

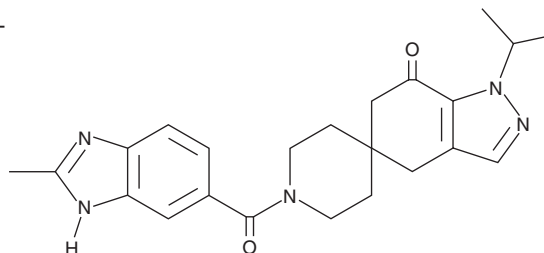
PRODUCT INFORMATION



PF-05175157

Item No. 21778

CAS Registry No.: 1301214-47-0
Formal Name: 1,4-dihydro-1'-[(2-methyl-1H-benzimidazol-6-yl)carbonyl]-1-(1-methylethyl)-spiro[5H-indazole-5,4'-piperidin]-7(6H)-one
MF: C₂₃H₂₇N₅O₂
FW: 405.5
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 249 nm
Supplied as: A crystalline 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

PF-05175157 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-05175157 in the solvent of choice, which should be purged with an inert gas. PF-05175157 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PF-05175157 in these solvents is approximately 25 mg/ml.

Description

PF-05175157 is an inhibitor of acetyl-CoA carboxylase 1 (ACC1) and ACC2 (IC₅₀s = 27, 33, 23.5, and 50.4 nM for human ACC1, human ACC2, rat ACC1, and rat ACC2, respectively).¹ It inhibits formation of malonyl-CoA in primary rat hepatocytes (EC₅₀ = 29.9 nM). *In vivo*, PF-05175157 inhibits formation of malonyl-CoA in the liver and skeletal muscle, inhibits hepatic *de novo* lipogenesis, and lowers the respiratory exchange rate in rats (ED₅₀s = 6, 11, 7.8, and 3.5 mg/kg, respectively).

Reference

1. Griffith, D.A., Kung, D.W., Esler, W.P., *et al.* Decreasing the rate of metabolic ketone reduction in the discovery of a clinical acetyl-CoA carboxylase inhibitor for the treatment of diabetes. *J. Med. Chem.* **57**(24), 10512-10526 (2014).

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, 12/01/2022

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