

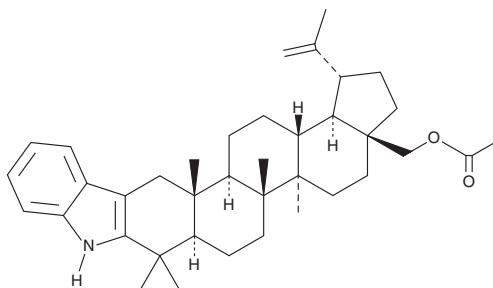
PRODUCT INFORMATION



28-acetyl-2,3-Indolobetulin

Item No. 33063

CAS Registry No.: 56855-69-7
Formal Name: 1'H-lupa-2,20(29)-dieno[3,2-b]indol-28-ol, acetate ester
MF: C₃₈H₅₃NO₂
FW: 555.8
Purity: ≥95%
UV/Vis.: λ_{max}: 229 nm
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-acetyl-2,3-Indolobetulin is supplied as a crystalline solid. A stock solution may be made by dissolving the 28-acetyl-2,3-indolobetulin in the solvent of choice, which should be purged with an inert gas. 28-acetyl-2,3-Indolobetulin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 28-acetyl-2,3-indolobetulin in ethanol and DMSO is approximately 1 mg/ml and approximately 10 mg/ml in DMF.

Description

28-acetyl-2,3-Indolobetulin is a 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