

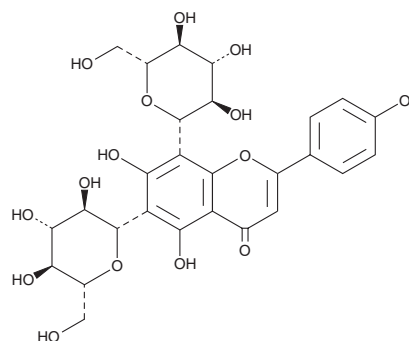
# PRODUCT INFORMATION



## Vicenin-2

Item No. 33403

**CAS Registry No.:** 23666-13-9  
**Formal Name:** 6,8-di-β-D-glucopyranosyl-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one  
**Synonyms:** Vicenin II, Violantin  
**MF:** C<sub>27</sub>H<sub>30</sub>O<sub>15</sub>  
**FW:** 594.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 218, 274, 335 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Desmodium styracifolium*



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

### Laboratory Procedures

Vicenin-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the vicenin-2 in the solvent of choice, which should be purged with an inert gas. Vicenin-2 is soluble in methanol.

### Description

Vicenin-2 is a polyketide synthase-derived flavonoid glycoside that has been found in *C. subternata* and has diverse biological activities.<sup>1-5</sup> It inhibits thrombin-induced aggregation of isolated mouse platelets when used at concentrations ranging from 2 to 30 μM, as well as increases tail bleeding time in mice when administered at a dose of 35.7 μg/animal.<sup>2</sup> Vicenin-2 (1 mg/kg) reduces tumor growth in a PC3 prostate cancer mouse xenograft model.<sup>3</sup> It decreases trabecular separation, as well as increases trabecular thickness and area, in an ovariectomized rat model of osteoporosis.<sup>4</sup> Vicenin-2 (50 mg/kg) decreases ulceration and increases colon weight and length in a mouse model of colitis induced by dextran sulfate (DSS; Item No. 23250).<sup>5</sup>

### References

1. Kitamura, K., Ando, Y., Matsumoto, T., *et al.* Total synthesis of aryl C-glycoside natural products: Strategies and tactics. *Chem. Rev.* **118**(4), 1495-1598 (2018).
2. Lee, W. and Bae, J.-S. Antithrombotic and antiplatelet activities of vicenin-2. *Blood Coagul. Fibrinolysis* **26**(6), 628-634 (2015).
3. Nagaprashantha, L.D., Vatsyayan, R., Singhal, J., *et al.* Anti-cancer effects of novel flavonoid vicenin-2 as a single agent and in synergistic combination with docetaxel in prostate cancer. *Biochem. Pharmacol.* **82**(9), 1100-1109 (2011).
4. Zhang, Z., Zhao, Q., Liu, T., *et al.* Effect of vicenin-2 on ovariectomy-induced osteoporosis in rats. *Biomed. Pharmacother.* **129**, 110474 (2020).
5. Yin, Y., Ye, L., Niu, Z., *et al.* Anti-inflammatory effects of vicenin-2 on dextran sulfate sodium-induced colitis in mice. *Drug Dev. Res.* **80**(5), 546-555 (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, 12/07/2022

#### 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](http://WWW.CAYMANCHEM.COM)