

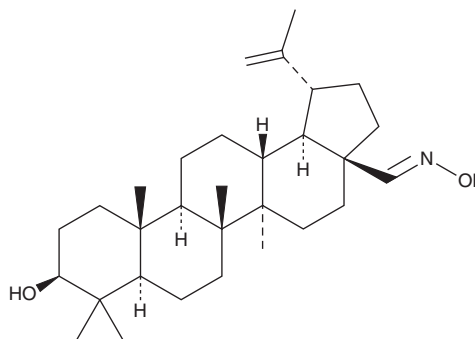
PRODUCT INFORMATION



Betulinic Aldehyde oxime

Item No. 33027

CAS Registry No.: 25613-12-1
Formal Name: 3β-hydroxy-lup-20(29)-en-28-al oxime
Synonym: Betulin 28-oxime
MF: C₃₀H₄₉NO₂
FW: 455.7
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Betulinic aldehyde oxime is supplied as a crystalline solid. A stock solution may be made by dissolving the betulinic aldehyde oxime in the solvent of choice, which should be purged with an inert gas. Betulinic aldehyde oxime is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of betulinic aldehyde oxime in these solvents is approximately 30 mg/ml.

Description

Betulinic aldehyde oxime is a pentacyclic triterpenoid and derivative of the cholesterol biosynthesis inhibitor betulin (Item No. 11041).¹ It is active against *E. aerogenes*, *E. coli*, *E. faecalis*, *P. aeruginosa*, *S. aureus*, and *C. albicans*, as well as *L. donovani* amastigotes, when used at a concentration of 50 μM.^{1,2} Betulinic aldehyde oxime (50 μM) is cytotoxic to Huh7 hepatocellular carcinoma cells.²

References

1. Alakurtti, S., Bergström, P., Sacerdoti-Sierra, N., et al. Anti-leishmanial activity of betulin derivatives. *J. Antibiot. (Tokyo)* **63(3)**, 123-126 (2010).
2. Haque, S., Nawrot, D.A., Alakurtti, S., et al. Screening and characterisation of antimicrobial properties of semisynthetic betulin derivatives. *PLoS One* **9(7)**, e102696 (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, 11/10/2022

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