

PRODUCT INFORMATION



Conagenin

Item No. 40222

CAS Registry No.: 134381-30-9
Formal Name: N-(3,5-dideoxy-3-methyl-D-xylonoyl)-2-methyl-L-serine

Synonym: CNG, (+)-Conagenin

MF: $C_{10}H_{19}NO_6$

FW: 249.3

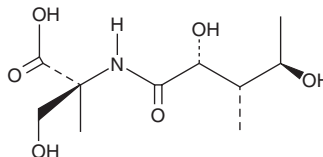
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Bacterium/*Streptomyces roseosporus* MI696-AF3



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

Laboratory Procedures

Conagenin (CNG) is supplied as a solid. A stock solution may be made by dissolving the CNG in the solvent of choice, which should be purged with an inert gas. CNG is soluble in organic solvents such as methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

CNG is a microbial metabolite originally isolated from *S. roseosporus* that has immunomodulatory activity.¹ It increases the proliferation of isolated rat splenic T cells induced by concanavalin A (Item No. 14951) when used at concentrations of 6.25 and 25 $\mu\text{g/ml}$. CNG (0.5 and 5 mg/kg) reduces tumor volume and increases the number of activated effector T cells in an IMC carcinoma mouse xenograft model.² It also increases the phagocytic activity of isolated rat alveolar macrophages when administered at a dose of 10 mg/kg.³

References

1. Yamashita, T., Iijima, M., Nakamura, H., *et al.* Conagenin, a low molecular weight immunomodulator produced by *Streptomyces roseosporus*. *J. Antibiot. (Tokyo)* **44**(5), 557-559 (1991).
2. Kawatsu, M., Yamashita, T., Osono, M., *et al.* Effect of conagenin in tumor bearing mice. Antitumor activity, generation of effector cells and cytokine production. *J. Antibiot. (Tokyo)* **46**(11), 1692-1698 (1993).
3. Hamada, M., Yamamoto, S., Moriguchi, S., *et al.* Phagocytosis of alveolar macrophages after conagenin injection to rats. *J. Antibiot. (Tokyo)* **54**(4), 349-353 (2001).

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, 07/25/2025

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