

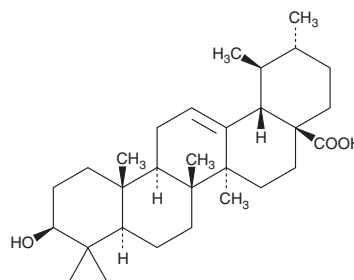
PRODUCT INFORMATION



Ursolic Acid

Item No. 10072

CAS Registry No.: 77-52-1
Formal Name: 3 β -hydroxy-urs-12-en-28-oic acid
Synonyms: Bungeolic Acid, Maerotaine, Malol, NSC 4060, NSC 167406, Prunol
MF: C₃₀H₄₈O₃
FW: 456.7
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

Ursolic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the ursolic acid in the solvent of choice. Ursolic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ursolic acid in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Ursolic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ursolic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ursolic acid has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ursolic acid is a pentacyclic triterpenoid that has been isolated from *M. pumila* and has diverse biological activities, including anticancer, hepatoprotective, anti-inflammatory, antioxidant, antimicrobial, and cardioprotective properties.¹⁻⁴ It inhibits the proliferation of HepG2 liver, MCF-7 breast, and Caco-2 colon cancer cells (EC₅₀s = 87.4, 14.4, and 34.4 μ M, respectively).² Ursolic acid reverses increases in hepatic steatosis, levels of hepatic triglycerides and free fatty acids, and hepatic TNF- α , IL-1 β , IL-6, and IL-8 mRNA expression in a dose-dependent manner in a high-fat diet-induced rat model of non-alcoholic fatty liver disease (NAFLD).⁴ Ursolic acid also scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH) radicals (IC₅₀ = 59.7 μ g/ml) and inhibits the growth of several strains of Gram-positive and Gram-negative bacteria *in vitro* including *S. aureus*, *E. coli*, *P. aeruginosa*, *K. pneumoniae*, and *S. flexneri* (MICs = 32-512 μ g/ml).³

References

1. Woźniak, Ł., Skąpska, S., and Marszałek, K. Ursolic acid--A pentacyclic triterpenoid with a wide spectrum of pharmacological activities. *Molecules* **20**(11), 20614-20641 (2015).
2. He, X. and Liu, R.H. Triterpenoids isolated from apple peels have potent antiproliferative activity and may be partially responsible for apple's anticancer activity. *J. Agric. Food Chem.* **55**(11), 4366-4370 (2007).
3. do Nascimento, P.G.G., Lemos, T.L.G., Bizerra, A.M.C., et al. Antibacterial and antioxidant activities of ursolic acid and derivatives. *Molecules* **19**(1), 1317-1327 (2014).
4. Li, S., Meng, F., Liao, X., et al. Therapeutic role of ursolic acid on ameliorating hepatic steatosis and improving metabolic disorders in high-fat diet-induced non-alcoholic fatty liver disease rats. *PLoS One* **9**(1):e86724, (2014).

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, 08/13/2018

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