

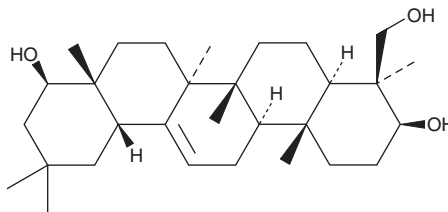
PRODUCT INFORMATION



Soyasapogenol B

Item No. 40270

CAS Registry No.: 595-15-3
Formal Name: (3 β ,4 β ,22 β)-olean-12-ene-3,22,23-triol
Synonyms: Soyasapogenin B, Soyasapogenol I
MF: C₃₀H₅₀O₃
FW: 458.7
Purity: $\geq$ 98%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years
Item Origin: Plant/*Glycine max*



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

Laboratory Procedures

Soyasapogenol B is supplied as a solid. A stock solution may be made by dissolving the soyasapogenol B in the solvent of choice, which should be purged with an inert gas. Soyasapogenol B is soluble in organic solvents such as DMSO. Soyasapogenol B is slightly soluble in chloroform and methanol.

Description

Soyasapogenol B is a triterpenoid that has been found in *A. membranaceus* and has diverse biological activities.¹⁻³ It inhibits LPS-induced increases in the levels of nitric oxide (NO) in RAW 264.7 macrophages when used at a concentration of 95.6 μ M.¹ Soyasapogenol B (12.5 μ M) decreases triglyceride levels, as well as increases glycerol and free fatty acids levels in, and decreases the levels of resistin secreted from, 3T3-L1 adipocytes.² *In vivo*, soyasapogenol B (80 mg/kg) reduces plasma levels of alanine transaminase (ALT) in a mouse model of autoimmune hepatitis induced by concanavalin A (Item No. 14951).³

References

1. Wang, Z.B., Zhai, Y.D., Ma, Z.P., *et al.* Triterpenoids and flavonoids from the leaves of *Astragalus membranaceus* and their inhibitory effects on nitric oxide production. *Chem. Biodivers.* **12**(10), 1575-1584 (2015).
2. Iwamoto, K., Kamo, S., Takada, Y., *et al.* Soyasapogenols reduce cellular triglyceride levels in 3T3-L1 mouse adipocyte cells by accelerating triglyceride lipolysis. *Biochem. Biophys. Rep.* **16**, 44-49 (2018).
3. Sasaki, K., Minowa, N., Kuzuhara, H., *et al.* Preventive effects of soyasapogenol B derivatives on liver injury in a concanavalin A-induced hepatitis model. *Bioorg. Med. Chem.* **13**(16), 4900-4911 (2005).

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, 02/08/2024

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