

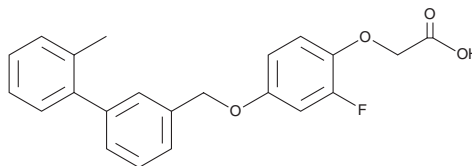
PRODUCT INFORMATION



HWL-088

Item No. 30227

CAS Registry No.: 2378617-96-8
Formal Name: 2-[2-fluoro-4-[(2'-methyl[1,1'-biphenyl]-3-yl)methoxy]phenoxy]-acetic acid
MF: C₂₂H₁₉FO₄
FW: 366.4
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

HWL-088 is supplied as a solid. A stock solution may be made by dissolving the HWL-088 in the solvent of choice, which should be purged with an inert gas. HWL-088 is soluble in acetonitrile.

Description

HWL-088 is an agonist of free fatty acid receptor 1 (FFAR1/GPR40) and peroxisome proliferator-activated receptor δ (PPAR δ).¹ It induces calcium mobilization in CHO cells expressing human FFAR1/GPR40 (EC₅₀ = 18.9 nM) and transactivation of PPAR δ in a reporter assay using HEK293 cells (EC₅₀ = 570.9 nM). HWL-088 is selective for FFAR1/GPR40 and PPAR δ over PPAR α and PPAR γ . It decreases plasma glucose, hemoglobin A1c (HbA1c), total cholesterol, triglyceride, and LDL levels, increases plasma insulin levels, and reduces hepatic lipid accumulation in *ob/ob* mice when administered at a dose of 40 mg/kg.

Reference

1. Chen, Y., Ren, Q., Zhou, Z., *et al.* HWL-088, a new potent free fatty acid receptor 1 (FFAR1) agonist, improves glucolipid metabolism and acts additively with metformin in *ob/ob* diabetic mice. *Br. J. Pharmacol.* **177**(10), 2286-2302 (2019).

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, 12/07/2023

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