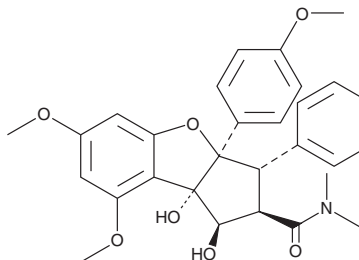


PRODUCT INFORMATION



Rocaglamide Item No. 14841

CAS Registry No.: 84573-16-0
Formal Name: (1R,2R,3S,3aR,8bS)-2,3,3a,8b-tetrahydro-1,8b-dihydroxy-6,8-dimethoxy-3a-(4-methoxyphenyl)-N,N-dimethyl-3-phenyl-1H-cyclopenta[b]benzofuran-2-carboxamide
Synonyms: NSC 326408, Roc-A, Rocaglamide A
MF: C₂₉H₃₁NO₇
FW: 505.6
Purity: ≥95%
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

Rocaglamide is supplied as a crystalline solid. A stock solution may be made by dissolving the rocaglamide in the solvent of choice, which should be purged with an inert gas. Rocaglamide is soluble in organic solvents such as ethanol and DMSO. The solubility of rocaglamide in these solvents is approximately 5 and 10 mg/ml, respectively.

Rocaglamide is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Rocaglamide is an anti-inflammatory, insecticidal, and anticancer tetrahydrobenzofuran isolated from *Aglaia* species. It has been shown to inhibit both TNF- α and the activation of NF- κ B in Jurkat T cells with IC₅₀ values in the nanomolar range.¹ At 25 nM, rocaglamide induces apoptosis in various human leukemia cell lines, activating p38 MAPK/JNK and suppressing ERK.² Rocaglamide also reduces IFN- γ , TNF- α , IL-2, and IL-4 production in peripheral blood T cells at a concentration of 50 nM.³ Furthermore, rocaglamide can inhibit the T cell expression of the immune response transcription factor, nuclear factor of activated T cells.³

References

1. Baumann, B., Bohnenstengel, F., Siegmund, D., *et al.* Rocaglamide derivatives are potent inhibitors of NF- κ B activation in T-cells. *J. Biol. Chem.* **277**(47), 44791-44800 (2002).
2. Zhu, J.Y., Lavrik, I.N., Mahlknecht, U., *et al.* The traditional Chinese herbal compound rocaglamide preferentially induces apoptosis in leukemia cells by modulation of mitogen-activated protein kinase activities. *Int. J. Cancer* **121**(8), 1839-1846 (2007).
3. Proksch, P., Giaisi, M., Treiber, M.K., *et al.* Rocaglamide derivatives are immunosuppressive phytochemicals that target NF-AT activity in T cells. *J. Immunol.* **174**(11), 7075-7084 (2005).

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, 12/12/2022

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