

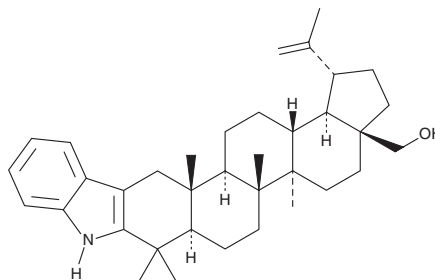
# PRODUCT INFORMATION



## 2,3-Indolobetulin

Item No. 33062

**CAS Registry No.:** 2365493-59-8  
**Formal Name:** 1'H-lupa-2,20(29)-dieno[3,2-b]indol-28-ol  
**MF:** C<sub>36</sub>H<sub>51</sub>NO  
**FW:** 513.8  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 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

2,3-Indolobetulin is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,3-indolobetulin in the solvent of choice, which should be purged with an inert gas. 2,3-Indolobetulin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2,3-indolobetulin in these solvents is approximately 10 and 25 mg/ml, respectively.

### Description

2,3-Indolobetulin is a pentacyclic triterpenoid and derivative of betulin (Item No. 11041).<sup>1</sup> It has been used in the synthesis of α-glucosidase inhibitors.

### Reference

1. Khusnutdinova, E.F., Smirnova, I.E., Kazakova, O.B., *et al.* Synthesis and evaluation of 2,3-indolotriterpenoids as new α-glucosidase inhibitors. *Med. Chem. Res.* **26**, 2732-2742 (2017).

**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