

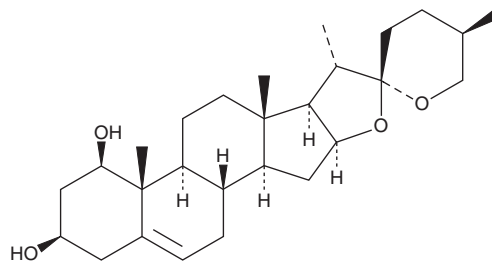
PRODUCT INFORMATION



Ruscogenin

Item No. 25217

CAS Registry No.: 472-11-7
Formal Name: (1 β ,3 β ,25R)-spirost-5-ene-1,3-diol
MF: C₂₇H₄₂O₄
FW: 430.6
Purity: $\geq 95\%$
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

Ruscogenin is supplied as a crystalline solid. A stock solution may be made by dissolving the ruscogenin in the solvent of choice, which should be purged with an inert gas. Ruscogenin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ruscogenin in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

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

Ruscogenin is a steroid sapogenin that has been found in *O. japonicus* with diverse biological activities.¹⁻⁵ It reduces adhesion of HL-60 cells to ECV304 cells induced by phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) with an IC₅₀ value of 7.76 nM and decreases peritoneal leukocyte migration induced by zymosan A (Item No. 21175) in mice when administered at a dose of 5 mg/kg.¹ Oral administration of ruscogenin (12.5 mg/kg) inhibits ADP-induced platelet aggregation in rats.² Ruscogenin prevents monocrotaline-induced increases in mean arterial pressure (MAP), expression of IL-1 β and TNF in plasma and lung, and endothelial cell apoptosis in pulmonary arterioles in a rat model of pulmonary hypertension.³ It also reduces proteinuria and renal macrophage infiltration in a rat model of diabetes induced by streptozotocin (Item No. 13104) and decreases gene expression of inflammatory cytokines, dyslipidemia, and liver steatosis in a hamster model of high-fat diet-induced nonalcoholic steatohepatitis (NASH).^{4,5}

References

1. Kou, J., Sun, Y., Lin, Y., et al. Anti-inflammatory activities of aqueous extract from *Radix Ophiopogon japonicus* and its two constituents. *Biol. Pharm. Bull.* **28(7)**, 1234-1238 (2005).
2. Kou, J., Tian, Y., Yan, J., et al. Antithrombotic activities of aqueous extract from *Radix Ophiopogon japonicus* and its two constituents. *Biol. Pharm. Bull.* **29(6)**, 1267-1270 (2006).
3. Bi, L.-Q., Zhu, R., Kong, H., et al. Ruscogenin attenuates monocrotaline-induced pulmonary hypertension in rats. *Int. Immunopharmacol.* **16(1)**, 7-16 (2013).
4. Lu, H.-H., Tzeng, T.-F., Liou, S.-S., et al. Ruscogenin ameliorates diabetic nephropathy by its anti-inflammatory and anti-fibrotic effects in streptozotocin-induced diabetic rat. *BMC Complement. Altern. Med.* **14:110**, (2014).
5. Lu, H.-J., Tzeng, T.-F., Liou, S.-S., et al. Ruscogenin ameliorates experimental nonalcoholic steatohepatitis via suppressing lipogenesis and inflammatory pathway. *Biomed. Res. Int.* 652680 (2014).

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, 10/17/2024

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