

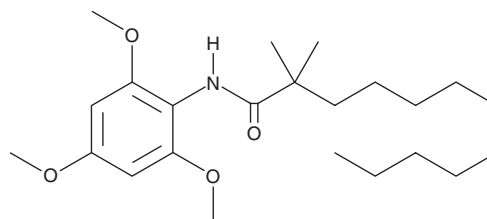
PRODUCT INFORMATION



CI-976

Item No. 17235

CAS Registry No.: 114289-47-3
Formal Name: 2,2-dimethyl-N-(2,4,6-trimethoxyphenyl)-dodecanamide
Synonym: PD 128042
MF: C₂₃H₃₉NO₄
FW: 393.6
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

CI-976 is supplied as a crystalline solid. A stock solution may be made by dissolving the CI-976 in the solvent of choice, which should be purged with an inert gas. CI-976 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CI-976 in these solvents is approximately 30 mg/ml.

CI-976 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CI-976 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CI-976 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Acyl-coenzyme A:cholesterol acyltransferase 1 (ACAT-1) is a sterol O-acyltransferase that catalyzes the formation of fatty acid-cholesterol esters, an important step in lipoprotein assembly and dietary cholesterol absorption.^{1,2} CI-976 is a potent, selective inhibitor of ACAT-1 (IC₅₀ = 73 nM).^{3,4} It is orally bioavailable and decreases plasma total cholesterol, very low density lipoprotein (VLDL) cholesterol, LDL cholesterol, apolipoprotein B, liver cholesteryl esters, and VLDL and LDL cholesteryl ester content in rabbits given a diet that induces hypercholesterolemia.⁵ CI-976 also diminishes atherosclerotic activity in hypercholesterolemic rabbits.⁶

References

1. Buhman, K.K., Chen, H.C., and Farese, R.V., Jr. The enzymes of neutral lipid synthesis. *J. Biol. Chem.* **276(44)**, 40369-40372 (2001).
2. Shi, Y. and Burn, P. Lipid metabolic enzymes: Emerging drug targets for the treatment of obesity. *Nat. Rev. Drug Discov.* **3(8)**, 695-710 (2004).
3. Field, F.J., Albright, E., and Mathur, S. Inhibition of acylcoenzyme A: cholesterol acyltransferase activity by PD128042: Effect on cholesterol metabolism and secretion in CaCo-2 cells. *Lipids* **26(1)**, 1-8 (1991).
4. O'Brien, P.M., Sliskovic, D.R., Blankley, C.J., et al. Inhibitors of acyl-CoA:cholesterol O-acyl transferase (ACAT) as hypocholesterolemic agents. 8. Incorporation of amide or amine functionalities into a series of disubstituted ureas and carbamates. Effects on ACAT inhibition in vitro and efficacy in vivo. *J. Med. Chem.* **37(12)**, 1810-1822 (1994).
5. Krause, B.R., Pape, M.E., Kieft, K., et al. ACAT inhibition decreases LDL cholesterol in rabbits fed a cholesterol-free diet. *Arterioscler. Thromb.* **14(4)**, 598-604 (1994).
6. Bocan, T.M.A., Mueller, S.B., Uhlendorf, P.D., et al. Comparison of CI-976, an ACAT inhibitor, and selected lipid-lowering agents for antiatherosclerotic activity in iliac-femoral and thoracic aortic lesions. *Arterioscler. Thromb.* **11(6)**, 1830-1843 (1991).

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, 11/04/2022

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