

PRODUCT INFORMATION

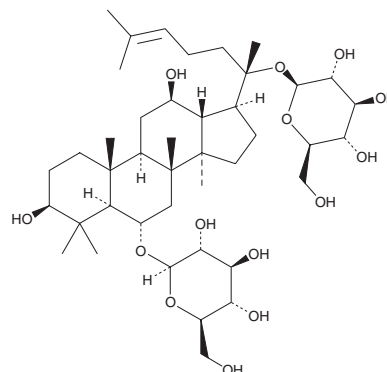


Ginsenoside Rg₁

Item No. 15315

CAS Registry No.: 22427-39-0
Formal Name: 3β,12β-dihydroxydammar-24-ene-6α,20-diyl bis-β-D-glucopyranoside
Synonyms: Ginsenoside A₂, Panaxoside A, Panaxoside Rg₁, Sanchinoside C1, Sanchinoside Rg₁

MF: C₄₂H₇₂O₁₄
FW: 801.0
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/Ginseng Radix et Rhizoma



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

Laboratory Procedures

Ginsenoside Rg₁ is supplied as a crystalline solid. A stock solution may be made by dissolving the ginsenoside Rg₁ in the solvent of choice, which should be purged with an inert gas. Ginsenoside Rg₁ is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ginsenoside Rg₁ in these solvents is approximately 10 mg/ml.

Ginsenoside Rg₁ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ginsenoside Rg₁ should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Ginsenoside Rg₁ has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ginsenoside Rg₁ is a steroid glycoside that has been found in *P. ginseng* and has diverse biological activities.¹⁻⁴ *In vivo*, ginsenoside Rg₁ (10, 20, and 40 mg/kg) decreases blood glucose, IL-1β, and IL-18 levels, increases insulin secretion, and inhibits hepatic and pancreatic activation of the NLRP3 inflammasome in a mouse model of diabetes induced by streptozotocin (STZ; Item No. 13104).¹ It facilitates fear extinction and decreases immobility time in the tail-suspension and forced swim tests in a mouse model of single-prolonged stress-induced post-traumatic stress disorder (PTSD).² Ginsenoside Rg₁ decreases hepatic alanine transaminase (ALT) and aspartate aminotransferase (AST) levels and attenuates hepatic steatosis in a mouse model of ethanol-induced hepatitis.³ It also decreases the abundance of *Bacteroidetes* in the large intestine, cortical neuron apoptosis, and latency to the platform in the Morris water maze in a tree shrew model of Alzheimer's disease induced by amyloid-β (25-35) (Aβ (25-35)).⁴

References

1. Gao, Y., Li, J., Chu, S., *et al.* Ginsenoside Rg₁ protects mice against streptozotocin-induced type 1 diabetic by modulating the NLRP3 and Keap1/Nrf2/HO-1 pathways. *Eur. J. Pharmacol.* **866**, 172801 (2020).
2. Zhang, Z., Song, Z., Shen, F., *et al.* Ginsenoside Rg₁ prevents PTSD-like behaviors in mice through promoting synaptic proteins, reducing Kir4.1 and TNF-α in the hippocampus. *Mol. Neurobiol.* **58(4)**, 1550-1563 (2021).
3. Yang, C., He, X., Zhao, J., *et al.* Hepatoprotection by Ginsenoside Rg₁ in alcoholic liver disease. *Int. Immunopharmacol.* **92**, 107327 (2021).
4. Guo, Y., Wang, L., Lu, J., *et al.* Ginsenoside Rg₁ improves cognitive capability and affects the microbiota of large intestine of tree shrew model for Alzheimer's disease. *Mol. Med. Rep.* **23(4)**, 291 (2021).

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