

PRODUCT INFORMATION



Ciwujianoside B

Item No. 42502

CAS Registry No.: 114902-16-8

Formal Name: (3β)-3-[[2-O-(6-deoxy-α-L-mannopyranosyl)-α-L-arabinopyranosyl]oxy]-30-noroleana-12,20(29)-dien-28-oic acid, O-6-deoxy-α-L-mannopyranosyl-(1→4)-O-β-D-glucopyranosyl-(1→6)-β-D-glucopyranosyl ester
Yemuoside YM₁₀

Synonym:

MF: C₅₈H₉₂O₂₅

FW: 1,189.3

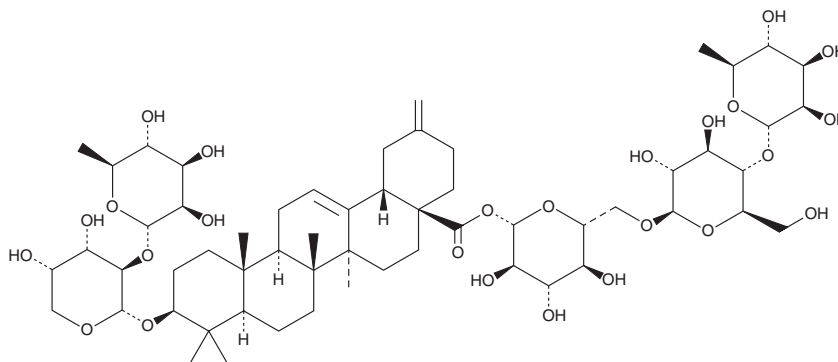
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Eleutherococcus senticosus*



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

Laboratory Procedures

Ciwujianoside B is supplied as a solid. A stock solution may be made by dissolving the ciwujianoside B in the solvent of choice, which should be purged with an inert gas. Ciwujianoside B is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Ciwujianoside B is sparingly soluble (1-10 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Ciwujianoside B is a triterpenoid saponin that has been found in *A. senticosus* and has diverse biological activities.¹⁻³ It inhibits α-glucosidase and scavenges peroxynitrite radicals in cell-free assays (IC₅₀s = 0.74 and 34.8 μM, respectively).¹ Ciwujianoside B (40 mg/kg) increases survival and reverses radiation-induced decreases in the number of peripheral blood leukocytes in irradiated mice.² It improves memory in the novel object recognition test in mice when administered at a dose of 0.5 mg/kg.³

References

1. Kwon, R.H., Na, H., Kim, J.H., *et al.* Comprehensive profiling of phenolic compounds and triterpenoid saponins from *Acanthopanax senticosus* and their antioxidant, α-glucosidase inhibitory activities. *Sci. Rep.* **14**(1), 26330 (2024).
2. Li, Y.-R., Cao, W., Guo, J., *et al.* Comparative investigations on the protective effects of rhodioside, ciwujianoside-B and astragaloside IV on radiation injuries of the hematopoietic system in mice. *Phytother. Res.* **25**(5), 644-653 (2011).
3. Yamauchi, Y., Ge, Y.-W., Yoshimatsu, K., *et al.* Memory enhancement by oral administration of extract of *Eleutherococcus senticosus* leaves and active compounds transferred in the brain. *Nutrients* **11**(5), 1142 (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, 01/14/2025

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