

PRODUCT INFORMATION

CP 640,186

Item No. 17691



CAS Registry No.: 591778-68-6

Formal Name: [(3R)-1'-(9-anthracenylcarbonyl)-
[1,4'-bipiperidin]-3-yl]-4-morpholinyl-
methanone

Synonym: ACC Inhibitor IV

MF: C₃₀H₃₅N₃O₃

FW: 485.6

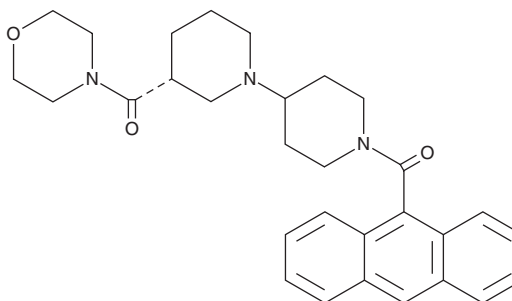
Purity: ≥95%

UV/Vis.: λ_{max}: 254 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

CP 640,186 is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 640,186 in the solvent of choice, which should be purged with an inert gas. CP 640,186 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CP 640,186 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CP 640,186 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of CP 640,186 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Acetyl-CoA carboxylase (ACC) is a key enzyme of fatty acid metabolism that enables the synthesis of malonyl-CoA, a precursor for fatty acids. CP 640,186 is an isozyme-nonspecific ACC inhibitor with IC₅₀ values of 53 and 61 nM for rat liver ACC1 and rat skeletal muscle ACC2, respectively.¹ It has been shown to lower malonyl-CoA levels and to inhibit hepatic fatty acid and triglyceride synthesis both *in vitro* and *in vivo*.¹ CP 640,186 can also stimulate muscle fatty acid oxidation (EC₅₀ = 1.3 μM in rat epitrochlearis muscle strips) and lower whole body respiratory quotient in rats (ED₅₀ = ~30 mg/kg).¹

Reference

1. Harwood, H.J., Jr., Petras, S.F., Shelly, L.D., *et al.* Isozyme-nonspecific N-substituted bipiperidylcarboxamide acetyl-CoA carboxylase inhibitors reduce tissue malonyl-CoA concentrations, inhibit fatty acid synthesis, and increase fatty acid oxidation in cultured cells and in experimental animals. *J. Biol. Chem.* **278**(39), 37099-37111 (2016).

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, 12/19/2022

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