

PRODUCT INFORMATION

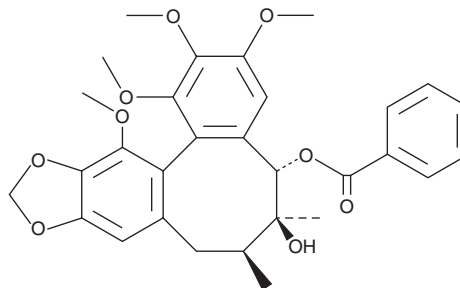


Gomisin C

Item No. 18643

CAS Registry No.: 58546-56-8
Formal Name: (5S,6S,7S,13aS)-5,6,7,8-tetrahydro-1,2,3,13-tetramethoxy-6,7-dimethylbenzo[3,4]cycloocta[1,2-f][1,3]benzodioxole-5,6-diol 5-benzoate

Synonym: Schisantherin A
MF: C₃₀H₃₂O₉
FW: 536.6
Purity: ≥98%
UV/Vis.: λ_{max}: 222 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

Gomisin C is supplied as a crystalline solid. A stock solution may be made by dissolving the gomisin C in the solvent of choice, which should be purged with an inert gas. Gomisin C is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of gomisin C in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Gomisin C is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, gomisin C should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Gomisin C has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Gomisin C, a lignan first isolated from the fruits of *S. chinensis*, scavenges oxygen radicals and reduces the production of reactive oxygen species by inhibiting neutrophil activity.¹ It has been shown to attenuate respiratory burst in rat peripheral neutrophils via suppression of NADPH oxidase (39% inhibition at 30 µg/ml) and inhibition of cytosolic calcium release from intracellular stores (36% inhibition at 50 µg/ml).² Gomisin C can induce liver microsomal cytochrome P450, thus facilitating the clearance of xenobiotics and protecting against hepatotoxicity in rodent models of liver injury.³

References

1. Lin, T.-J., Liu, G.-T., Li, X.-J., et al. Detection of free radical scavenging activity of schisanhenol by electron spin resonance. *Acta. Pharmacol. Sin.* **11(6)**, 534-539 (1990).
2. Wang, J.P., Raung, S.L., Hsu, M.f., et al. Inhibition by gomisin C (a lignan from *Schizandra chinensis*) of the respiratory burst of rat neutrophils. *Br. J. Pharmacol.* **113(3)**, 945-953 (1994).
3. Li, X.-Y. Bioactivity of neolignans from *fructus Schizandrae*. *Mem. Inst. Oswaldo Cruz* **86 Suppl 2**, 31-37 (1991).

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, 11/11/2022

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