

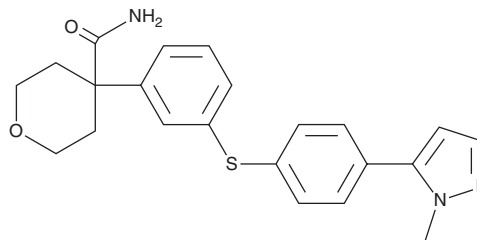
PRODUCT INFORMATION



PF-4191834

Item No. 46096

CAS Registry No.: 1029317-21-2
Formal Name: tetrahydro-4-[[3-[[4-(1-methyl-1H-pyrazol-5-yl)phenyl]thio]phenyl]-2H-pyran-4-carboxamide
Synonym: PF-04191834
MF: C₂₂H₂₃N₃O₂S
FW: 393.5
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

PF-4191834 is supplied as a solid. A stock solution may be made by dissolving the PF-4191834 in the solvent of choice, which should be purged with an inert gas. PF-4191834 is soluble (≥10 mg/ml) in ethanol. PF-4191834 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

PF-4191834 is a non-redox inhibitor of 5-lipoxygenase (5-LO; IC₅₀ = 229 nM).¹ It is selective for 5-LO over 12-LO, 15-LO, COX-1, and COX-2 at 30 μM. PF-4191834 inhibits leukotriene B₄ (LTB₄; Item No. 20110) production in whole blood isolated from patients with asthma (IC₅₀ = 180 nM). It reverses mechanical hyperalgesia in a rat model of inflammatory pain at 1-10 mg/kg.

Reference

1. Masferrer, J.L., Zweifel, B.S., Hardy, M., et al. Pharmacology of PF-4191834, a novel, selective non-redox 5-lipoxygenase inhibitor effective in inflammation and pain. *J. Pharmacol. Exp. Ther.* **334**(1), 294-301 (2010).

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, 08/07/2026

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