

# PRODUCT INFORMATION

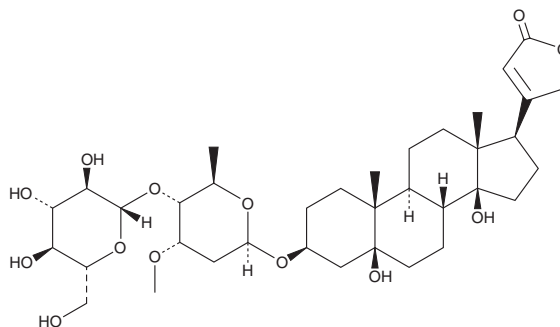


## Periplocin

Item No. 25216

**CAS Registry No.:** 13137-64-9  
**Formal Name:** (3 $\beta$ ,5 $\beta$ )-3-[(2,6-dideoxy-4-O- $\beta$ -D-glucopyranosyl-3-O-methyl- $\beta$ -D-ribohexopyranosyl)oxy]-5,14-dihydroxycard-20(22)-enolide

**MF:** C<sub>36</sub>H<sub>56</sub>O<sub>13</sub>  
**FW:** 696.8  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 219 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

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

Periplocin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, periplocin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Periplocin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Periplocin is a cardiac glycoside that has been isolated from *P. sepium* and has cardiac and anticancer activity.<sup>1</sup> It increases viability and proliferation of mouse cardiac microvascular endothelial cells (CMECs) when used at concentrations ranging from 2 to 50  $\mu$ M and improves left ventricular structure and function in a rat model of chronic heart failure.<sup>2,3</sup> Periplocin inhibits cell proliferation in nine lung cancer cell lines in a time- and dose-dependent manner with IC<sub>50</sub> values ranging from 0.12 to 53  $\mu$ M.<sup>4</sup> It induces apoptosis in SGC-7901 and MGC-803 gastric cancer cells and activates the ERK1/2-EGR1 pathway.<sup>5</sup> Periplocin (5 and 20 mg/kg) reduces tumor growth in a hepatocellular carcinoma (HCC) mouse xenograft model.<sup>1</sup> It also inhibits AKT and ERK autophosphorylation and tumor growth in an A549 lung cancer mouse xenograft model when administered at doses of 50 and 100  $\mu$ g.<sup>4</sup>

### References

1. Cheng, C.-F., Lu, I.-H., Tseng, H.-W., et al. *Evid. Based Complement. Alternat. Med.* **2013**, 958025 (2013).
2. Wang, X.-y., Gao, X.-m., Liu, H., et al. *Chin. J. Integr. Med.* **16**(1), 33-40 (2010).
3. Li, Y., Li, J., Zhou, K., et al. *Molecules* **21**(12), E1702 (2016).
4. Lu, Z.J., Zhou, Y., Song, Q., et al. *Cell Physiol. Biochem.* **26**(4-5), 609-618 (2010).
5. Li, L., Zhao, L.-M., Dai, S.-I., et al. *Cell Physiol. Biochem.* **38**(5), 1939-1951 (2016).

#### 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, 12/20/2022

#### 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](http://WWW.CAYMANCHEM.COM)