

PRODUCT INFORMATION



Dipsacoside B

Item No. 35624

CAS Registry No.: 33289-85-9
Formal Name: (3 β ,4 α)-3-[[2-O-(6-deoxy- α -L-mannopyranosyl)- α -L-arabinopyranosyl]oxy]-23-hydroxy-olean-12-en-28-oic acid, 6-O- β -D-glucopyranosyl- β -D-glucopyranosyl

Synonyms: Glycoside L-G₃, Tauroside G₃

MF: C₅₃H₈₆O₂₂

FW: 1,075.3

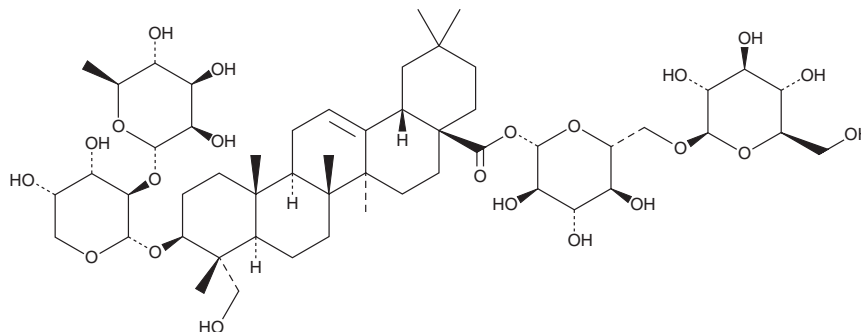
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Lonicera confusa* DC



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

Laboratory Procedures

Dipsacoside B is supplied as a solid. A stock solution may be made by dissolving the dipsacoside B in the solvent of choice, which should be purged with an inert gas. Dipsacoside B is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of dipsacoside B in these solvents is approximately 5, 16, and 12 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of dipsacoside B can be prepared by directly dissolving the solid in aqueous buffers. The solubility of dipsacoside B in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Dipsacoside B is a triterpenoid saponin that has been found in *Lonicera* and has hepatoprotective activity.^{1,2} It enhances acetaminophen-induced autophagy and inhibits acetaminophen-induced apoptosis in HepG2 cells when used at a concentration of 0.5 μ M.² Dipsacoside B (50 mg/kg) increases survival in a mouse model of acute liver failure induced by acetaminophen (Item No. 10024).

References

1. Yang, R., Min, K., Wang, Y., *et al.* Rapid semi-quantitative analysis of hemolytic triterpenoid saponins in *Lonicerae Flos* crude drugs and preparations by paper spray mass spectrometry. *Talanta* **239**, 123148 (2022).
2. Chen, S., Li, M., Jiang, W., *et al.* The role of Neu1 in the protective effect of dipsacoside B on acetaminophen-induced liver injury. *Ann. Transl. Med.* **8(13)**, 823 (2020).

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, 12/12/2022

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