

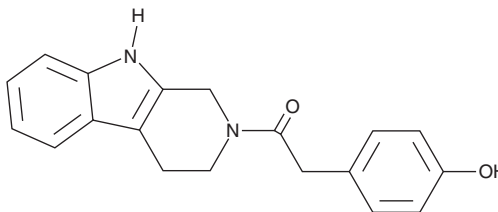
PRODUCT INFORMATION



YH-306

Item No. 41874

CAS Registry No.: 1373764-75-0
Formal Name: 2-(4-hydroxyphenyl)-1-(1,3,4,9-tetrahydro-2H-pyrido[3,4-b]indol-2-yl)-ethanone
MF: C₁₉H₁₈N₂O₂
FW: 306.4
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

YH-306 is supplied as a solid. A stock solution may be made by dissolving the YH-306 in the solvent of choice, which should be purged with an inert gas. YH-306 is soluble (≥10 mg/ml) in DMSO.

Description

YH-306 is an anticancer agent.¹ It reduces TGF-β-induced transcriptional activity by 72.4% in a reporter assay using HaCaT cells when used at a concentration of 1 μM. YH-306 inhibits the proliferation of PC-9, H1299, and A549 lung cancer cells (IC₅₀s = 5.87, 8.07, and 7.9 μM, respectively). It decreases migration of H1299 cells in a wound healing assay when used at concentrations of 0.1, 0.5, or 1 μM. YH-306 (50 μM) inhibits actin polymerization in HT-29 colon cancer cells.² It inhibits fibronectin-induced increases in the levels of matrix metalloproteinase-2 (MMP-2), MMP-9, and phosphorylated focal adhesion kinase (FAK), PI3K, and ERK in CT26 colon cancer cells when used at a concentration of 50 μM. YH-306 (5 mg/kg per day) decreases tumor volume without reducing body weight in an A549 mouse xenograft model.¹ It also reduces tumor growth and the number of hepatic metastases in a CT26-luc colon cancer mouse xenograft model with hepatic metastases when administered at a dose of 50 mg/kg per day.²

References

1. Zheng, C., Fang, Y., Tong, W., *et al.* Synthesis and biological evaluation of novel tetrahydro-β-carboline derivatives as antitumor growth and metastasis agents through inhibiting the transforming growth factor-β signaling pathway. *J. Med. Chem.* **57**(3), 600-612 (2014).
2. Dai, F., Chen, Y., Huang, L., *et al.* A novel synthetic small molecule YH-306 suppresses colorectal tumour growth and metastasis via FAK pathway. *J. Cell. Mol. Med.* **19**(2), 383-395 (2015).

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, 10/24/2024

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