

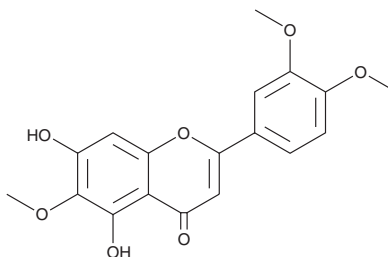
# PRODUCT INFORMATION



## Eupatilin

Item No. 19601

**CAS Registry No.:** 22368-21-4  
**Formal Name:** 2-(3,4-dimethoxyphenyl)-5,7-dihydroxy-6-methoxy-4H-1-benzopyran-4-one  
**Synonym:** NSC 122413  
**MF:** C<sub>18</sub>H<sub>16</sub>O<sub>7</sub>  
**FW:** 344.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 214, 242, 274, 337 nm  
**Supplied as:** A powder  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Eupatilin is supplied as a powder. A stock solution may be made by dissolving the eupatilin in the solvent of choice, which should be purged with an inert gas. Eupatilin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of eupatilin in these solvents is approximately 33 and 50 mg/ml, respectively.

### Description

Eupatilin is a flavonoid that has been found in *Artemisia* and has diverse biological activities.<sup>1-5</sup> It inhibits cell growth in A375 melanoma cells when used at concentrations ranging from 25 to 800 μM.<sup>1</sup> Eupatilin binds to peroxisome proliferator-activated receptor α (PPARα; IC<sub>50</sub> = 1.18 μM) and induces reporter gene expression in CV-1 cells expressing human PPARα, but not PPARγ, when used at concentrations ranging from 10 to 300 μM.<sup>2</sup> It prevents hydrogen peroxide-induced disruption of the F-actin cytoskeleton and inhibition of cell migration in AGS gastric epithelial cells.<sup>6</sup> Eupatilin also inhibits 5-lipoxygenase (5-LO) activity in cultured mastocytoma cells (IC<sub>50</sub> = 14 μM).<sup>3</sup> *In vivo*, eupatilin (5, 10, and 15 mg/kg) reduces pulmonary edema and levels of the lung injury markers surfactant protein A (SPA) and SPD in a rat model of LPS-induced acute lung injury.<sup>4</sup> It reduces serum levels of histamine and increases survival in a mouse model of allergic anaphylactic shock induced by compound 48/80 (Item No. 22173).<sup>5</sup>

### References

1. Shawi, A.A., Rasul, A., Khan, M., et al. *Afr. J. Pharm. and Pharmacol.* **5(5)**, 582-588 (2011).
2. Choi, Y., Jung, Y. and Kim, S.N. *Molecules* **20(8)**, 13753-13763 (2015).
3. Koshihara, Y., Neichi, T., Murota, S.i., et al. *FEBS Letters* **158(1)**, 41-44 (1983).
4. Liu, H., Hao, J., Wu, C., et al. *Med. Sci Monit.* **25**, 8289-8296 (2019).
5. Song, E.-H., Chung, K.-S., Kang, Y.-M., et al. *Phytomedicine* **42**, 1-8 (2018).
6. Choi, E.J., Oh, H.M., Na, B.R., et al. *Pharmaceutical Research* **25(6)**, 1355-1364 (2008).

#### 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, 11/14/2022

#### 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