

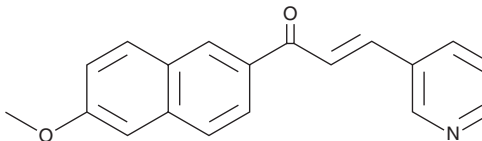
PRODUCT INFORMATION



DMU2139

Item No. 37301

CAS Registry No.: 104890-70-2
Formal Name: 1-(6-methoxy-2-naphthalenyl)-3-(3-pyridinyl)-2-propen-1-one
MF: C₁₉H₁₅NO₂
FW: 289.3
Purity: ≥98%
UV/Vis.: λ_{max}: 232 nm
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

DMU2139 is supplied as a solid. A stock solution may be made by dissolving the DMU2139 in the solvent of choice, which should be purged with an inert gas. DMU2139 is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

DMU2139 is an inhibitor of the cytochrome P450 (CYP) isoform CYP1B1 (IC₅₀ = 0.009 μM).¹ It is selective for CYP1B1 over CYP1A1, CYP1A2, CYP3A4, and CYP2D6 (IC₅₀s = 0.795, 1.2, 10.6, and 5 μM, respectively). DMU2139 enhances cisplatin-induced cell death in HEK293 cells overexpressing CYP1B1.

Reference

1. Horley, N.J., Beresford, K.J.M., Chawla, T., *et al.* Discovery and characterization of novel CYP1B1 inhibitors based on heterocyclic chalcones: Overcoming cisplatin resistance in CYP1B1-overexpressing lines. *Eur. J. Med. Chem.* **129**, 159-174 (2017).

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, 10/04/2023

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