

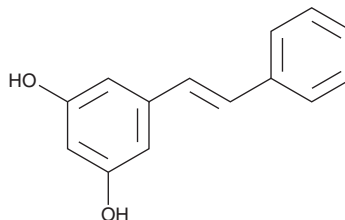
# PRODUCT INFORMATION



## Pinosylvin

Item No. 27612

**CAS Registry No.:** 22139-77-1  
**Formal Name:** 5-[(1E)-2-phenylethenyl]-1,3-benzenediol  
**Synonym:** *trans*-3,5-Dihydroxystilbene  
**MF:** C<sub>14</sub>H<sub>12</sub>O<sub>2</sub>  
**FW:** 212.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 211, 301, 309 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

Pinosylvin is supplied as a solid. A stock solution may be made by dissolving the pinosylvin in the solvent of choice, which should be purged with an inert gas. Pinosylvin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pinosylvin in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Pinosylvin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pinosylvin should first be dissolved in ethanol or DMF and then diluted with the aqueous buffer of choice. Pinosylvin has a solubility of approximately 0.01 mg/ml in a 1:50 solution of ethanol:PBS (pH 7.2) or DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pinosylvin is a stilbene originally isolated from pine heartwood that has diverse biological activities.<sup>1-6</sup> It activates sirtuin 1 (SIRT1) and induces glucose uptake in isolated rat L6 skeletal muscle myotubes.<sup>2</sup> Pinosylvin reduces radial growth in a panel of 28 plant pathogenic fungi when used at a concentration of 100 µg/ml.<sup>3</sup> It reduces expression of matrix metalloproteinase-2 (MMP-2), MMP-9, and membrane type 1-MMP in and inhibits migration of HT-1080 cells.<sup>4</sup> Pinosylvin (10 mg/kg, i.p.) reduces the number of tumor nodules and lung tumor weight in a CT26 mouse xenograft model of metastatic colon cancer. It decreases hind paw volume and myeloperoxidase (MPO) activity in a rat model of adjuvant-induced arthritis.<sup>5</sup> Pinosylvin also exhibits plant antifeedant activity against *L. americanus* (snowshoe hares).<sup>6</sup>

### References

1. Alvarez-Novoa, J.C., Erdtman, H., and Lindstedt, G. *Acta Chemica Scandinavica* **4(3)**, 444-447 (1950).
2. Modi, S., Yaluri, N., Kokkola, T., et al. *Sci. Rep.* **7(1)**, 17606 (2017).
3. Seppänen, S.K., Syrjälä, L., von Weissenberg, K., et al. *Plant Cell. Rep.* **22(8)**, 584-593 (2004).
4. Park, E.-J., Park, H.-J., Chung, H.-J., et al. *J. Nutr. Biochem.* **23(8)**, 946-952 (2012).
5. Mačičková, T., Drábiková, K., Nosál, R., et al. *Neuro. Endocrinol. Lett.* **31(Suppl 2)**, 91-95 (2013).
6. Sullivan, T.P., Crump, D.R., Wieser, H., et al. *J. Chem. Ecol.* **18(7)**, 1151-1164 (1992).

#### 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, 12/13/2022

#### 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