

PRODUCT INFORMATION



Saikosaponin B₂

Item No. 39936

CAS Registry No.: 58316-41-9
Formal Name: (3 β ,4 α ,16 α)-16,23,28-trihydroxyoleana-11,13(18)-dien-3-yl 6-deoxy-3-O- β -D-glucopyranosyl- β -D-galactopyranoside

MF: C₄₂H₆₈O₁₃

FW: 781.0

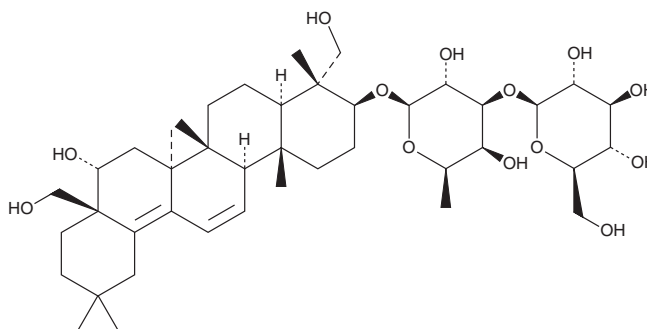
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Bupleurum falcatum*



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

Laboratory Procedures

Saikosaponin B₂ is supplied as a solid. A stock solution may be made by dissolving the saikosaponin B₂ in the solvent of choice, which should be purged with an inert gas. Saikosaponin B₂ is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

Saikosaponin B₂ is a triterpenoid saponin that has been found in *B. chinensis* and has diverse biological activities.¹⁻³ It is an antagonist of transient receptor potential ankyrin 1 (TRPA1) and inhibits calcium flux induced by polygodial (Item No. 14979) or allyl isothiocyanate in HEK293 cells expressing human TRPA1 when used at concentrations ranging from 0.5 to 1.5 μ M.¹ Saikosaponin B₂ inhibits hepatitis C virus (HCV) viral entry into, but not replication in, Huh7.5 cells (EC_{50} = 16.13 μ M) with a 50% cytotoxic concentration (CC_{50}) value of 740.4 μ M.² It protects SH-SY5Y cells from hydrogen peroxide-induced cell death when used at a concentration of 50 μ g/ml.³

References

1. Lee, G., Choi, J., Nam, Y.-J., *et al.* Identification and characterization of saikosaponins as antagonists of transient receptor potential A1 channel. *Phytother Res.* **34**(4), 788-795 (2020).
2. Lin, L.T., Chung, C.Y., Hsu, W.C., *et al.* Saikosaponin b₂ is a naturally occurring terpenoid that efficiently inhibits hepatitis C virus entry. *J. Hepatol.* **62**(3), 541-548 (2015).
3. Zhang, Y., Liu, F., Zhang, X., *et al.* Recognition and identification of active components from Radix Bupleuri using human neuroblastoma SH-SY5Y cells. *Biomed. Chromatogr.* **30**(3), 440-446 (2015).

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, 08/20/2024

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