

PRODUCT INFORMATION



Ginsenoside Rg₂

Item No. 15331

CAS Registry No.: 52286-74-5
Formal Name: (6 α)-3 β ,12 β ,20-trihydroxydammar-24-en-6-yl 2-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranoside

Synonyms: Chikusetsusaponin I, Panaxoside Rg₂, Prosapogenin C₂

MF: C₄₂H₇₂O₁₃

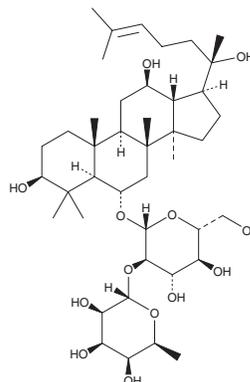
FW: 785.0

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 Rg₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the ginsenoside Rg₂ in the solvent of choice. Ginsenoside Rg₂ is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of ginsenoside Rg₂ in these solvents is approximately 10 mg/ml.

Ginsenoside Rg₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ginsenoside Rg₂ should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ginsenoside Rg₂ has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ginsenosides are pharmacologically active natural compounds from ginseng and other plants of the genus *Panax*.¹ Structurally, they are steroid glycosides derived from the triterpene squalene.¹ Ginsenoside Rg₂ is a protopanaxatriol that is more abundant in some *Panax* species (e.g., white and red *P. ginseng*) than others.² This ginsenoside and its metabolites have diverse *in vitro* and *in vivo* effects, including neuroprotective, anti-inflammatory, and anti-diabetic actions.³⁻⁵ It also protects against DNA damage and apoptosis induced by ultraviolet light.⁶ Notably, this ginsenoside is increased by the metabolism of other bioactive ginsenosides during the steaming or heating of plant materials, particularly in *P. quinquefolium*.^{2,7}

References

1. Liang, Y. and Zhao, S. *Plant Biol. (Stuttg)* **10(4)**, 415-421 (2008).
2. Chen, C.-F., Chiou, W.-F., and Zhang, J.-T. *Acta. Pharmacol. Sin.* **29(9)**, 1103-1108 (2008).
3. Zhang, G., Liu, A., Zhou, Y., et al. *J. Ethnopharmacol.* **115(3)**, 441-448 (2008).
4. Cho, Y.S., Kim, C.H., Ha, T.S., et al. *Korean J. Physiol. Pharmacol.* **17(2)**, 133-137 (2013).
5. Yuan, H.D., Kim, D.Y., Quan, H.Y., et al. *Chem. Biol. Interact.* **195(1)**, 35-42 (2012).
6. Ha, S.E., Shin, D.H., Kim, H.D., et al. *Naunyn Schmiedebergs Arch. Pharmacol.* **382(1)**, 89-101 (2010).
7. Kim, I.-W., Cha, K.-M., Wee, J.J., et al. *J. Ginseng Res.* **37(4)**, 475-482 (2013).

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, 11/28/2022

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