

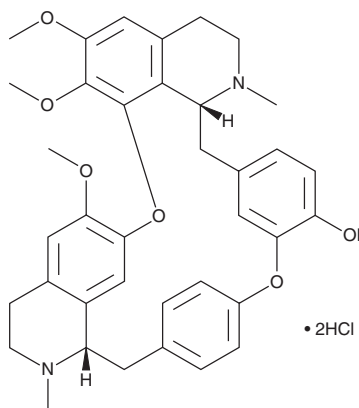
# PRODUCT INFORMATION



## Berberamine (hydrochloride)

Item No. 35593

**CAS Registry No.:** 6078-17-7  
**Formal Name:** 3,4,4a,5,16a,17,18,19-octahydro-21,22,26-trimethoxy-4,17-dimethyl-16H-1,24:6,9-dietheno-11,15-metheno-2H-pyrido[2',3':17,18][1,11]dioxacycloeicosino[2,3,4-ij]isoquinolin-12-ol, dihydrochloride  
**MF:**  $C_{37}H_{40}N_2O_6 \cdot 2HCl$   
**FW:** 681.7  
**Purity:**  $\geq 95\%$   
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years  
**Item Origin:** Plant/*Berberis julianae*



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

### Laboratory Procedures

Berberamine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the berberamine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Berberamine (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of berberamine (hydrochloride) in these solvents is approximately 12 mg/ml. Berberamine (hydrochloride) is also slightly soluble in ethanol.

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 berberamine (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of berberamine (hydrochloride) in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Berberamine is a bisbenzylisoquinoline alkaloid that has been found in *B. aristata* and has diverse biological activities.<sup>1-4</sup> It induces apoptosis and autophagy in, and is cytotoxic to, HT-29 cells when used at concentrations ranging from 3 to 100  $\mu M$ .<sup>1</sup> Berberamine reduces viral infectivity in Vero E6 cells infected with severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2;  $IC_{50} = 2.35 \mu M$ ) and decreases cell surface levels of angiotensin-converting enzyme 2 (ACE2) in Huh7 cells when used at a concentration of 10  $\mu M$ .<sup>2</sup> *In vivo*, berberamine (25 and 50 mg/kg) reduces hepatic lipid accumulation and inflammatory cytokine levels, serum and hepatic triglyceride and total cholesterol levels, and serum aspartate aminotransferase (AST) and alanine transaminase (ALT) levels in a mouse model of ethanol-induced liver injury.<sup>3</sup> It also protects against myocardial apoptosis, mitochondrial dysfunction, and lipid peroxidation in a rat model of myocardial infarction induced by isoproterenol (Item No. 15592).<sup>4</sup>

### References

1. Mou, L., Liang, B., Liu, G., et al. *J. BUON* **24**(5), 1870-1875 (2019).
2. Huang, L., Yuen, T.T.-T., Ye, Z., et al. *Signal Transduct. Target. Ther.* **6**(1), 168 (2021).
3. Liu, X.-Y., Chen, G.-N., Du, G.-M., et al. *Chin. J. Nat. Med.* **18**(3), 186-195 (2020).
4. Saranya, S., Baskaran, R., Poornima, P., et al. *J. Cell. Biochem.* **120**(3), 3101-3113 (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, 12/07/2022

#### 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)