

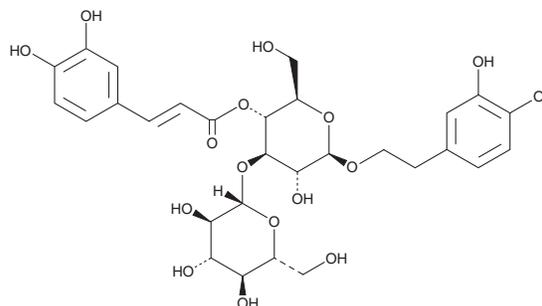
PRODUCT INFORMATION



Plantamajoside

Item No. 27296

CAS Registry No.: 104777-68-6
Formal Name: 2-(3,4-dihydroxyphenyl)ethyl 3-O-β-D-glucopyranosyl-β-D-glucopyranoside 4-[(2E)-3-(3,4-dihydroxyphenyl)-2-propenoate]
MF: C₂₉H₃₆O₁₆
FW: 640.6
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 334 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Plantago asiatica* L.



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

Laboratory Procedures

Plantamajoside is supplied as a crystalline solid. A stock solution may be made by dissolving the plantamajoside in the solvent of choice, which should be purged with an inert gas. Plantamajoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of plantamajoside in these solvents is approximately 10, 33, and 50 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 plantamajoside can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of plantamajoside in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Plantamajoside is a phenylethanoid that has been found in *P. lanceolata* and has diverse biological activities, including antioxidative, anticancer, and anti-inflammatory properties.¹⁻³ It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in a cell-free assay (IC₅₀ = 11.8 μM). It also inhibits 5-lipoxygenase (5-LO) and 15-LO (IC₅₀s = 0.375 and 96 μM, respectively).^{1,2} Plantamajoside decreases viability of TE-1 esophageal squamous cell carcinoma (ESCC) cells when used at concentrations ranging from 20 to 320 μg/ml, as well as inhibits LPS-induced epithelial-to-mesenchymal transition (EMT) in Eca-109 and TE-1 cells and decreases phosphorylated NF-κB levels in TE-1 cells.³ Plantamajoside (3 mg/ear) inhibits mouse ear edema induced by arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) by 25%.¹

References

1. Murai, M., Tamayama, Y., and Nishibe, S. *Planta. Med.* **61**(5), 479-480 (1995).
2. Skari, K.P., Malterud, K.E., and Haugli, T. *Natural antioxidants and anticarcinogens in nutrition, health and disease.* Kumpulainen, J.T. and Salonen, J.T., editors, *The Royal Society of Chemistry* (1999).
3. Li, X., Chen, D., Li, M., et al. *Biomed. Pharmacother.* **102**, 1045-1051 (2018).

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, 12/19/2022

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