

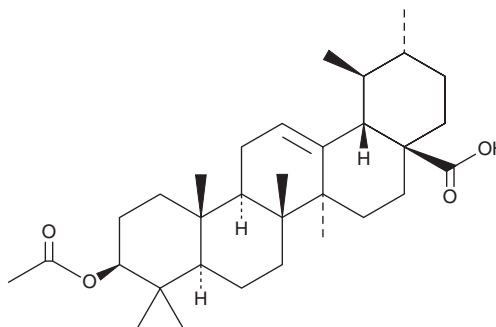
# PRODUCT INFORMATION



## 3-Acetylursolic Acid

Item No. 30220

**CAS Registry No.:** 7372-30-7  
**Formal Name:** 3 $\beta$ -(acetyloxy)-urs-12-en-28-oic acid  
**Synonym:** Ursolic Acid Acetate  
**MF:** C<sub>32</sub>H<sub>50</sub>O<sub>4</sub>  
**FW:** 498.7  
**Purity:**  $\geq$ 95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

3-Acetylursolic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-acetylursolic acid in the solvent of choice, which should be purged with an inert gas. 3-Acetylursolic acid is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 3-acetylursolic acid in these solvents is approximately 1 and 15 mg/ml, respectively.

### Description

3-Acetylursolic acid is a triterpene and derivative of ursolic acid (Item No. 10072) with diverse biological activities.<sup>1-3</sup> It inhibits ADP-, thrombin-, or epinephrine-induced aggregation of isolated rat platelets (IC<sub>50</sub>s = <1, 0.8, and <1 mg/ml, respectively).<sup>1</sup> 3-Acetylursolic acid is active against *P. falciparum* *in vitro* (IC<sub>50</sub> = 4  $\mu$ M).<sup>2</sup> It is also active against *B. cereus*, *S. aureus*, and *S. pneumoniae* (MICs = 39.1, 19.6, and 4.81  $\mu$ M, respectively).<sup>3</sup> It is cytotoxic to HepG2 cells (LC<sub>50</sub> = 351  $\mu$ M).

### References

1. Habila, J.D., Shode, F.O., and Opoku, A.R. Triterpenoids from *Eucalyptus grandis* hill ex maiden inhibits platelet aggregation. *Afr. J. Microbiol. Res.* **5(26)**, 4646-4651 (2011).
2. Innocente, A.M., Silva, G.N.S., Cruz, L.N., et al. Synthesis and antiplasmodial activity of betulinic acid and ursolic acid analogues. *Molecules* **17(10)**, 12003-12014 (2012).
3. Setzer, W.N., Setzer, M.C., Bates, R.B., et al. Biologically active triterpenoids of *Syncarpia glomulifera* bark extract from Paluma, North Queensland, Australia. *Planta Med.* **66(2)**, 176-177 (2000).

#### 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, 10/26/2022

#### 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