

PRODUCT INFORMATION



Lauric Acid Leelamide

Item No. 10008618

Formal Name: 1R,2,3,4,4aS,9,10,10aR-octahydro-1,4a-dimethyl-7-(1-methylethyl)-1-phenanthrene-dodecanamide

MF: C₃₂H₅₃NO

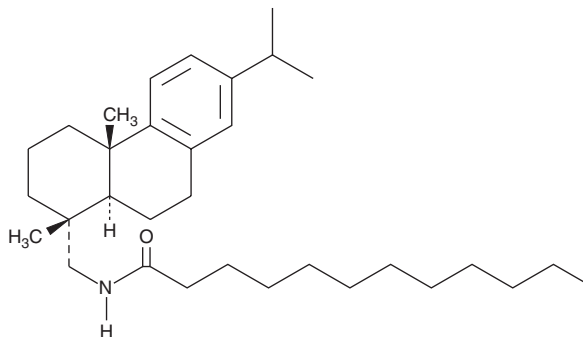
FW: 467.8

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥1 year



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

Laboratory Procedures

Lauric acid leelamide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of lauric acid leelamide in these solvents is approximately 20 mg/ml.

Description

Lauric acid leelamide is the lauric (C-12) amide analog of leelamine. Leelamine has weak affinity for the human central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) receptors, exhibiting 20% displacement of [³H]-CP55940 at a concentration of 10 μM.¹ Leelamine inhibits pyruvate dehydrogenase kinase (PDHK) with an IC₅₀ value of 9.5 μM.² Derivatives of leelamine exhibit anti-inflammatory activity and show moderate inhibition of phospholipase A₂ activity from a variety of sources.³ There are no published studies of the pharmacological properties of lauric acid leelamide.

References

1. Martin, B.R. Leelamine. Personal Communication (2006).
2. Aicher, T.D., Damon, R.E., Koletar, J., *et al.* Triterpene and diterpene inhibitors of pyruvate dehydrogenase kinase (PDK). *Bioorg. Med. Chem. Lett.* **9(15)**, 2223-2228 (1999).
3. Wilkerson, W., DeLucca, I., Galbraith, W., *et al.* Antiinflammatory phospholipase-A2 inhibitors. I. *Eur. J. Med. Chem.* **26**, 667-676 (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, 06/10/2021

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