

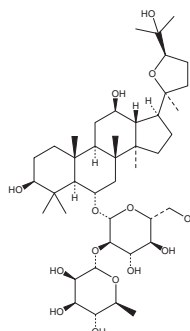
PRODUCT INFORMATION



Pseudoginsenoside F₁₁

Item No. 30222

CAS Registry No.: 69884-00-0
Formal Name: (3 β ,6 α ,12 β ,24R)-20,24-epoxy-3,12,25-trihydroxydammaran-6-yl 2-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranoside
Synonym: Ginsenoside A₁
MF: C₄₂H₇₂O₁₄
FW: 801.0
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Panax ginseng*



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

Laboratory Procedures

Pseudoginsenoside F₁₁ is supplied as a solid. A stock solution may be made by dissolving the pseudoginsenoside F₁₁ in the solvent of choice, which should be purged with an inert gas. Pseudoginsenoside F₁₁ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of pseudoginsenoside F₁₁ in these solvents is approximately 0.1, 10, and 15 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of pseudoginsenoside F₁₁ can be prepared by directly dissolving the solid in aqueous buffers. The solubility of pseudoginsenoside F₁₁ in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pseudoginsenoside F₁₁ is an ocotillol-type ginsenoside that has been found in *P. ginseng* and has diverse biological activities.¹⁻⁴ *In vivo*, pseudoginsenoside F₁₁ (10 mg/kg) prevents tubular cell apoptosis, decreases in renal glutathione peroxidase (GPX) and superoxide dismutase (SOD) levels, and increases in renal lipid peroxide levels in a rat model of nephrotoxicity induced by cisplatin (Item No. 13119).¹ It reduces infarct size, brain water content, and cortical accumulation of autophagosomes in a rat model of ischemic stroke induced by permanent middle cerebral artery occlusion.² Pseudoginsenoside F₁₁ (4 and 8 mg/kg) inhibits morphine-induced memory impairment in the Morris water maze and development of morphine-induced conditioned place preference in mice.³ It also reduces hippocampal advanced glycation end product (AGE) and malondialdehyde (MDA) levels, increases hippocampal SOD activity and glutathione (GSH) levels, and attenuates cognitive impairment in the Morris water maze in a mouse model of D-galactose-induced mild cognitive impairment.⁴

References

1. Wang, H., Kong, L., Zhang, J., et al. *Sci. Rep.* **4**, 4986 (2014).
2. Liu, Y.Y., Zhang, T.Y., Xue, X., et al. *CNS Neurosci. Ther.* **23(7)**, 567-579 (2017).
3. Li, Z., Wu, C.F., Pei, G., et al. *Pharmacol. Biochem. Behav.* **66(3)**, 595-601 (2000).
4. Zhang, Z., Yang, H., Yang, J., et al. *Int. Immunopharmacol.* **67**, 78-86 (2019).

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, 11/07/2022

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