

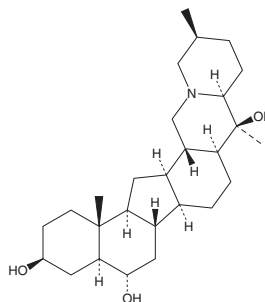
# PRODUCT INFORMATION



## Peimine

Item No. 29963

**CAS Registry No.:** 23496-41-5  
**Formal Name:** (5 $\alpha$ )-cevine-3 $\beta$ ,6 $\alpha$ ,20-triol  
**MF:** C<sub>27</sub>H<sub>45</sub>NO<sub>3</sub>  
**FW:** 431.7  
**Purity:**  $\geq$ 98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 2 years  
**Item Origin:** Plant/*Fritillaria thunbergii*



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

### Laboratory Procedures

Peimine is supplied as a crystalline solid. A stock solution may be made by dissolving the peimine in the solvent of choice, which should be purged with an inert gas. Peimine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of peimine in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Peimine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, peimine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Peimine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Peimine is an alkaloid that has been found in *Fritillaria* and has diverse biological activities, including ion channel inhibitory, anti-tussive, anticancer, and anti-inflammatory properties.<sup>1-5</sup> It is an inhibitor of voltage-gated sodium channel 1.7 (Na<sub>v</sub>1.7; IC<sub>50</sub> = 47.2  $\mu$ M) and human-ether-a-go-go (hERG), also known as K<sub>v</sub>11.1 (IC<sub>50</sub> = 43.7  $\mu$ M).<sup>2,3</sup> It increases the latent period of coughs, decreases the number of coughs, and has expectorant activity in a mouse model of ammonia liquor-induced cough when administered at a dose of 3 mg/kg.<sup>1</sup> Peimine (2.5, 5, and 10  $\mu$ M) inhibits the growth of DU145, LNCaP, and PC3 prostate cancer, but not RWPE-1 non-cancerous, cells.<sup>4</sup> It also induces phosphorylation of calcium/calmodulin-dependent protein kinase II (CaMKII) in, and apoptosis of, PC3 cells and inhibits PC3 cell invasion, migration, and epithelial-to-mesenchymal transition. Peimine (10, 20, or 50 mg/animal) reduces tumor growth in a PC3 mouse xenograft model. It decreases IL-1 $\beta$ -induced nitric oxide (NO) and prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) production and reduces inducible NO synthase (iNOS) and COX-2 protein levels in primary mouse articular chondrocytes *in vitro*.<sup>5</sup> It also decreases the severity of osteoarthritis-associated histopathological alterations in a mouse model of osteoarthritis when administered at a dose of 20 mg/kg.

### References

1. Wang, D., Zhu, J., Wang, S., et al. *Fitoterapia* **82**(8), 1290-1294 (2011).
2. Xu, J., Zhao, W., Pan, L., et al. *Fitoterapia* **111**, 1-6 (2016).
3. Kan, L., Zhao, W., Pan, L., et al. *Biomed. Pharmacother.* **89**, 838-844 (2017).
4. Tan, H., Zhang, G., Yang, X., et al. *J. Cell. Biochem.* **121**(1), 81-92 (2020).
5. Chen, K., Lv, Z.-T., Zhou, C.-H., et al. *Int. J. Mol. Med.* **43**(5), 2241-2251 (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, 05/07/2020

#### 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