

# PRODUCT INFORMATION



## Peimisine (hydrochloride)

Item No. 26438

CAS Registry No.: 900498-44-4

Formal Name: (2'R,3S,3'R,3'aS,4aS,6'S,6aR,6bS,7'aR,11aS,11bR)-1,2,3,3'a,4,4'a,5',6',6a,6b,7,7',7'a,8,11,11a,11b-octadecahydro-3-hydroxy-3',6',10,11b-tetramethyl-spiro[9H-benzo[a]fluorene-9,2'(3'H)-furo[3,2-b]pyridin]-5(6H)-one, monohydrochloride

MF:  $C_{27}H_{41}NO_3 \cdot HCl$

FW: 464.1

Purity:  $\geq 95\%$

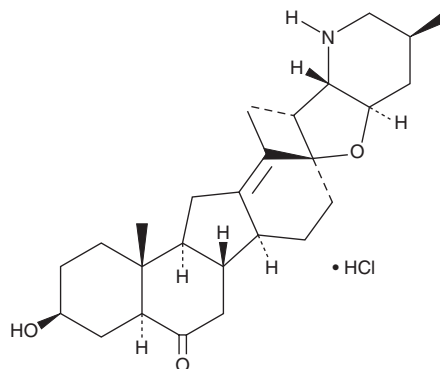
UV/Vis.:  $\lambda_{max}$ : 204 nm

Supplied as: A crystalline solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years

Item Origin: Plant/*Fritillaria hupehensis* Hsiao et K.C. Hsia



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

### Laboratory Procedures

Peimisine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the peimisine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Peimisine (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of peimisine (hydrochloride) in these solvents is approximately 1 and 16 mg/ml, respectively.

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 peimisine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of peimisine (hydrochloride) is slightly soluble in PBS, pH 7.2. We do not recommend storing the aqueous solution for more than one day.

### Description

Peimisine is a steroidal alkaloid originally isolated from *Fritillaria*.<sup>1</sup> It inhibits rabbit lung acetylcholinesterase (AChE) by 20.2% when used at a concentration of 200  $\mu M$ .

### Reference

1. An, J.-J., Zhou, J.-L., Li, H.-J., et al. Puqienine E: An angiotensin converting enzyme inhibitory steroidal alkaloid from *Fritillaria puqiensis*. *Fitoterapia* **81**(3), 149-152 (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, 11/28/2022

#### 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