

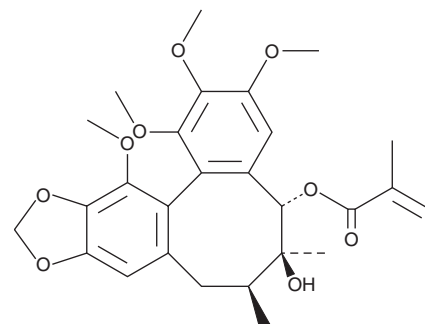
# PRODUCT INFORMATION



## Schisantherin B

Item No. 26716

**CAS Registry No.:** 58546-55-7  
**Formal Name:** (2Z)-2-methyl-2-butenoic acid, (5S,6S,7S,13aS)-5,6,7,8-tetrahydro-6-hydroxy-1,2,3,13-tetramethoxy-6,7-dimethylbenzo[3,4]cycloocta[1,2-f][1,3]benzodioxol-5-yl ester  
**Synonyms:** Gomisin B, Schisantherin B, Wuweizi ester B  
**MF:** C<sub>28</sub>H<sub>34</sub>O<sub>9</sub>  
**FW:** 514.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219 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

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

Schisantherin B is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, schisantherin B should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Schisantherin 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

Schisantherin B (STB) is a lignan originally isolated from *S. sphenanthera* fruit.<sup>1</sup> It inhibits formation of advanced glycation end products (AGEs) *in vitro*.<sup>2</sup> STB (0.1 mg/kg) decreases escape latency in the Morris water maze and increases spontaneous alternation in the Y-maze in a mouse model of Alzheimer's disease induced by amyloid-β (1-42) (Item No. 20574).<sup>3</sup> It increases glial expression of glutamate transporter 1 (GLP-1) and decreases immobility time in the forced swim test (FST) in mice, indicating antidepressant-like activity.<sup>4</sup> STB also increases spontaneous alternation in the Y-maze in a mouse model of stress-induced short-term learning impairment induced by the FST.

### References

1. Liu, C.S., Fang, S.D., Huang, M.F., *et al.* Studies on the active principles of *Schisandra sphenanthera* Rehd. *et Wils.* The structures of schisantherin A, B, C, D, E, and the related compounds. *Sci. Sin.* **21(4)**, 483-502 (1978).
2. 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).
3. Xu, M., Dong, Y., Wan, S., *et al.* Schisantherin B ameliorates Aβ<sub>1-42</sub>-induced cognitive decline via restoration of GLT-1 in a mouse model of Alzheimer's disease. *Physiol. Behav.* **167**, 265-273 (2016).
4. Xu, M., Xiao, F., Wang, M., *et al.* Schisantherin B improves the pathological manifestations of mice caused by behavior desperation in different ages-depression with cognitive impairment. *Biomol. Ther. (Seoul)* **27(2)**, 160-167 (2019).

#### 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, 12/21/2022

#### 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