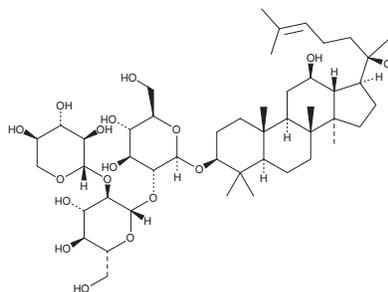


PRODUCT INFORMATION



Notoginsenoside Ft₁ Item No. 24976

CAS Registry No.: 155683-00-4
Formal Name: (3β,12β,20R)-12,20-dihydroxydammar-24-en-3-yl O-β-D-xylopyranosyl-(1→2)-O-β-D-glucopyranosyl-(1→2)-β-D-glucopyranoside
MF: C₄₇H₈₀O₁₇
FW: 917.1
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

Notoginsenoside Ft₁ is supplied as a crystalline solid. A stock solution may be made by dissolving the notoginsenoside Ft₁ in the solvent of choice, which should be purged with an inert gas. Notoginsenoside Ft₁ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of notoginsenoside Ft₁ in these solvents is approximately 5, 15, and 20 mg/ml, respectively.

Notoginsenoside Ft₁ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, notoginsenoside Ft₁ should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Notoginsenoside Ft₁ has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Notoginsenoside Ft₁ is a saponin originally isolated from *P. notoginseng* with diverse biological activities.¹⁻⁵ It induces proliferation, migration, and tube formation of human umbilical vein endothelial cells (HUVECs) via nuclear translocation of hypoxia-inducible factor-1α (HIF-1α) and activation of the PI3K/AKT and Raf/MEK/ERK signaling pathways in a manner dependent on mammalian target of rapamycin (mTOR).¹ Notoginsenoside Ft₁ (45 μM) induces cell cycle arrest at the S and G₂/M phases and promotes apoptosis of SH-SY5Y cells.² It increases cGMP levels and induces relaxation of isolated precontracted rat mesenteric arteries, effects that are reversed by the nitric oxide synthase inhibitor L-NAME (Item No. 80210) and ODQ (Item No. 81410), an inhibitor of soluble guanylyl cyclase.³ *In vivo*, notoginsenoside Ft₁ (0.25, 2.5, and 25 mg/kg) promotes angiogenesis and decreases wound diameter in a mouse model of punched-hole ear injury.¹ Notoginsenoside Ft₁ (1.25 mg/kg) decreases tail bleeding time and increases thrombus weight in a rat tail bleeding assay.⁴ Topical administration of notoginsenoside Ft₁ increases mRNA expression of the collagen expression, fibroblast proliferation, and scar formation genes COL1A1, COL3A1, TGF-β1, TGF-β3, and fibronectin, promotes neovascularization, reduces monocyte infiltration, and shortens wound closure time in a *db/db* mouse model of diabetic foot ulcers.⁵

References

1. Shen, K., Ji, L., Gong, C., et al. *Biochem. Pharmacol.* **84**(6), 784-792 (2012).
2. Gao, B., Shi, H.-L., Li, X., et al. *Life Sci.* **108**(2), 63-70 (2014).
3. Shen, K., Leung, S.W.S., Ji, L., et al. *Biochem. Pharmacol.* **88**(1), 66-74 (2014).
4. Gao, B., Huang, L., Liu, H., et al. *Br. J. Pharmacol.* **171**(1), 214-223 (2014).
5. Zhang, E., Gao, B., Yang, L., et al. *J. Pharmacol. Exp. Ther.* **356**(2), 324-332 (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, 11/18/2022

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