

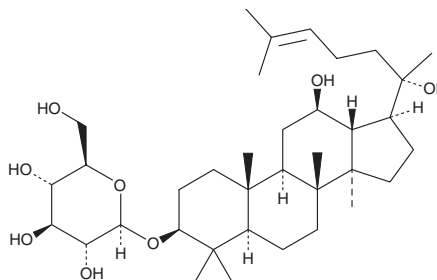
PRODUCT INFORMATION



20(S)-Ginsenoside Rh₂

Item No. 23390

CAS Registry No.: 78214-33-2
Formal Name: 12 β ,20-dihydroxydammar-24-en-3 β -yl, β -D-glucopyranoside
Synonym: 20(S)-Ginsenoside Rh₂
MF: C₃₆H₆₂O₈
FW: 622.9
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 Rh₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the ginsenoside Rh₂ in the solvent of choice, which should be purged with an inert gas. Ginsenoside Rh₂ is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ginsenoside Rh₂ in these solvents is approximately 10 mg/ml.

Ginsenoside Rh₂ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ginsenoside Rh₂ should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ginsenoside Rh₂ 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

Ginsenoside Rh₂ is a steroid glycoside found in plants of the genus *Panax* that has diverse biological activities.¹⁻⁵ It inhibits release of β -hexosaminidase from RBL-2H3 cells (IC₅₀ = 100 μ M) and inhibits the IgE-dependent passive cutaneous anaphylaxis reaction in mice at a dose of 25 mg/kg.¹ It inhibits growth of HRA ovarian and BxPC-3 pancreatic cancer cells in a dose-dependent manner.^{3,4} *In vivo*, ginsenoside Rh₂ decreases immobility time in the forced swim test in a mouse model of colorectal carcinoma.⁵ Ginsenoside Rh₂ also reduces infarct volume in a rat model of ischemia-reperfusion-induced brain injury.²

References

1. Park, E.-K., Choo, M.-K., Kim, E.-J., *et al.* Antiallergic activity of ginsenoside Rh₂. *Biol. Pharm. Bull.* **26**(11), 1581-1584 (2003).
2. Park, E.-K., Choo, M.-K., Oh, J.K., *et al.* Ginsenoside Rh₂ reduces ischemic brain injury in rats. *Biol. Pharm. Bull.* **27**(3), 433-436 (2004).
3. Kikuchi, Y., Sasa, H., Kita, T., *et al.* Inhibition of human ovarian cancer cell proliferation *in vitro* by ginsenoside Rh₂ and adjuvant effects to cisplatin *in vivo*. *Anticancer Drugs* **2**(1), 63-67 (1991).
4. Tang, X.-P., Tang, G.-D., Fang, C.-Y., *et al.* Effects of ginsenoside Rh₂ on growth and migration of pancreatic cancer cells. *World J. Gastroenterol.* **19**(10), 1582-1592 (2013).
5. Wang, J., Chen, Y., Dai, C.L., *et al.* Ginsenoside Rh₂ alleviates tumor-associated depression in a mouse model of colorectal carcinoma. *Am. J. Transl. Res.* **8**(5), 2189-2195 (2016).

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, 12/05/2022

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