, L., An, L.-B., *et al.* Inhibition of UDP-glucuronosyltransferases (UGTs) activity by constituents of *Schisandra chinensis*. *Phytother. Res.* **29**(10), 1658-1664 (2015).
3. Poornima, B., Kumar, D.A., Siva, B., *et al.* Advanced glycation end-products inhibitors isolated from *Schisandra grandiflora*. *Nat. Prod. Res.* **30**(4), 493-496 (2016).

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, 11/03/2022

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