PRODUCT INFORMATION



Isorhynchophylline

Item No. 29650

CAS Registry No.:	6859-01-4	
Formal Name:	(αΕ,1′S,6′R,7′S,8′aS)-6′-ethyl-	
	1,2,2',3',6',7',8',8'a-octahydro-α-	0,
	(methoxymethylene)-2-oxo-spiro[3H-indole-	H
	3,1'(5'H)-indolizine]-7'-acetic acid, methyl ester	
Synonym:	(+)-lsorhynchophylline	
MF:	C ₂₂ H ₂₈ N ₂ O ₄	
FW:	384.5	
Purity:	≥95%	•
UV/Vis.:	λ _{max} : 242 nm	
Supplied as:	A solid	
Storage:	-20°C	Ō
Stability:	≥4 years	
Item Origin:	Plant/Uncaria rhynchophylla	
Informer attack managements		

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

Laboratory Procedures

Isorhynchophylline is supplied as a solid. A stock solution may be made by dissolving the isorhynchophylline in the solvent of choice, which should be purged with an inert gas. Isorhynchophylline is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of isorhynchophylline in ethanol is approximately 1 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Isorhynchophylline is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isorhynchophylline should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Isorhynchophylline has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Isorhynchophylline is an oxindole alkaloid that has been found in Uncaria and has diverse biological activities, including neuroprotective, anticancer, and antihypertensive properties.¹⁻⁴ It inhibits decreases in cell viability and increases in intracellular levels of reactive oxygen species (ROS) and malondialdehyde (MDA) induced by amyloid-β (25-35) (Aβ (25-35); Item No. 24155) in PC12 cells in a concentration-dependent manner.¹ Isorhynchophylline (30 and 100 μ M) attenuates decreases in population spike amplitudes in the CA1 region of a rat hippocampal slice model of ischemia induced by oxygen and D-glucose deprivation.² It decreases the viability of HepG2, A549, BxPC-3, Caki-1, RPMI-8226, and 786-O cells (IC₅₀s = 130-292 μ M).³ In HepG2 cells, isorhynchophylline induces apoptosis, inhibits cell migration in a wound healing assay, and decreases invasion in a Matrigel™ assay. Isorhynchophylline (0.1% in the diet) decreases right ventricular end systolic pressure, hypertrophy, and fibrosis in a rat model of pulmonary arterial hypertension induced by monocrotaline (Item No. 16666).⁴

References

- 1. Xian, Y.-F., Lin, Z.-X., Mao, Q.-Q., et al. Cell Mol. Neurobiol. 32(3), 353-360 (2012).
- 2. Kang, T.-H., Murakami, Y., Takayama, H., et al. Life Sci. 76(3), 331-343 (2004).
- 3. Lee, H., Baek, S.H., Lee, J.H., et al. Int. J. Mol. Sci. 18(5), E1095 (2017).
- 4. Guo, H., Zhang, H., Cui, Y., et al. Biochem. Biophys. Res. Commun. 450(1), 729-734 (2014).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 10/20/2022

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