PRODUCT INFORMATION



Vinpocetine

Item No. 19578

CAS Registry No.: Formal Name:	42971-09-5 (3α,16α)-eburnamenine-14-	0,0
Synonyms:	carboxylic acid, ethyl ester Apovincaminic Acid ethyl ester, AY 27255, RGH-4405	
MF:	$C_{22}H_{26}N_2O_2$	
FW:	350.5	
Purity:	≥98%	
UV/Vis.:	λ _{may} : 229, 274, 314 nm	L N J
Supplied as:	A crystalline solid	\sim
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

Vinpocetine is supplied as a crystalline solid. A stock solution may be made by dissolving the vinpocetine in the solvent of choice, which should be purged with an inert gas. Vinpocetine is soluble in the organic solvent dimethyl formamide at a concentration of approximately 0.25 mg/ml.

Description

Vinpocetine is a semisynthetic alkaloid commonly used as an inhibitor of phosphodiesterase 1 (PDE1), blocking the hydrolysis of both cAMP and cGMP (K_i = 14 μ M).^{1,2} In addition to its use as a selective PDE1 inhibitor in basic research, vinpocetine has diverse cerebral and neurological effects in vivo.² Vinpocetine also directly inhibits the kinase activity of IKK β (IC₅₀ = 17 μ M in a cell-free system) and blocks TNF- α - and LPS-mediated activation of NF- κ B in cells and *in vivo*.³

References

- 1. Manallack, D.T., Hughes, R.A., and Thompson, P.E. The next generation of phosphodiesterase inhibitors: Structural clues to ligand and substrate selectivity of phosphodiesterases. J. Med. Chem. 48(10), 3449-3462 (2005).
- 2. Medina, A.E. Therapeutic utility of phosphodiesterase type I inhibitors in neurological conditions. Front. Neurosci. 5(21), (2011).
- 3. Jeon, K.-I., Xu, X., Aizawa, T., et al. Vinpocetine inhibits NF-кB-dependent inflammation via an IKK-dependent but PDE-independent mechanism. Proc. Natl. Acad. Sci. USA 107(21), 9795-9800 (2016).

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.

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, 11/16/2022

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