

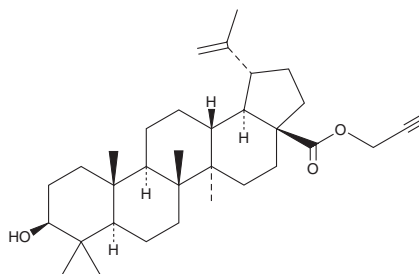
# PRODUCT INFORMATION



## Betulinic Acid propargyl ester

Item No. 33030

**CAS Registry No.:** 1400262-27-2  
**Formal Name:** (3β)-3-hydroxy-lup-20(29)-en-28-oic acid, 2-propyn-1-yl ester  
**Synonym:** Click Tag™ Betulinic Acid propargyl ester  
**MF:** C<sub>33</sub>H<sub>50</sub>O<sub>3</sub>  
**FW:** 494.8  
**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 acid propargyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the betulinic acid propargyl ester in the solvent of choice, which should be purged with an inert gas. Betulinic acid propargyl ester is soluble in the organic solvent dimethyl formamide at a concentration of approximately 30 mg/ml.

### Description

Betulinic acid propargyl ester is a pentacyclic lupane triterpene and alkyne derivative of betulinic acid (Item No. 11686).<sup>1,2</sup> It has been used in click chemistry reactions for the synthesis of cytotoxic anticancer betulinic acid derivatives.

### References

1. Thi, T.A.D., Tuyet, N.T.K., The, C.P., *et al.* Synthesis and cytotoxic evaluation of novel ester-triazole-linked triterpenoid-AZT conjugates. *Bioorg. Med. Chem. Lett.* **24**(22), 5190-5194 (2014).
2. Bębenek, E., Chrobaka, E., Wietrzyk, J., *et al.* Synthesis, structure and cytotoxic activity of acetylenic derivatives of betulonic and betulinic acids. *J. Mol. Struct.* **1106**, 210-216 (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, 08/24/2023

#### 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