

PRODUCT INFORMATION

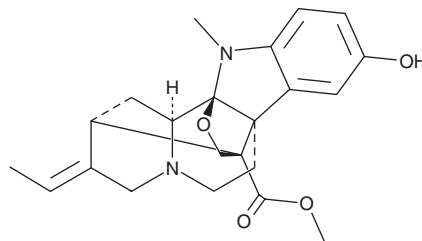


Akuammine

Item No. 35149

CAS Registry No.: 3512-87-6
Formal Name: (2S,3E,7aS,12aS,12bS,15R)-3-ethylidene-1,3,4,6,7,12b-hexahydro-9-hydroxy-12-methyl-2H,12H-12a,2,7a-(epoxyethanylylidene)indolo[2,3-a]quinolizine-15-carboxylic acid, methyl ester
Synonyms: (-)-Akuammine, Vincamajoridine

MF: C₂₂H₂₆N₂O₄
FW: 382.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Picralima nitida*



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

Laboratory Procedures

Akuammine is supplied as a crystalline solid. A stock solution may be made by dissolving the akuammine in the solvent of choice, which should be purged with an inert gas. Akuammine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of akuammine in these solvents is approximately 1 and 2 mg/ml, respectively.

Description

Akuammine is an indole alkaloid that has been found in *Picralima nitida* and has analgesic activity.¹ It selectively binds to the μ - and κ -opioid receptors over the δ -opioid receptor (K_i s = 0.3, 1.68, and 10.4 μ M for the human receptors, respectively). Akuammine inhibits forskolin-induced cAMP production in HEK293 cells expressing human μ - or κ -opioid receptors (IC_{50} s = 2.6 and 0.073 μ M, respectively). It increases the latency to withdrawal in the tail-flick or hot plate test in mice when administered at a dose of 60 mg/kg.

Reference

1. Creed, S.M., Gutridge, A.M., Argade, M.D., *et al.* Isolation and pharmacological characterization of six opioidergic *Picralima nitida* alkaloids. *J. Nat. Prod.* **84**(1), 71-80 (2021).

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, 04/22/2025

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