

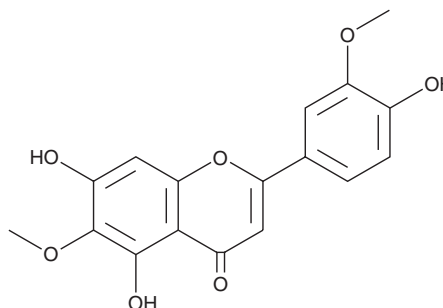
# PRODUCT INFORMATION



## Jaceosidin

Item No. 19758

CAS Registry No.: 18085-97-7  
Formal Name: 5,7-dihydroxy-2-(4-hydroxy-3-methoxyphenyl)-6-methoxy-4H-1-benzopyran-4-one  
MF: C<sub>17</sub>H<sub>14</sub>O<sub>7</sub>  
FW: 330.3  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 215, 275, 346 nm  
Supplied as: A crystalline solid  
Storage: -20°C  
Stability: ≥4 years



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

### Laboratory Procedures

Jaceosidin is supplied as a crystalline solid. A stock solution may be made by dissolving the jaceosidin in the solvent of choice, which should be purged with an inert gas. Jaceosidin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of jaceosidin in ethanol is approximately 0.2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

Jaceosidin is a natural methylated trihydroxyflavone first isolated from plants of the Compositae family. Like many flavones, jaceosidin has antioxidative actions on cells in culture.<sup>1</sup> It induces cell cycle arrest and apoptosis in cancer cells, suggesting potential roles in cancer therapy.<sup>2-4</sup> Jaceosidin also blocks cell signaling related to inflammation, including activation of NF-κB and induced expression of cyclooxygenase 1, inducible nitric oxide synthase, and matrix metalloproteinase-9.<sup>1,5</sup>

### References

1. Kim, M.-J., Han, J.-M., Jin, Y.-Y., *et al.* *In vitro* antioxidant and anti-inflammatory activities of Jaceosidin from *Artemisia princeps* Pampanini cv. Sajabal. *Arch. Pharm. Res.* **31**(4), 42-437 (2008).
2. Lee, J.-G., Kim, J.-H., Ahn, J.-H., *et al.* Jaceosidin, isolated from dietary mugwort (*Artemisia princeps*), induces G2/M cell cycle arrest by inactivating cdc25C-cdc2 via ATM-Chk1/2 activation. *Food Chem. Toxicol.* **55**, 214-221 (2013).
3. Lv, W., Sheng, X., Chen, T., *et al.* Jaceosidin induces apoptosis in human ovary cancer cells through mitochondrial pathway. *J. Biomed. Biotechnol.* 394802 (2008).
4. Woo, S.M., and Kwon, T.K. Jaceosidin induces apoptosis through Bax activation and down-regulation of Mcl-1 and c-FLIP expression in human renal carcinoma Caki cells. *Chem. Biol. Interact.* **260**, 168-175 (2016).
5. Jeong, M.A., Lee, K.W., Yoon, D.-Y., *et al.* Jaceosidin, a pharmacologically active flavone derived from *Artemisia argyi*, inhibits phorbol-ester-induced upregulation of COX-2 and MMP-9 by blocking phosphorylation of ERK-1 and -2 in cultured human mammary epithelial cells. *Ann. N.Y. Acad. Sci.* **1095**, 458-466 (2007).

#### 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, 12/22/2022

#### 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