PRODUCT INFORMATION



Leelamine (hydrochloride)

Item No. 10008614

CAS Registry No.: Formal Name:	16496-99-4 1R,2,3,4,4aS,9,10,10aR-octahydro- 1,4a-dimethyl-7-(1-methylethyl)-1- phenanthrenemethanamine, monohydrochloride	CHa
Synonym:	Dehydroabietylamine	
MF:	$C_{20}H_{31}N \bullet HCI$	ſ Ť Ť
FW:	321.9	
Purity:	≥98%	• HCI
Supplied as:	A crystalline solid	H ₃ C H
Storage:	-20°C	
Stability:	≥4 years	H ₂ N

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

Laboratory Procedures

Leelamine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the leelamine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Leelamine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of leelamine (hydrochloride) in ethanol and DMF is approximately 20 mg/ml and approximately 25 mg/ml in DMSO.

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.

Description

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 and IC₅₀ value of 9.5 μ M.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). 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(7), 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, 04/30/2024

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