

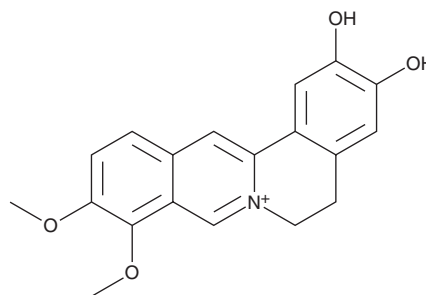
PRODUCT INFORMATION



Demethyleneberberine

Item No. 43099

CAS Registry No.: 25459-91-0
Formal Name: 5,6-dihydro-2,3-dihydroxy-9,10-dimethoxy-dibenzo[a,g]quinolizinium
Synonyms: DMB, DM-BBR, 2-Hydroxyjatrorrhizine
MF: C₁₉H₁₈NO₄
FW: 324.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Thalictrum javanicum* Blume



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

Laboratory Procedures

Demethyleneberberine is supplied as a solid. A stock solution may be made by dissolving the demethyleneberberine in the solvent of choice, which should be purged with an inert gas. Demethyleneberberine is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

Demethyleneberberine is an isoquinoline alkaloid that has been found in *P. chinensis* and has diverse biological activities.¹⁻⁵ It is also an active metabolite of berberine (Item No. 10006427).^{1,2} Demethyleneberberine is formed from berberine by cytochrome P450 (CYP) isoforms CYP2D6, CYP1A2, and CYP3A4.² It reduces viability, colony formation, and migration and invasion in A549 and H1299 lung cancer cells when used at a concentration of 40 μM.³ Demethyleneberberine (40 mg/kg per day) increases hepatic mitochondrial levels of glutathione (GSH) and glutathione peroxidase 4 (Gpx4), reduces hepatic mitochondrial levels of thiobarbituric acid reactive substances (TBARS), decreases hepatic swelling of mitochondria, and increases serum activity of alanine transaminase (ALT) in a mouse model of ethanol-induced oxidative liver damage.⁴ It decreases hepatic levels of inducible nitric oxide synthase (iNOS) and triglycerides in the same model at the same dose. Demethyleneberberine (10 mg/kg per day) reduces hepatic collagen levels and increases survival in a mouse model of thioacetamide-induced hepatic fibrosis.⁵

References

1. Cao, S., Zhou, Y., Xu, P., et al. Berberine metabolites exhibit triglyceride-lowering effects via activation of AMP-activated protein kinase in Hep G2 cells. *J. Ethnopharmacol.* **149**(2), 576-582 (2013).
2. Li, Y., Ren, G., Wang, Y.X., et al. Bioactivities of berberine metabolites after transformation through CYP450 isoenzymes. *J. Transl. Med.* **9**, 62 (2011).
3. Liu, J., Huang, X., Liu, D., et al. Demethyleneberberine induces cell cycle arrest and cellular senescence of NSCLC cells via c-Myc/HIF-1α pathway. *Phytomedicine* **91**:153678, (2021).
4. Zhang, P., Qiang, X., Zhang, M., et al. Demethyleneberberine, a natural mitochondria-targeted antioxidant, inhibits mitochondrial dysfunction, oxidative stress, and steatosis in alcoholic liver disease mouse model. *J. Pharmacol. Exp. Ther.* **352**(1), 139-147 (2015).
5. Wang, Y., Zhao, Z., Yan, Y., et al. Demethyleneberberine protects against hepatic fibrosis in mice by modulating NF-κB signaling. *Int. J. Mol. Sci.* **17**(7), 1036 (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, 04/08/2025

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