

# PRODUCT INFORMATION

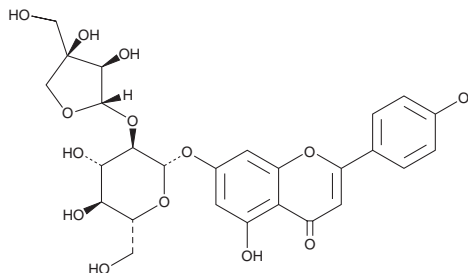


## Apiin

Item No. 34596

**CAS Registry No.:** 26544-34-3  
**Formal Name:** 7-[(2-O-D-apio-β-D-furanosyl-β-D-glucopyranosyl)oxy]-5-hydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one  
**Synonyms:** Apigenin 7-apisoylglucoside, Apigenin 7-O-apisoylglucoside

**MF:** C<sub>26</sub>H<sub>28</sub>O<sub>14</sub>  
**FW:** 564.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 269, 338 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Apium graveolens*



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

### Laboratory Procedures

Apiin is supplied as a solid. A stock solution may be made by dissolving the apiin in the solvent of choice, which should be purged with an inert gas. Apiin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of apiin in these solvents is approximately 30 mg/ml.

Apiin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, apiin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Apiin has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Apiin is a polyketide synthase-derived flavonoid glycoside that has been found in *A. graveolens* and has diverse biological activities.<sup>1-4</sup> It inhibits Aurora B kinase in a cell-free assay using the human enzyme (IC<sub>50</sub> = 12.14 μM).<sup>2</sup> Apiin (10 and 50 μg/ml) inhibits LPS-induced increases in nitric oxide (NO) release and inducible nitric oxide synthase (iNOS) levels in J774.A1 macrophages.<sup>3</sup> It scavenges DPPH (Item No. 14805), superoxide, and hydroxyl radicals (IC<sub>50</sub>s = 68, 390, and 48 μg/ml, respectively), as well as increases the levels of superoxide dismutase (SOD) and catalase in mouse serum, brain, heart, liver, and kidney when administered at a dose of 50 mg/kg.<sup>4</sup>

### References

- Cheung, V.W.N., Xue, B., Hernandez-Valladares, M., *et al.* Identification of polyketide inhibitors targeting 3-dehydroquininate dehydratase in the shikimate pathway of *Enterococcus faecalis*. *PLoS One* **9**(7), e103598 (2014).
- Jung, Y., Shin, S.Y., Yong, Y., *et al.* Plant-derived flavones as inhibitors of aurora B kinase and their quantitative structure-activity relationships. *Chem. Biol. Drug Des.* **85**(5), 574-585 (2015).
- Mencherini, T., Cau, A., Bianco, G., *et al.* An extract of *Apium graveolens* var. dulce leaves: Structure of the major constituent, apiin, and its anti-inflammatory properties. *J. Pharm. Pharmacol.* **59**(6), 891-897 (2007).
- Li, P., Jia, J., Zhang, D., *et al.* *In vitro* and *in vivo* antioxidant activities of a flavonoid isolated from celery (*Apium graveolens* L. var. dulce). *Food Funct.* **5**(1), 50-56 (2014).

#### 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, 12/06/2022

#### 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