

# PRODUCT INFORMATION



## Chrysosplenetin

Item No. 30625

**CAS Registry No.:** 603-56-5  
**Formal Name:** 5-hydroxy-2-(4-hydroxy-3-methoxyphenyl)-3,6,7-trimethoxy-4H-1-benzopyran-4-one  
**Synonyms:** Chrysosplenol B, Chrysosptertin B, EMD 20940, Polycladin

**MF:** C<sub>19</sub>H<sub>18</sub>O<sub>8</sub>

**FW:** 374.3

**Purity:** ≥98%

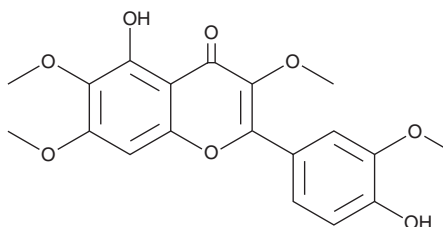
**UV/Vis.:** λ<sub>max</sub>: 257, 271, 350 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Plant/*Laggera pterodonta*



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

### Laboratory Procedures

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

### Description

Chrysosplenetin is a flavonoid that has been found in *A. annua* and has diverse biological activities.<sup>1-3</sup> It is active against *P. falciparum* *in vitro* (IC<sub>50</sub> = 23 μM).<sup>1</sup> Chrysosplenetin inhibits the cytopathic effect of enterovirus 71 (EV71) in Vero cells (EC<sub>50</sub> = 0.68 μM) and increases survival in a neonatal mouse model of EV71 infection when administered at doses of 1 and 5 mg/kg.<sup>2</sup> It also increases proliferation and osteogenic differentiation of isolated human bone marrow stromal cells (BMSCs) and prevents estrogen deficiency-induced bone loss in ovariectomized mice.<sup>3</sup>

### References

1. Liu, K.C., Yang, S.L., Roberts, M.F., *et al.* Antimalarial activity of *Artemisia annua* flavonoids from whole plants and cell cultures. *Plant Cell Rep.* **11(12)**, 637-640 (1992).
2. Dai, W., Bi, J., Li, F., *et al.* Antiviral efficacy of flavonoids against enterovirus 71 infection *in vitro* and in newborn mice. *Viruses* **11(7)**, 625 (2019).
3. Hong, G., He, X., Shen, Y., *et al.* Chrysosplenetin promotes osteoblastogenesis of bone marrow stromal cells via Wnt/β-catenin pathway and enhances osteogenesis in estrogen deficiency-induced bone loss. *Stem Cell Res. Ther.* **10(1)**, 277 (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, 10/21/20222

#### 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