

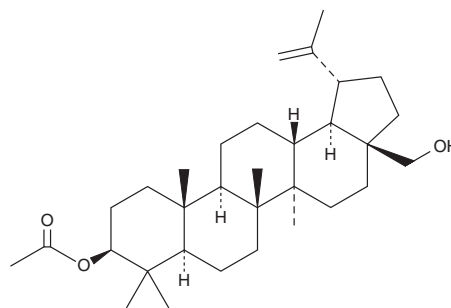
PRODUCT INFORMATION



3-Acetyl Betulin

Item No. 33049

CAS Registry No.: 27570-20-3
Formal Name: (3 β)-3-acetate-lup-20(29)-ene-3,28-diol
Synonym: 3-O-Acetyl Betulin
MF: C₃₂H₅₂O₃
FW: 484.8
Purity: $\geq 95\%$
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

3-Acetyl betulin is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-acetyl betulin in the solvent of choice, which should be purged with an inert gas. 3-Acetyl betulin is soluble in the organic solvent dimethyl formamide at a concentration of approximately 5 mg/ml.

Description

3-Acetyl betulin is a pentacyclic triterpenoid and derivative of the cholesterol biosynthesis inhibitor betulin (Item No. 11041).^{1,2} It is an intermediate in the synthesis of various lupane and germanicane triterpenes and saponins with anticancer activity.

References

1. Thibeault, D., Gauthier, C., Legault, J., *et al.* Synthesis and structure-activity relationship study of cytotoxic germanicane- and lupane-type 3 β -O-monodesmosidic saponins starting from betulin. *Bioorg. Med. Chem.* **15**(18), 6144-6157 (2007).
2. Bębenek, E., Kadela-Tomanek, M., Chrobak, E., *et al.* Novel triazoles of 3-acetylbetulin and betulone as anticancer agents. *Med. Chem. Res.* **27**(9), 2051-2061 (2018).

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, 11/03/2022

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