

PRODUCT INFORMATION



Blinin

Item No. 30945

CAS Registry No.: 125675-09-4
Formal Name: 4-[2-[(1S,2R,4aS,7S,8aR)-4a-[(acetyloxy)methyl]-1,2,3,4,4a,7,8,8a-octahydro-7-hydroxy-5-(hydroxymethyl)-1,2-dimethyl-1-naphthalenyl]ethyl]-2(5H)-furanone

MF: C₂₂H₃₂O₆

FW: 392.5

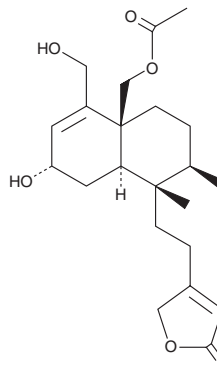
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Conyza blinii*



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

Laboratory Procedures

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

Blinin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, blinin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Blinin has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Blinin is a diterpene originally isolated from *C. blinii* with antiulcerogenic activity.^{1,2} *In vivo*, blinin decreases gastric mucosal levels of malondialdehyde (MDA) and ulcer area in a rat model of pyloric ligation-induced ulcers.²

References

1. Dogan, M., Erol, M., Cesur, Y., *et al.* The effect of 25-hydroxyvitamin D₃ on the immune system. *J. Pediatr. Endocrinol. Metab.* **22**(10), 929-935 (2009).
2. Su, Y., Chen, L., Luo, Y., *et al.* Chemical constituents and their antiulcerogenic studies on whole herb of *Conyza blinii*. *Zhongcaoyao* **38**(3), 332-334 (2007).

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, 12/09/2022

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