

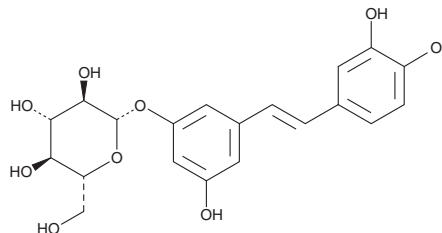
PRODUCT INFORMATION



Astringin

Item No. 27979

CAS Registry No.: 29884-49-9
Formal Name: 3-[(1E)-2-(3,4-dihydroxyphenyl)ethenyl]-5-hydroxyphenyl β-D-glucopyranoside
Synonyms: (E)-Astringin, *trans*-Astringin, 3,4,3',5'-Tetrahydroxystilbene 3'-glucoside
MF: C₂₀H₂₂O₉
FW: 406.4
Purity: ≥90%
UV/Vis.: λ_{max}: 219, 322 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Rheum hotaoense*



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

Laboratory Procedures

Astringin is supplied as a crystalline solid. A stock solution may be made by dissolving the astringin in the solvent of choice, which should be purged with an inert gas. Astringin is soluble in DMSO.

Description

Astringin is a phenolic stilbene glucoside that has been found in *V. vinifera* and has antioxidant and antineoplastic activities.^{1,2} It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals (IC₅₀ = 30.2 μM) in a cell-free assay and inhibits cupric ion-induced lipid peroxidation of human LDL (IC₅₀ = 3 μM).¹ Astringin (10 μg/ml) inhibits development of preneoplastic lesions induced by 7,12-dimethylbenz(a)anthracene (DMBA) in mouse mammary gland organ cultures by 68.8%.²

References

1. Tegeo, P.W., Fauconneau, B., Deffieux, G., *et al.* Isolation, identification, and antioxidant activity of three stilbene glucosides newly extracted from *Vitis vinifera* cell cultures. *J. Nat. Prod.* **61**(5), 655-657 (1998).
2. Waffo-Tégou, P., Hawthorne, M.E., Cuendet, M., *et al.* Potential cancer-chemopreventive activities of wine stilbenoids and flavans extracted from grape (*Vitis vinifera*) cell cultures. *Nutr. Cancer* **40**(2), 173-179 (2001).

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, 09/30/2022

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