

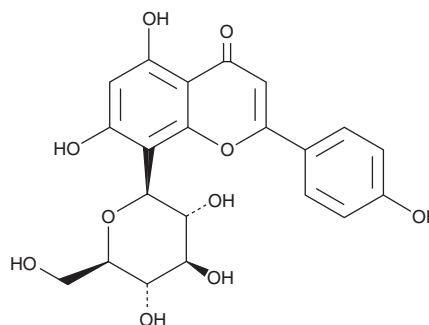
PRODUCT INFORMATION



Vitexin

Item No. 15116

CAS Registry No.: 3681-93-4
Formal Name: 8-β-D-glucopyranosyl-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
Synonym: Orientoside
MF: C₂₁H₂₀O₁₀
FW: 432.4
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 270, 332 nm
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

Vitexin is supplied as a crystalline solid. A stock solution may be made by dissolving the vitexin in the solvent of choice, which should be purged with an inert gas. Vitexin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of vitexin in these solvents is approximately 16.6 and 14.3 mg/ml, respectively.

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

Vitexin is an apigenin flavone C-glycoside isolated from *Crataegus* spp. (hawthorn), *Passiflora* spp. (passionflower), *P. glaucum* R. Br. (pearl millet), *P. nigra* (bamboo leaves), *V. agnuscastus* L. (chasteberry), and various other tropical plant species.¹⁻³ It demonstrates antioxidant activity in a DPPH bleaching assay (IC₅₀ = 24.2 μM) and also inhibits the activity of the carbohydrate-hydrolyzing enzyme α-glucosidase (IC₅₀ = 244 μM).^{1,3} It has been shown to have broad anti-tumor activity, activating caspases and inducing apoptosis in cancer xenograft models.⁴ Additionally, vitexin at 10 mg/kg exhibits an analgesic effect in several different models of inflammatory pain by targeting TRPV1 channel activity, by preventing the decrease of reduced glutathione levels, and by modulating the production of cytokines.⁵

References

1. Simirgiotis, M.J., Schmeda-Hirschmann, G., Bórquez, J., *et al.* The *Passiflora tripartita* (banana passion) fruit: A source of bioactive flavonoid C-glycosides isolated by HSCCC and characterized by HPLC-DAD-ESI/MS/MS. *Molecules* **18**(2), 1672-1692 (2013).
2. Wang, J., Yue, Y.-D., Jiang, H., *et al.* Rapid screening for flavone C-glycosides in the leaves of different species of bamboo and simultaneous quantitation of four marker compounds by HPLC-UV/DAD. *Int. J. Anal. Chem.* 205101 (2012).
3. Chen, Y.-G., Ping, L., Peng, L., *et al.* α-Glucosidase inhibitory effect and simultaneous quantification of three major flavonoid glycosides in *Microctis folium*. *Molecules* **18**(4), 4221-4232 (2013).
4. Zhou, Y., Liu, Y.E., Cao, J., *et al.* Vitexins, nature-derived lignan compounds, induce apoptosis and suppress tumor growth. *Clin. Cancer Res.* **15**(16), 5161-5169 (2009).
5. Borghi, S.M., Carvalho, T.T., Staurengo-Ferrari, L., *et al.* Vitexin inhibits inflammatory pain in mice by targeting TRPV1, oxidative stress, and cytokines. *J. Nat. Prod.* **76**(6), 1141-1149 (2013).

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, 12/20/2022

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