

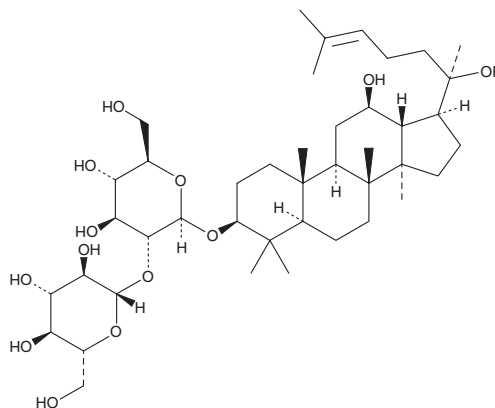
PRODUCT INFORMATION



20(R)-Ginsenoside Rg₃

Item No. 24978

CAS Registry No.: 38243-03-7
Formal Name: 12β,20R-dihydroxydammar-24-en-3β-yl 2-O-β-D-glucopyranosyl-β-D-glucopyranoside
Synonym: 20(R)-Propanaxadiol
MF: C₄₂H₇₂O₁₃
FW: 785.0
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

20(R)-Ginsenoside Rg₃ is supplied as a crystalline solid. A stock solution may be made by dissolving the 20(R)-ginsenoside Rg₃ in the solvent of choice, which should be purged with an inert gas. 20(R)-Ginsenoside Rg₃ is slightly soluble in ethanol.

Description

20(R)-Ginsenoside Rg₃ is a saponin that has been found in *P. ginseng* and inhibits angiogenesis.¹ 20(R)-Ginsenoside Rg₃ (0.001-1 μM) dose-dependently reduces proliferation, capillary tube branching, and the expression of matrix metalloproteinase-9 (MMP-9) and MMP-2 in human umbilical vein endothelial cells (HUVECs). It reduces tumor cell invasion of MM1 rat hepatoma, B16FE7 mouse melanoma, OC-10 human lung carcinoma, and PSN-1 human pancreatic adenocarcinoma cells *in vitro* by 89, 73.7, 84.2, and 59.1%, respectively, when used at a concentration of 25 μM.² However, it does not inhibit the proliferation of MM1 cells when used at concentrations ranging from 1.6 to 64 μM. 20(R)-Ginsenoside Rg₃ (20 mg/kg) reduces tumor volume, increases the tumor necrosis rate by approximately 16%, and increases survival in a human Lewis lung carcinoma mouse xenograft model.³ It also reduces the total number of B16-BL6 melanoma cell and 26-M3.1 colon carcinoma cell lung colonies in mice by 27 and 51%, respectively, when administered at a dose of 100 μg.⁴ Topical administration of 20(R)-ginsenoside Rg₃ (0.05% w/v) reduces ear thickness by 47.5% compared with control animals in a mouse model of oxazolone-induced contact dermatitis and reduces expression of COX-2, IL-4, IL-1, IFN, and TNF mRNA by greater than 25%.⁵

References

1. Yue, P.Y., Wong, D.Y., Wu, P.K., *et al.* The angiosuppressive effects of 20(R)- ginsenoside Rg₃. *Biochem. Pharmacol.* **72**(4), 437-445 (2006).
2. Shinkai, K., Akedo, H., Mukai, M., *et al.* Inhibition of *in vitro* tumor cell invasion by ginsenoside Rg₃. *Jpn. J. Cancer Res.* **87**(4), 357-362 (1996).
3. Liu, T.G., Huang, Y., Cui, D.D., *et al.* Inhibitory effect of ginsenoside Rg₃ combined with gemcitabine on angiogenesis and growth of lung cancer in mice. *BMC Cancer* **9**, 250 (2009).
4. Mochizuki, M., Yoo, Y.C., Matsuzawa, K., *et al.* Inhibitory effect of tumor metastasis in mice by saponins, ginsenoside-Rb2, 20(R)- and 20(S)-ginsenoside-Rg3, of red ginseng. *Biol. Pharm. Bull.* **18**(9), 1197-1202 (1995).
5. Bae, E.A., Han, M.J., Shin, Y.W., *et al.* Inhibitory effects of Korean red ginseng and its genuine constituents ginsenosides Rg₃, Rf, and Rh2 in mouse passive cutaneous anaphylaxis reaction and contact dermatitis models. *Biol. Pharm. Bull.* **29**(9), 1862-1867 (2006).

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, 12/21/2022

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