

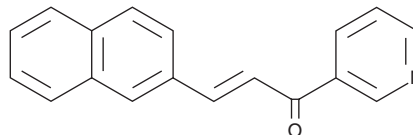
PRODUCT INFORMATION



DMU2105

Item No. 38522

CAS Registry No.: 1031063-36-1
Formal Name: 3-(2-naphthalenyl)-1-(3-pyridinyl)-
2-propen-1-one
MF: C₁₈H₁₃NO
FW: 259.3
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

DMU2105 is supplied as a solid. A stock solution may be made by dissolving the DMU2105 in the solvent of choice, which should be purged with an inert gas. DMU2105 is soluble in the organic solvent methanol.

Description

DMU2105 is an inhibitor of the cytochrome P450 (CYP) isoform CYP1B1 (IC₅₀ = 0.01 μM).¹ It is selective for CYP1B1 over CYP1A1, CYP1A2, CYP3A4, and CYP2D6 (IC₅₀s = 0.742, 1.2, 18, and 10 μM, respectively). DMU2105 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, 11/01/2023

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