

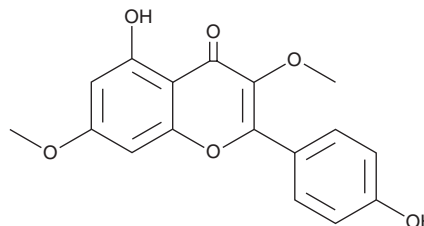
# PRODUCT INFORMATION



## Kumatakenin

Item No. 35514

**CAS Registry No.:** 3301-49-3  
**Formal Name:** 5-hydroxy-2-(4-hydroxyphenyl)-3,7-dimethoxy-4H-1-benzopyran-4-one  
**Synonyms:** Jaranol, Kaempferol 3,7-O-dimethyl ether  
**MF:** C<sub>17</sub>H<sub>14</sub>O<sub>6</sub>  
**FW:** 314.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 267, 350 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Astragalus membranaceus*



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

### Laboratory Procedures

Kumatakenin is supplied as a solid. A stock solution may be made by dissolving the kumatakenin in the solvent of choice, which should be purged with an inert gas. Kumatakenin is soluble in acetone, chloroform, dichloromethane, DMSO, and ethyl acetate.

### Description

Kumatakenin is a flavonoid that has been found in *S. aromaticum* and has diverse biological activities.<sup>1-3</sup> It inhibits the growth of SKOV3 and A2780 ovarian cancer cells in a concentration-dependent manner.<sup>1</sup> It induces apoptosis and decreases the protein levels of the tumor-associated macrophage (TAM) recruitment factors MCP-1 and RANTES in SKOV3 cells when used at a concentration of 30 μM. Kumatakenin (10 μM) inhibits neutrophil elastase activity and inhibits antigen-induced RBL-2H3 mast cell degranulation (IC<sub>50</sub> = 80.4 μM).<sup>2</sup> It is active against *S. aureus*, *E. coli*, *P. aeruginosa*, and *L. monocytogenes* (MICs = 64, 128, 128, and 512 μg/ml, respectively) and acts synergistically with ampicillin (Item No. 14417), gentamicin, or erythromycin (Item No. 16486) to reduce *L. monocytogenes* biofilm formation.<sup>3</sup>

### References

1. Woo, J.-H., Ahn, J.-H., Jang, D.S., *et al.* Effect of kumatakenin isolated from cloves on the apoptosis of cancer cells and the alternative activation of tumor-associated macrophages. *J. Agric. Food Chem.* **65**(36), 7893-7899 (2017).
2. Ebada, S.S., Al-Jawabri, N.A., Youssef, F.S., *et al.* Anti-inflammatory, antiallergic and COVID-19 protease inhibitory activities of phytochemicals from the Jordanian hawksbeard: Identification, structure-activity relationships, molecular modeling and impact on its folk medicinal uses. *RSC Adv.* **10**(62), 38128-38141 (2020).
3. Zhang, X., Wang, L., Mu, H., *et al.* Synergistic antibacterial effects of *Buddleja albiflora* metabolites with antibiotics against *Listeria monocytogenes*. *Lett. Appl. Microbiol.* **68**(1), 38-47 (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, 09/29/2022

#### 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