

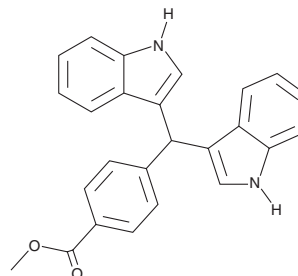
PRODUCT INFORMATION



DIM-C-pPhCO₂Me

Item No. 37540

CAS Registry No.: 151358-48-4
Formal Name: 4-(di-1H-indol-3-ylmethyl)-benzoic acid, methyl ester
MF: C₂₅H₂₀N₂O₂
FW: 380.4
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

DIM-C-pPhCO₂Me is supplied as a solid. A stock solution may be made by dissolving the DIM-C-pPhCO₂Me in the solvent of choice, which should be purged with an inert gas. DIM-C-pPhCO₂Me is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of DIM-C-pPhCO₂Me in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

DIM-C-pPhCO₂Me is a nerve growth factor IB (NR4A1) antagonist.¹ It decreases expression of the mRNA encoding integrin β1 (ITGB1) in MIA PaCa-2 and PANC-1 pancreatic and RKO and SW480 colon cancer cells when used at concentrations of 15 and 20 μM. DIM-C-pPhCO₂Me (15 and 20 μM) decreases the migration and adhesion of MIA PaCa-2 and SW480 cells, effects that can be prevented by siRNA knockdown of NR4A1 or ITGB1.

Reference

1. Hedrick, E., Lee, S.-O., and Safe, S. The nuclear orphan receptor NR4A1 regulates β1-integrin expression in pancreatic and colon cancer cells and can be targeted by NR4A1 antagonists. *Mol. Carcinog.* **56(9)**, 2066-2075 (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, 10/27/2023

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