

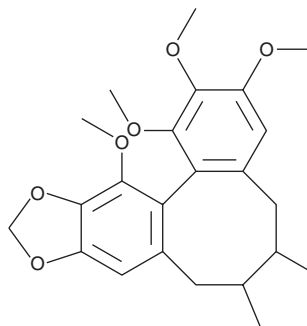
PRODUCT INFORMATION



Schizandrin B

Item No. 24968

CAS Registry No.: 61281-37-6
Formal Name: (6R,7S,13aR)-rel-5,6,7,8-tetrahydro-1,2,3,13-tetramethoxy-6,7-dimethyl-benzo[3,4]cycloocta[1,2-f][1,3]benzodioxole
Synonym: Wuweizisu B
MF: C₂₃H₂₈O₆
FW: 400.5
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

Schizandrin B is supplied as a crystalline solid. A stock solution may be made by dissolving the schizandrin B in the solvent of choice, which should be purged with an inert gas. Schizandrin B is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of schizandrin B in these solvents is approximately 2, 5, and 30 mg/ml, respectively.

Schizandrin B is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, schizandrin B should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Schizandrin B has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Schizandrin B is a dibenzocyclooctadiene that has been found in *F. schisandrae* and *S. chinensis* with diverse biological activities.¹⁻³ Schizandrin B (50 μM) completely inhibits anti-CD3- and anti-CD28-induced release of the cytokines IL-2, IL-3, IL-4, IL-6, and IFN-γ in T cells.⁴ In A549 adenocarcinoma cells, schizandrin B (1-30 μM) dose-dependently reduces cell survival and UV-induced phosphorylation of p53 and checkpoint kinase 1 (Chk1).¹ Schizandrin B reduces kinase activity of the serine/threonine-protein kinase ATR with an IC₅₀ value of 7.25 μM for the recombinant enzyme. *In vivo*, schizandrin B (25-100 mg/kg, p.o.) dose-dependently inhibits lipid peroxidation, formation of DNA strand breaks, and NADPH oxidase-dependent oxygen production induced by doxorubicin (Item No. 15007) in mouse heart.² It also reduces the expression of p53, 3-nitrotyrosine, phosphorylated p38 MAPK and MAPKAPK-2, and matrix metalloproteinase-2 (MMP-2) and MMP-9. Schizandrin B (1 or 2 mmol/kg per day, p.o.) pretreatment reduces *tert*-butylhydroperoxide-induced mortality and malondialdehyde formation and increases glutathione (GSH) levels and Se-glutathione peroxidase (GSH-Px) activity in brain in mice.³

References

1. Nishida, H., Tatewaki, N., Nakajima, Y., et al. *Nucleic Acids Res.* **37(17)**, 5678-5689 (2009).
2. Thandavarayan, R.A., Giridharan, V.V., Arumugam, S., et al. *PLoS One* **10(3)**, e0119214 (2015).
3. Ko, K.M. and Lam, B.Y. *Mol. Cell. Biochem.* **238(1-2)**, 181-186 (2002).
4. Checker, R., Patwardhan, R.S., Sharma, D., et al. *Free Radic. Biol. Med.* **53(7)**, 1421-1430 (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, 12/05/2022

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