

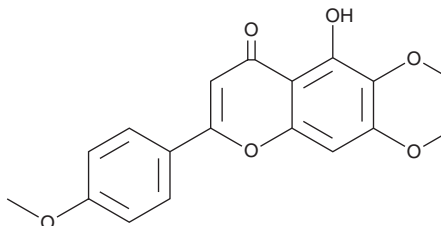
PRODUCT INFORMATION



Salvigenin

Item No. 35632

CAS Registry No.: 19103-54-9
Formal Name: 5-hydroxy-6,7-dimethoxy-2-(4-methoxyphenyl)-4H-1-benzopyran-4-one
Synonyms: 5-hydroxy-4',6,7-Trimethoxyflavone, 7-O-Methylpectolarigenin, Psathyrotin
MF: C₁₈H₁₆O₆
FW: 328.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Salvia japonica*



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

Laboratory Procedures

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

Description

Salvigenin is a polyphenol flavonoid originally found in *Alnus japonica* with diverse biological activities.¹⁻³ It inhibits hydrogen peroxide-induced apoptosis and reduces the generation of reactive oxygen species (ROS) in SH-SY5Y human neuroblastoma cells when used at a concentration of 25 μM.² Salvigenin (25 and 50 μM) decreases cleaved caspase-3 levels and the Bax/B cell lymphoma 2 (Bcl-2) ratio in SH-SY5Y cells. It reduces the viability of MCF-7 human breast cancer cells in a concentration-dependent manner.³ Salvigenin (3.65-9.68 μg/mouse per day) enhances the delayed-type hypersensitivity (DTH) response to sheep red blood cells in mice. It increases the production of IFN-γ induced by lysate antigens in isolated mouse splenocytes. Salvigenin (9.68 μg/mouse per day) reduces tumor growth in a spontaneous mouse mammary tumor (SMMT) model.

References

1. Wollenweber, E. Flavonoids from *Alnus crispa*, *A. japonica*, *A. koehnei* and *A. sinuata*. *Phytochemistry* **13**(10), 2318-2319 (1974).
2. Rafatian, G., Khodaghali, F., Farimani, M.M., et al. Increase of autophagy and attenuation of apoptosis by Salvigenin promote survival of SH-SY5Y cells following treatment with H₂O₂. *Mol. Cell. Biochem.* **371**(1-2), 9-22 (2012).
3. Noori, S., Hassan, Z.M., Yaghmaei, B., et al. Antitumor and immunomodulatory effects of salvigenin on tumor bearing mice. *Cell. Immunol.* **286**(1-2), 16-21 (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, 12/12/2022

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