

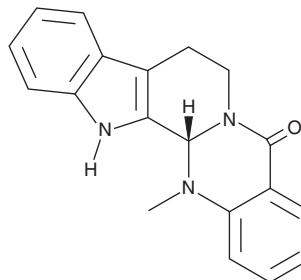
PRODUCT INFORMATION



(+)-Evodiamine

Item No. 16885

CAS Registry No.: 518-17-2
Formal Name: 8,13,13b,14-tetrahydro-14-methyl-indolo[2',3':3,4]pyrido[2,1-b]quinazolin-5(7H)-one
Synonym: D-Evodiamine
MF: C₁₉H₁₇N₃O
FW: 303.4
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 270 nm
Supplied as: A crystalline 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

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

Description

Evodiamine is a natural indole alkaloid found in the fruits of Wu Zhu Yu, a plant used in traditional Chinese medicine. Studies involving evodiamine have demonstrated beneficial effects in cancer, obesity, inflammation, and many other conditions.¹⁻³ It can act as an aryl hydrocarbon antagonist (K_i = 28 nM), activator of the transient receptor potential vanilloid 1 channel (EC₅₀ = 45 nM), and inhibit signaling through NF-κB.^{1,4,5}

References

1. Yu, H., Gong, W., Wang, Z., *et al.* Pharmacological actions of multi-target-directed evodiamine. *Molecules* **18**(2), 1826-1843 (2013).
2. Wang, T., Wang, X., and Yamashita, H. Evodiamine inhibits adipogenesis via the EGFR-PKCα-ERK signaling pathway. *FEBS Lett.* **583**(22), 3655-3659 (2009).
3. Ogasawara, M., Matsubara, T., and Suzuki, H. Screening of natural compounds for inhibitory activity on colon cancer cell migration. *Biol. Pharm. Bull.* **24**(6), 720-723 (2001).
4. Pearce, L.V., Petukhov, P.A., Szabo, T., *et al.* Evodiamine functions as an agonist for the vanilloid receptor TRPV1. *Org. Biomol. Chem.* **2**, 2281-2286 (2004).
5. Takada, Y., Kobayashi, Y., and Aggarwal, B.B. Evodiamine abolishes constitutive and inducible NF-κB activation by inhibiting IκBα kinase activation, thereby suppressing NF-κB-regulated antiapoptotic and metastatic gene expression, up-regulating apoptosis, and inhibiting invasion. *J. Biol. Chem.* **280**(17), 17203-17212 (2005).

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, 12/15/2022

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