

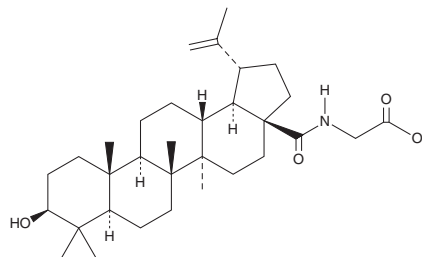
PRODUCT INFORMATION



Betulinic glycine amide

Item No. 33031

CAS Registry No.: 174740-40-0
Formal Name: N-[(3β)-3-hydroxy-28-oxolup-20(29)-en-28-yl]-glycine
MF: C₃₂H₅₁NO₄
FW: 513.8
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 glycine amide is supplied as a crystalline solid. A stock solution may be made by dissolving the betulinic glycine amide in the solvent of choice, which should be purged with an inert gas. Betulinic glycine amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of betulinic glycine amide in these solvents is approximately 15, 10, and 30 mg/ml, respectively.

Description

Betulinic glycine amide is a pentacyclic triterpenoid and derivative of betulinic acid (Item No. 11686).^{1,2} It has been used as a precursor in the synthesis of betulinic acid derivatives with anticancer and antiviral activity.

References

1. Soler, F., Poujade, C., Evers, M., *et al.* Betulinic acid derivatives: A new class of specific inhibitors of human immunodeficiency virus type 1 entry. *J. Med. Chem.* **39(5)**, 1069-1083 (1996).
2. Jeong, H.J., Chai, H.B., Park, S.Y., *et al.* Preparation of amino acid conjugates of betulinic acid with activity against human melanoma. *Bioorg. Med. Chem. Lett.* **9(8)**, 1201-1204 (1999).

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