

PRODUCT INFORMATION



Soyasaponin Ba

Item No. 42534

CAS Registry No.: 114590-20-4

Formal Name: (3 β ,4 β ,22 β)-22,23-dihydroxyolean-12-en-3-yl O- β -D-glucopyranosyl-(1 $\rightarrow$ 2)-O- β -D-galactopyranosyl-(1 $\rightarrow$ 2)- β -D-glucopyranosiduronic acid

Synonyms: Soyasaponin M₂, Soyasaponin V

MF: C₄₈H₇₈O₁₉

FW: 959.1

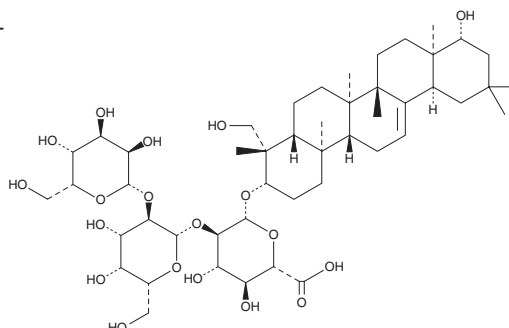
Purity: $\geq$ 98%

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

Soyasaponin Ba is supplied as a solid. A stock solution may be made by dissolving the soyasaponin Ba in the solvent of choice, which should be purged with an inert gas. Soyasaponin Ba is soluble ($\geq$ 10 mg/ml) in DMSO. Soyasaponin Ba is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Soyasaponin Ba is a saponin that has been found in *P. vulgaris* and has enzyme inhibitory and immune stimulatory activities.^{1,2} It is an inhibitor of aldose reductase 1 (K_i = 66 μ M for the recombinant human enzyme).¹ Soyasaponin Ba (50 μ g/mouse) potentiates ovalbumin-induced increases in IgG and IgG1 serum levels in mice.²

References

1. Balestri, F., De Leo, M., Sorce, C., *et al.* Soyasaponins from Zolfino bean as aldose reductase differential inhibitors. *J. Enzyme Inhib. Med. Chem.* **34**(1), 350-360 (2019).
2. Qiao, N., Liu, Q., Meng, H., *et al.* Haemolytic activity and adjuvant effect of soyasaponins and some of their derivatives on the immune responses to ovalbumin in mice. *Int. Immunopharmacol.* **18**(2), 333-339 (2014).

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, 01/24/2025

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