

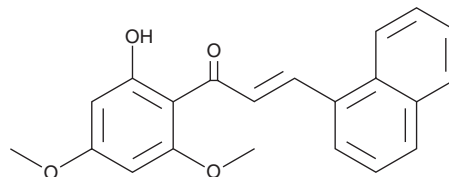
PRODUCT INFORMATION



Flavokawain 1i

Item No. 39394

CAS Registry No.: 1098176-51-2
Formal Name: (2E)-1-(2-hydroxy-4,6-dimethoxyphenyl)-3-(1-naphthalenyl)-2-propen-1-one
Synonym: Flavokavain 1i
MF: C₂₁H₁₈O₄
FW: 334.4
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

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

Description

Flavokawain 1i is a derivative of the chalcones flavokawain A (Item No. 36083), flavokawain B (Item No. 19652), and flavokawain C (Item No. 39393) and has anticancer and antiviral activities.^{1,2} It inhibits the proliferation of gefitinib-resistant H1975 non-small cell lung cancer (NSCLC) cells by 36%, as well as decreases protein levels of the heat shock protein 90 (Hsp90) client proteins EGFR, c-Met, HER2, Akt, and cyclin-dependent kinase 4 (Cdk4) and increases protein levels of Hsp70, markers of Hsp90 inhibition, when used at a concentration of 30 μM.¹ Intramuscular administration of flavokawain 1i reduces viral titers in pigs infected with porcine reproductive and respiratory syndrome virus (PRRSV).²

References

1. Seo, Y.H. and Park, S.Y. Synthesis of flavokawain analogues and their anti-neoplastic effects on drug-resistant cancer cells through Hsp90 inhibition. *Bull. Korean Chem. Soc.* **35**(4), 1154-1158 (2014).
2. Khatun, A., Park, S.Y., Shabir, N., et al. Evaluation of the inhibitory effects of (E)-1-(2-hydroxy-4,6-dimethoxyphenyl)-3-(naphthalen-1-yl)prop-2-en-1-one (DiNap), a natural product analog, on the replication of type 2 PRRSV in vitro and in vivo. *Molecules* **24**(5), 887 (2019).

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, 08/11/2023

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