

PRODUCT INFORMATION



PF-06424439

Item No. 17680

CAS Registry No.: 1469284-79-4

Formal Name: [(3R)-1-[2-[1-(4-chloro-1H-pyrazol-1-yl)cyclopropyl]-3H-imidazo[4,5-b]pyridin-5-yl]-3-piperidinyl]-1-pyrrolidinyl-methanone, monomethanesulfonate

MF: C₂₂H₂₆ClN₇O • CH₃SO₃H

FW: 536.0

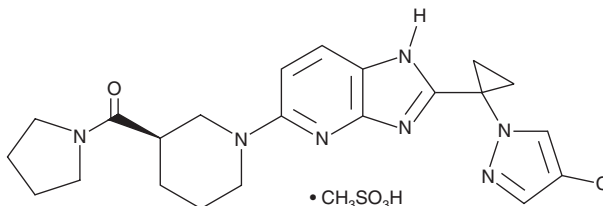
Purity: ≥98%

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.



Description

PF-06424439 is an orally bioavailable inhibitor of diacylglycerol acyltransferase 2 (DGAT-2; IC₅₀ = 14 nM).¹ It is selective for DGAT-2 over the related acyltransferases DGAT-1, MGAT-2, and MGAT-3 (IC₅₀s = >50 μM). PF-06424439 (0.1-10 mg/kg) reduces plasma triglyceride levels in sucrose-fed rats in a dose-dependent manner. It also reduces plasma levels of cholesterol and triglycerides as well as hepatic triglycerides in LDL receptor knockout mice fed a high-fat, high-cholesterol diet when administered at a dose of 60 mg/kg per day.

Reference

1. Futatsugi, K., Kung, D.W., Orr, S.T., *et al.* Discovery and optimization of imidazopyridine-based inhibitors of diacylglycerol acyltransferase 2 (DGAT2). *J. Med. Chem.* **58**(18), 7173-7185 (2015).

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, 12/05/2022

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