

PRODUCT INFORMATION



28-(Poc-amino)betulin

Item No. 33069

Formal Name: prop-2-yn-1-yl (((1R,3aS,5aR,5bR,7aR,9S,11aR,11bR,13aR,13bR)-9-hydroxy-5a,5b,8,8,11a-pentamethyl-1-(prop-1-en-2-yl)icosahydro-3aH-cyclopenta[a]chrysen-3a-yl)methyl)carbamate

Synonym: Click Tag™ 28-(Poc-amino)betulin

MF: C₃₄H₅₃NO₃

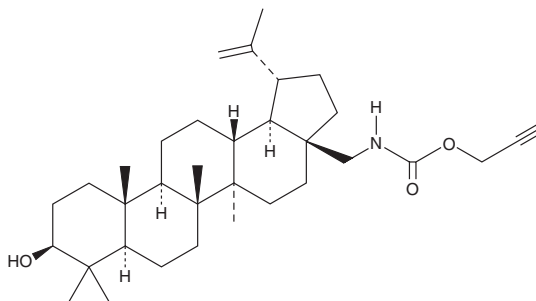
FW: 523.8

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

28-(Poc-amino)betulin is supplied as a crystalline solid. A stock solution may be made by dissolving the 28-(Poc-amino)betulin in the solvent of choice, which should be purged with an inert gas. 28-(Poc-amino)betulin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 28-(Poc-amino)betulin in these solvents is approximately 1 mg/ml in ethanol and approximately 30 mg/ml in DMSO and DMF.

28-(Poc-amino)betulin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 28-(Poc-amino)betulin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 28-(Poc-amino)betulin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

28-(Poc-amino)betulin is a terpene and an alkyne derivative of the cholesterol biosynthesis inhibitor betulin (Item No. 11041). 28-(Poc-amino)betulin can be used for click chemistry applications.

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, 08/25/2023

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