

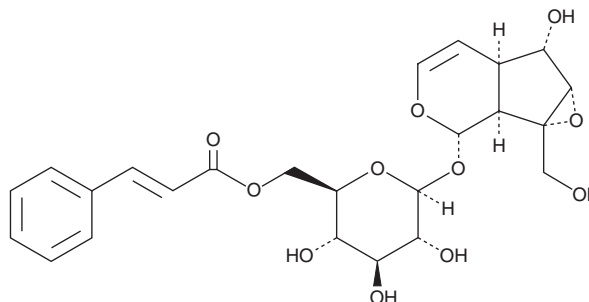
# PRODUCT INFORMATION



## Picoside I

Item No. 15431

**CAS Registry No.:** 27409-30-9  
**Formal Name:** 6-[(2E)-3-phenyl-2-propenoate]-1aS,1bS,2S,5aR,6S,6aS-hexahydro-6-hydroxy-1a-(hydroxymethyl)oxireno[4,5]cyclopenta[1,2-c]pyran-2-yl-β-D-glucopyranoside  
**Synonym:** 6'-Cinnamoylcatalpol  
**MF:** C<sub>24</sub>H<sub>28</sub>O<sub>11</sub>  
**FW:** 492.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216, 222, 276 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

Picoside I is supplied as a crystalline solid. A stock solution may be made by dissolving the picoside I in the solvent of choice, which should be purged with an inert gas. Picoside I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of picoside I in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of picoside I can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of picoside I in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Picoliv is a hepatoprotective mixture of compounds that is isolated from an herb native to the Himalayas.<sup>1</sup> Picoside I is an iridoid glycoside found in picroliv.<sup>1</sup> By itself, picoside I blocks changes in acid phosphatase activity, phospholipid levels, and lipid peroxide production induced by D-galactosamine in rat liver.<sup>2</sup> Picoside I also enhances neurite outgrowth in PC12D cells induced by basic fibroblast growth factor and 7S nerve growth factor when given at 60 μM.<sup>3,4</sup> At concentrations as low as 5 μM, picoside I enhances the ATPase activity of the efflux transporter P-glycoprotein.<sup>5</sup>

### References

1. Girish, C. and Pradhan, S.C. *Fundam. Clin. Pharmacol.* **22**, 623-632 (2008).
2. Dwivedi, Y., Rastogi, R., Garg, N.K., *et al. Pharmacol. Toxicol.* **71**, 383-387 (1992).
3. Li, P., Matsunaga, K., Yamakuni, T., *et al. Eur. J. Pharmacol.* **406(2)**, 203-208 (2000).
4. Li, P., Matsunaga, K., Yamakuni, T., *et al. Life Sci.* **71(15)**, 1821-1835 (2002).
5. Najar, I.A., Sachin, B.S., Sharma, S.C., *et al. Phytother. Res.* **24(3)**, 454-458 (2010).

#### 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, 10/11/2022

#### 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