

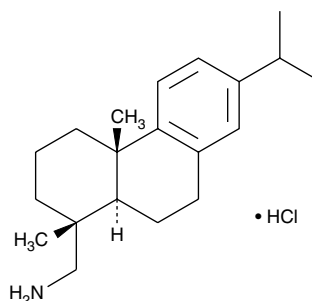
Product Information



Leelamine (hydrochloride)

Item No. 10008614

CAS Registry No.: 16496-99-4
Formal Name: 1R,2,3,4,4aS,9,10,10aR-octahydro-1,4a-dimethyl-7-(1-methylethyl)-1-phenanthrenemethanamine, monohydrochloride
Synonym: Dehydroabietylamine
MF: C₂₀H₃₁N • HCl
FW: 321.9
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that leelamine (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

Leelamine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the leelamine (hydrochloride) in an organic solvent purged with an inert gas. Leelamine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of leelamine (hydrochloride) in these solvents is approximately 20 mg/ml.

Leelamine (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, leelamine (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Leelamine (hydrochloride) has a solubility of approximately 0.15 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Leelamine is a diterpene molecule whose name derives from the Sanskrit word *leela* which means “play”. It 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 (PDK) 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

References

1. Martin, B.R. Personal Communication.
2. Aicher, T.D., Damon, R.E., Koletar, J., *et al.* Triterpene and diterpene inhibitors of pyruvate dehydrogenase kinase (PDK). *Bioorganic & Medicinal Chemistry Letters* **9**, 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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008614

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

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