

PRODUCT INFORMATION



28-Acetalbetulin

Item No. 33053

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

MF: $C_{33}H_{54}O_3$

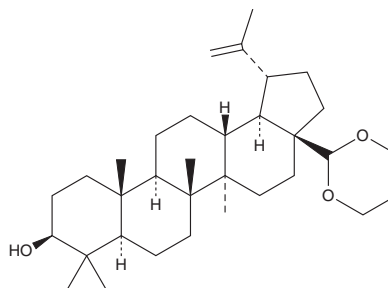
FW: 498.8

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

28-Acetalbetulin is supplied as a crystalline solid. A stock solution may be made by dissolving the 28-acetalbetulin in the solvent of choice, which should be purged with an inert gas. 28-Acetalbetulin is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml. 28-Acetalbetulin is also slightly soluble in DMSO.

Description

28-Acetalbetulin is a pentacyclic triterpenoid and derivative of the cholesterol biosynthesis inhibitor betulin (Item No. 11041).

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, 11/11/2022

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