

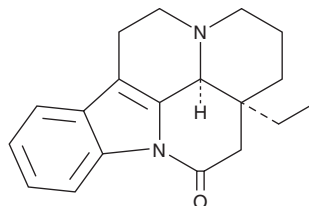
PRODUCT INFORMATION



(-)-Eburnamonine

Item No. 37569

CAS Registry No.: 4880-88-0
Formal Name: (3 α ,16 α)-eburnamenin-14(15H)-one
Synonyms: l-Eburnamonine, Vinburnine, cis-Vincamone, VNB
MF: C₁₉H₂₂N₂O
FW: 294.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 243, 268 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

(-)-Eburnamonine is supplied as a solid. A stock solution may be made by dissolving the (-)-eburnamonine in the solvent of choice, which should be purged with an inert gas. (-)-Eburnamonine is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Description

(-)-Eburnamonine is an indole alkaloid that has been found in *K. jasminiflora* and has vasodilator and neuroprotective activities.¹⁻³ It increases cerebral blood flow and capillary circulation in anesthetized dogs when administered at a dose of 2.5 mg/kg.³ (-)-Eburnamonine (20 mg/kg) prevents scopolamine-induced amnesia in a passive avoidance test in mice.² It enhances acquisition of the conditioned avoidance response induced by hypobaric hypoxia in rats when administered at a dose of 20 mg/kg. (-)-Eburnamonine (0.25 mg/kg per minute) increases the onset time to hypoxia-induced brain death in rats.³ It increases survival time and inhibits decreases in body temperature induced by hypobaric hypoxia in rats when administered at a dose of 45 mg/kg.

References

1. Kitajima, M., Anbe, M., Kogure, N., *et al.* Indole alkaloids from *Kopsia jasminiflora*. *Tetrahedron* **70**(47), 9099-9106 (2014).
2. Drago, F., Grassi, M., Valerio, C., *et al.* Effects of vinburnine on experimental models of learning and memory impairments. *Pharmacol. Biochem. Behav.* **37**(1), 53-57 (1990).
3. Linée, P., Lacroix, P., Le Polles, J.B., *et al.* Cerebral metabolic, hemodynamic and antihypoxic properties of l-eburnamonine. *Eur. Neurol.* **17**(Suppl 1), 113-120 (1978).

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, 08/08/2023

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