

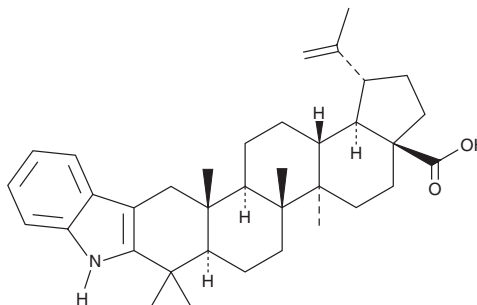
PRODUCT INFORMATION



2,3-Indolobetulonic Acid

Item No. 33065

CAS Registry No.: 905837-93-6
Formal Name: 1'H-lupa-2,20(29)-dieno[3,2-b]
indol-28-oic acid
MF: $C_{36}H_{49}NO_2$
FW: 527.8
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 228 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

2,3-Indolobetulonic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,3-indolobetulonic acid in the solvent of choice, which should be purged with an inert gas. 2,3-Indolobetulonic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2,3-indolobetulonic acid in ethanol and DMSO is approximately 5 mg/ml and approximately 30 mg/ml in DMF.

Description

2,3-Indolobetulonic acid is pentacyclic triterpenoid derivative of betulonic acid (Item No. 29765) and an inhibitor of α -glucosidase ($IC_{50} = 1.8 \mu M$).¹

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