

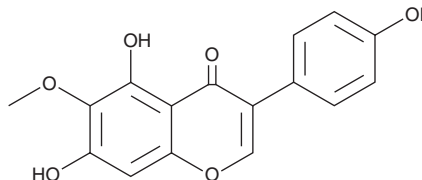
# PRODUCT INFORMATION



## Tectorigenin

Item No. 35860

**CAS Registry No.:** 548-77-6  
**Formal Name:** 5,7-dihydroxy-3-(4-hydroxyphenyl)-6-methoxy-4H-1-benzopyran-4-one  
**Synonym:** Tectorigenine  
**MF:** C<sub>16</sub>H<sub>12</sub>O<sub>6</sub>  
**FW:** 300.3  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Belamcanda chinensis*



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

### Laboratory Procedures

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

### Description

Tectorigenin is an isoflavonoid that has been found in *B. chinensis* and has diverse biological activities and is an active metabolite of the flavonoid glycoside and phytoestrogen tectoridin (Item No. 33969).<sup>1-5</sup> It is formed from tectoridin by gut microbiota.<sup>1</sup> Tectorigenin scavenges hydroxyl and superoxide radicals in cell-free assays (IC<sub>50</sub>s = 87 and 46.62 μg/ml, respectively) and inhibits the formation of thiobarbituric acid reactive substances (TBARS) in egg yolk (IC<sub>50</sub> = 23 μg/ml).<sup>3</sup> It inhibits the production of prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) induced by the PKC activator 12-O-tetradecanoylphorbol 13-acetate (TPA; Item No. 10008014) or the sarcoplasmic/endoplasmic reticulum Ca<sup>2+</sup>-ATPase (SERCA) and endoplasmic reticulum (ER) stress inducer thapsigargin (Item No. 10522) in isolated rat peritoneal macrophages when used at concentrations of 3, 10, or 30 μM.<sup>3</sup> Tectorigenin (10 μg/ml) reduces cell death induced by hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) and increases catalase levels and activity in V79-4 Chinese hamster lung fibroblasts.<sup>4</sup> *In vivo*, tectorigenin (100 mg/kg per day) prevents increases in hepatic levels of malondialdehyde (MDA) and blood levels of alanine transaminase (ALT) and aspartate aminotransferase (AST) and decreases in hepatic superoxide dismutase (SOD), catalase, and glutathione peroxidase (GPX) activities in a rat model of liver injury induced by carbon tetrachloride (CCl<sub>4</sub>).<sup>5</sup>

### References

1. Chen, Y., Song, W., Peng, Z.H., *et al.* Identification of metabolites of tectoridin in-vivo and in-vitro by liquid chromatography-tandem mass spectrometry. *J. Pharm. Pharmacol.* **60(6)**, 709-716 (2008).
2. Han, T., Cheng, G., Liu, Y., *et al.* *In vitro* evaluation of tectoridin, tectorigenin and tectorigenin sodium sulfonate on antioxidant properties. *Food Chem. Toxicol.* **50(2)**, 409-414 (2012).
3. Kim, Y.P., Yamada, M., Lim, S.S., *et al.* Inhibition by tectorigenin and tectoridin of prostaglandin E<sub>2</sub> production and cyclooxygenase-2 induction in rat peritoneal macrophages. *Biochim. Biophys. Acta* **1438(3)**, 399-407 (1999).
4. Zhang, R., Piao, M.J., Oh, M.C., *et al.* Protective effect of an isoflavone, tectorigenin, against oxidative stress-induced cell death via catalase activation. *J. Cancer Prev.* **21(4)**, 257-263 (2016).
5. Jung, S.H., Lee, Y.S., Lim, S.S., *et al.* Antioxidant activities of isoflavones from the rhizomes of *Belamcanda chinensis* on carbon tetrachloride-induced hepatic injury in rats. *Arch. Pharm. Res.* **27(2)**, 184-188 (2004).

#### 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, 01/17/2024

#### 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