

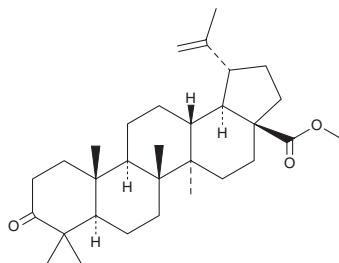
PRODUCT INFORMATION



Betulonic Acid methyl ester

Item No. 33054

CAS Registry No.: 4356-31-4
Formal Name: 3-oxo-lup-20(29)-en-28-oic acid, methyl ester
Synonyms: Betulonic Acid 28-O-methyl ester, Methyl Betulonate, NSC 152535
MF: C₃₁H₄₈O₃
FW: 468.7
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

Betulonic acid methyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the betulonic acid methyl ester in the solvent of choice, which should be purged with an inert gas. Betulonic acid methyl ester is soluble in the organic solvent dimethyl formamide at a concentration of approximately 5 mg/ml.

Description

Betulonic acid methyl ester is a pentacyclic lupane triterpenoid and a methyl ester form of betulonic acid (Item No. 29765).

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