

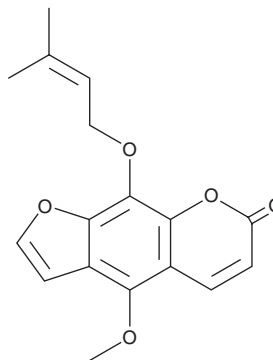
# PRODUCT INFORMATION



## Phellopterin

Item No. 38965

**CAS Registry No.:** 2543-94-4  
**Formal Name:** 7-butyl-5-hydroxy-2-methyl-2-(4-methyl-3-penten-1-yl)-2H-1-benzopyran-6-carboxylic acid  
**Synonym:** NSC 152469  
**MF:** C<sub>17</sub>H<sub>16</sub>O<sub>5</sub>  
**FW:** 300.3  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Saposhnikovia divaricata*



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

### Laboratory Procedures

Phellopterin is supplied as a solid. A stock solution may be made by dissolving the phellopterin in the solvent of choice, which should be purged with an inert gas. Phellopterin is soluble in chloroform and methanol.

### Description

Phellopterin is a furanocoumarin that has been found in *A. dahurica* and has diverse biological activities.<sup>1-3</sup> It is an antagonist of GABA<sub>A</sub> receptors (IC<sub>50</sub> = 360 nM for the rat receptor).<sup>1</sup> Phellopterin (20 μM) decreases LPS-induced secretion of prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) from primary rat peritoneal macrophages.<sup>2</sup> Topical application of phellopterin (1.2 μg/g) accelerates wound healing in a mouse model of diabetes induced by streptozotocin (Item No. 13104).<sup>3</sup> It increases the levels of mRNA encoding sirtuin 1 (Sirt1) and decreases mRNA encoding intracellular adhesion molecule-1 (Icam-1) in epithelium isolated from mice in the same model.

### References

1. Bergendorff, O., Dekermendjian, K., Nielsen, M., *et al.* Furanocoumarins with affinity to brain benzodiazepine receptors *in vitro*. *Phytochemistry* **44**(6), 1121-1124 (1997).
2. Ban, H.S., Lim, S.S., Suzuki, K., *et al.* Inhibitory effects of furanocoumarins isolated from the roots of *Angelica dahurica* on prostaglandin E<sub>2</sub> production. *Planta Med.* **69**(5), 408-412 (2003).
3. Zou, J., Duan, Y., Wang, Y., *et al.* Phellopterin cream exerts an anti-inflammatory effect that facilitates diabetes-associated cutaneous wound healing *via* SIRT1. *Phytomedicine* **107**, 154447 (2022).

#### 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, 11/15/2023

#### 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