

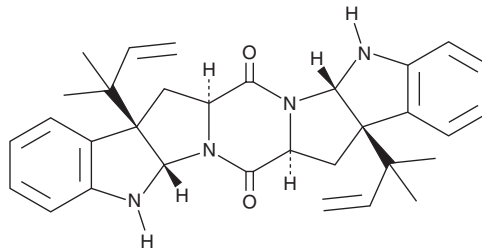
PRODUCT INFORMATION



Amauromine

Item No. 23886

CAS Registry No.: 88360-87-6
Formal Name: (5aS,7aS,8aR,13aS,15aS,16aR)-8a,16a-bis(1,1-dimethyl-2-propen-1-yl)-5a,8,8a,13,13a,15a,16,16a-octahydro-pyrazino[1'',2'':1,5;4'',5'':1',5']dipyrrolo[2,3-b:2',3'-b']diindole-7,15(5H,7aH)-dione
MF: C₃₂H₃₆N₄O₂
FW: 508.7
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Amauromine is supplied as a solid. A stock solution may be made by dissolving the amauroamine in the solvent of choice. Amauroamine is soluble in organic solvents such as ethanol, DMSO, and methanol, which should be purged with an inert gas.

Description

Amauroamine is a neutral antagonist of the cannabinoid (CB) receptor CB₁ that is selective for CB₁ (K_i = 178 nM; K_b = 66.6 nM) over CB₂, with no activity at CB₂ receptors at concentrations up to 10 μM.¹ It is also an antagonist of GPR18 (IC₅₀ = 3.74 μM).² Amauroamine has vasodilatory activity.³

References

1. Elsebai, M.F., Rempel, V., Schnakenburg, G., *et al.* Identification of a potent and selective cannabinoid CB₁ receptor antagonist from *Auxarthron reticulatum*. *ACS Med. Chem. Lett.* **2(11)**, 866-869 (2011).
2. Nazir, M., Harms, H., Loef, I., *et al.* GPR18 inhibiting amauroamine and the novel triterpene glycoside auxarthonoside from the sponge-derived fungus *Auxarthron reticulatum*. *Planta. Med.* **81(12-13)**, 1141-1145 (2015).
3. Takase, S., Kawai, Y., Uchida, I., *et al.* Structure of amauroamine, a new alkaloid with vasodilating activity produced by *Amauroascus*. sp. *Tetrahedron Lett.* **25(41)**, 4673-4676 (1984).

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, 12/22/2022

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