

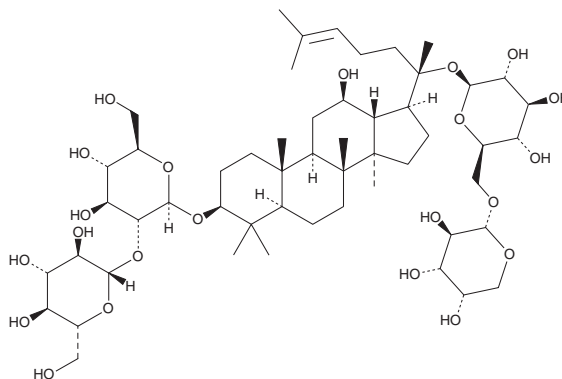
PRODUCT INFORMATION



Ginsenoside Rb₂

Item No. 22265

CAS Registry No.: 11021-13-9
Formal Name: (3 β ,12 β)-20-[(6-O- α -L-arabinopyranosyl- β -D-glucopyranosyl)oxy]-12-hydroxydammar-24-en-3-yl 2-O- β -D-glucopyranosyl- β -D-glucopyranoside
Synonyms: Ginsenoside C, NSC 308878
MF: C₅₃H₉₀O₂₂
FW: 1,079.3
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Ginsenoside Rb₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the ginsenoside Rb₂ in the solvent of choice. Ginsenoside Rb₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ginsenoside Rb₂ in these solvents is approximately 0.1, 10, and 15 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 ginsenoside Rb₂ can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ginsenoside Rb₂ in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ginsenoside Rb₂ is a steroid glycoside found in plants of the genus *Panax* that has diverse biological activities.¹ It inhibits LPS-stimulated increases in vascular cell adhesion molecule 1 (VCAM-1) and intercellular adhesion molecule 1 (ICAM-1) expression *in vitro*. It dose-dependently reverses LPS-induced decreases in the expression of I κ B α in human umbilical vein endothelial cells (HUVECs).² Ginsenoside Rb₂ reduces adhesion of THP-1 cells to LPS-stimulated HUVEC cells, also in a dose-dependent manner. Ginsenoside Rb₂ reduces UVB-induced cytotoxicity and apoptotic nuclear fragmentation in HaCaT cells.³ It also lowers total cholesterol and triacylglycerol in 3T3-L1 adipocytes cultured under high fatty acid conditions to levels comparable to lovastatin (Item No. 10010338).⁴

References

1. Liang, Y. and Zhao, S. *Plant Biol.(Stuttg)* **10**(4), 415-421 (2008).
2. Cho, Y.S., Kim, C.H., Ha, T.S., *et al.* *Korean J.Physiol.Pharmacol.* **17**(2), 133-137 (2013).
3. Ha, S.E., Shin, D.H., Kim, H.D., *et al.* *Naunyn Schmiedebergs Arch. Pharmacol.* **382**(1), 89-101 (2010).
4. Kim, E.J., Lee, H.I., Chung, K.J., *et al.* *BMB Rep.* **42**(4), 194-199 (2009).

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, 12/02/2022

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