

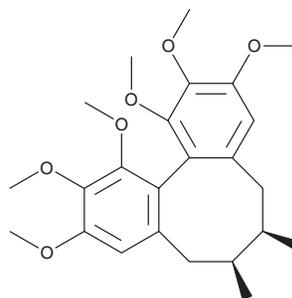
PRODUCT INFORMATION



Schisandrin A

Item No. 19849

CAS Registry No.: 61281-38-7
Formal Name: (6R,7S,12aR)-5,6,7,8-tetrahydro-1,2,3,10,11,12-hexamethoxy-6,7-dimethyl-dibenzo[a,c]cyclooctene
Synonyms: Deoxyschizandrin, Wuweizisu A
MF: C₂₄H₃₂O₆
FW: 416.5
Purity: ≥98%
UV/Vis.: λ_{max}: 217 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

Schisandrin A is supplied as a crystalline solid. A stock solution may be made by dissolving the schisandrin A in the solvent of choice, which should be purged with an inert gas. Schisandrin A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of schisandrin A in ethanol and DMSO is approximately 20 mg/ml and approximately 25 mg/ml in DMF.

Schisandrin A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, schisandrin A should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Schisandrin A has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Schisandrin A is a lignan that has been found in *Schisandra* and has diverse biological activities.¹⁻⁴ It binds to adiponectin receptor 2 (IC₅₀ = 3.2 μM).¹ Schisandrin A (10, 20, and 50 μM) reduces LPS-induced production of nitric oxide (NO), IL-6, and TNF-α and apoptosis in primary mouse microglial cells.² It reverses P-glycoprotein-mediated doxorubicin resistance in MCF-7/dox cells when used at a concentration of 20 μM.³ Schisandrin A (4, 12, and 36 mg/kg) reverses short-term and spatial memory deficits, as well as decreases in cerebral superoxide dismutase (SOD) and glutathione peroxidase (GPX) activities induced by amyloid β (1-42) (Aβ₄₂; Item No. 20574) in a mouse model of Alzheimer's disease.⁴

References

1. Sun, Y., Zang, Z., Zhong, L., *et al.* Identification of adiponectin receptor agonist utilizing a fluorescence polarization based high throughput assay. *PLoS One* **8(5)**, e63354 (2013).
2. Song, F., Zeng, K., Liao, L., *et al.* Schisandrin A inhibits microglia-mediated neuroninflammation through inhibiting TRAF6-NF-κB and Jak2-stat3 signaling pathways. *PLoS One* **11(2)**, (2016).
3. Zhang, Z.-L., Jiang, Q.-C., and Wang, S.-R. Schisandrin A reverses doxorubicin-resistant human breast cancer cell line by the inhibition of P65 and Stat3 phosphorylation. *Breast Cancer* **25(2)**, 233-242 (2018).
4. Hu, D., Li, C., Han, N., *et al.* Deoxyschizandrin isolated from the fruits of *Schisandra chinensis* ameliorates Aβ₁₋₄₂-induced memory impairment in mice. *Planta Med.* **78(12)**, 1332-1336 (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, 12/19/2022

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