

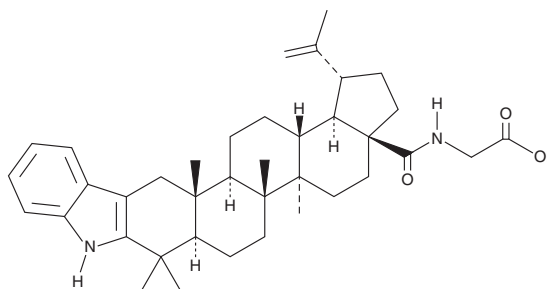
PRODUCT INFORMATION



2,3-Indolobetulinic glycine amide

Item No. 33064

CAS Registry No.: 905838-14-4
Formal Name: N-(28-oxo-1'H-lupa-2,20(29)-dieno[3,2-b]indol-28-yl)-glycine
MF: $C_{38}H_{52}N_2O_3$
FW: 584.8
Purity: $\geq 95\%$
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

2,3-Indolobetulinic glycine amide is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,3-indolobetulinic glycine amide in the solvent of choice, which should be purged with an inert gas. 2,3-Indolobetulinic glycine amide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2,3-indolobetulinic glycine amide in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

Description

2,3-Indolobetulinic glycine amide is a pentacyclic triterpene and derivative of betulinic acid (Item No. 11686).¹ It has been used in the synthesis of betulinic acid derivatives with anticancer activity.

Reference

1. Kumar, V., Rani, N., Aggarwal, P., *et al.* Synthesis and cytotoxic activity of heterocyclic ring-substituted betulinic acid derivatives. *Bioorg. Med. Chem. Lett.* **18(18)**, 5058-5062 (2008).

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