

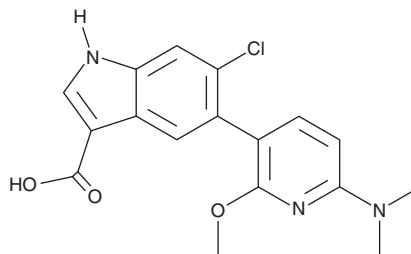
# PRODUCT INFORMATION



## PF-249

Item No. 38416

**CAS Registry No.:** 1467059-70-6  
**Formal Name:** 6-chloro-5-[6-(dimethylamino)-2-methoxy-3-pyridinyl]-1H-indole-3-carboxylic acid  
**Synonym:** PF-06885249  
**MF:** C<sub>17</sub>H<sub>16</sub>ClN<sub>3</sub>O<sub>3</sub>  
**FW:** 345.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 255, 301 nm  
**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-249 is supplied as a solid. A stock solution may be made by dissolving the PF-249 in the solvent of choice, which should be purged with an inert gas. PF-249 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PF-249 in these solvents is approximately 5, 1, and 10 mg/ml, respectively.

### Description

PF-249 is an allosteric activator of AMP-activated protein kinase β1 (AMPKβ1) heterotrimers (EC<sub>50</sub>s = 8 and 6 nM for human AMPK α1β1γ1 and -α2β1γ1, respectively).<sup>1,2</sup> It is selective for AMPK α1β1γ1 and -α2β1γ1 over AMPK α1β2γ1, -α2β2γ1, and -α2β2γ3 at 40 μM and a panel of receptors, channels, and phosphodiesterases at 10 μM.<sup>1,2</sup> PF-249 decreases proteinuria and improves kidney function in an obese ZSF1 rat model of diabetic nephropathy in a dose-dependent manner.<sup>1</sup>

### References

1. Salatto, C.T., Miller, R.A., Cameron, K.O., *et al.* Selective activation of AMPK β1-containing isoforms improves kidney function in a rat model of diabetic nephropathy. *J. Pharmacol. Exp. Ther.* **361**(2), 303-311 (2017).
2. Ryder, T.F., Calabrese, M.F., Walker, G.S., *et al.* Acyl glucuronide metabolites of 6-chloro-5-[4-(1-hydroxycyclobutyl)phenyl]-1H-indole-3-carboxylic acid (PF-06409577) and related indole-3-carboxylic acid derivatives are direct activators of adenosine monophosphate-activated protein kinase (AMPK). *J. Med. Chem.* **61**(16), 7273-7288 (2018).

#### 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, 06/02/2023

#### 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