

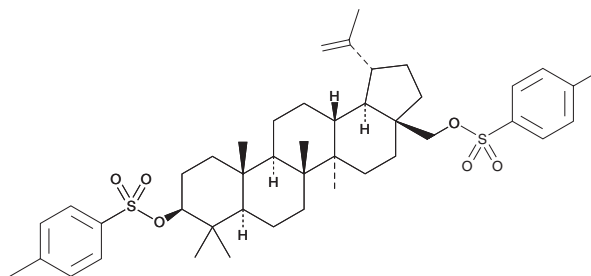
PRODUCT INFORMATION



Betulin ditosylate

Item No. 33029

CAS Registry No.: 77265-02-2
Formal Name: (3 β)-lup-20(29)-ene-3,28-diol, 3,28-bis(4-methylbenzenesulfonate)
Synonym: Betulin 3,28-di-O-tosylate
MF: C₄₄H₆₂O₆S₂
FW: 751.1
Purity: \geq 95%
UV/Vis.: λ_{max} : 225 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Betulin ditosylate is supplied as a crystalline solid. A stock solution may be made by dissolving the betulin ditosylate in the solvent of choice, which should be purged with an inert gas. Betulin ditosylate is soluble in the organic solvent chloroform at a concentration of approximately 1 mg/ml. Betulin ditosylate is also slightly soluble in dimethyl formamide.

Description

Betulin ditosylate is a derivative of betulin (Item No. 11041).¹

Reference

1. Peipiņš, U., Freimanis, N., Stepanovs, D., *et al.* Betulin 3,28-di-O-tosyl-ate. *Acta Crystallogr. Sect. E Struct. Rep. Online* **70(Pt 8)**, o879-o880 (2014).

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