

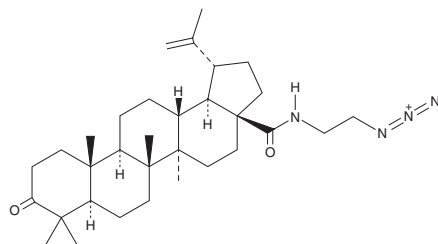
PRODUCT INFORMATION



N-(2-Azidoethyl)betulonamide

Item No. 33061

CAS Registry No.: 2055270-64-7
Formal Name: N-(2-azidoethyl)-3-oxo-lup-20(29)-en-28-amide
Synonym: Click Tag™ N-(2-Azidoethyl)betulonamide
MF: C₃₂H₅₀N₄O₂
FW: 522.8
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

N-(2-Azidoethyl)betulonamide is supplied as a crystalline solid. A stock solution may be made by dissolving the N-(2-azidoethyl)betulonamide in the solvent of choice, which should be purged with an inert gas. N-(2-Azidoethyl)betulonamide is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

N-(2-Azidoethyl)betulonamide is a pentacyclic triterpenoid, a derivative of betulonic acid (Item No. 29765), and an intermediate in the synthesis of betulonic acid derivatives with *in vitro* cancer cell cytotoxicity.¹

Reference

1. Suman, P., Patel, A., Solano, L., *et al.* Synthesis and cytotoxicity of Baylis-Hillman template derived betulinic acid-triazole conjugates. *Tetrahedron* **73**(29), 4214-4226 (2017).

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