

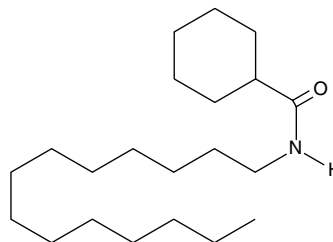
Product Information



N-Cyclohexanecarbonyltetradecylamine

Item No. 10008317

CAS Registry No.: 1215071-05-8
Formal Name: N-tetradecyl-cyclohexanecarboxamide
MF: C₂₁H₄₁NO
FW: 323.6
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that N-cyclohexanecarbonyltetradecylamine be stored as supplied at -20°C. It will be stable for at least one year.

N-Cyclohexanecarbonyltetradecylamine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-cyclohexanecarbonyltetradecylamine in an organic solvent purged with an inert gas. N-cyclohexanecarbonyltetradecylamine is soluble in ethanol. The solubility of N-cyclohexanecarbonyltetradecylamine in this solvent is approximately 2 mg/ml.

N-cyclohexanecarbonyltetradecylamine is sparingly soluble in aqueous buffers. If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that 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.

Numerous analogs of fatty acyl ethanolamides potentiate the intrinsic biological activity of endocannabinoids.¹ This potentiation is ascribed either to inhibition of AEA reuptake into neurons, or inhibition of fatty acid amide hydrolase (FAAH) within the neurons.² However, Ueda, *et al.* has recently cloned another amidase, the acidic PEAase that promotes the hydrolysis of palmitoylethanolamide.³ N-Cyclohexanecarbonyltetradecylamine is an analog of N-cyclohexanecarbonylpentadecylamine, a selective inhibitor of acidic PEAase with an IC₅₀ value of 4.5 μM, that contains one less carbon in the alkyl chain.⁴ The biological activity of N-cyclohexanecarbonyltetradecylamine has not been documented.

References

1. Khanolkar, A.D. and Makriyannis, A. Structure-activity relationships of anandamide, an endogenous cannabinoid ligand. *Life Sci.* **65**, 607-616 (1999).
2. Deutsch, D.G., Glaser, S.T., Howell, J.M., *et al.* The cellular uptake of anandamide is coupled to its breakdown by fatty-acid amide hydrolase. *J. Biol. Chem.* **276**(10), 6967-6973 (2001).
3. Ueda, N., Yamanaka, K., and Yamamoto, S. Purification and characterization of an acid amidase selective for N-palmitoylethanolamine, a putative endogenous anti-inflammatory substance. *J. Biol. Chem.* **276**(38), 35552-35557 (2001).
4. Tsuboi, K., Hilligsmann, C., Vandevoorde, S., *et al.* N-cyclohexanecarbonylpentadecylamine: A selective inhibitor of the acid amidase hydrolysing N-acyl ethanolamines, as a tool to distinguish acid amidase from fatty acid amide hydrolase. *Biochem J.* **379**, 99-106 (2004).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008317

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/12/2013