

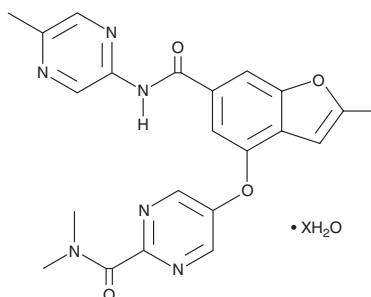
# PRODUCT INFORMATION



## PF-04937319 (hydrate)

Item No. 30926

**Formal Name:** N,N-dimethyl-5-[[[2-methyl-6-[[[(5-methyl-2-pyrazinyl)amino]carbonyl]-4-benzofuranyl]oxy]-2-pyrimidinecarboxamide (hydrate)  
**MF:** C<sub>22</sub>H<sub>20</sub>N<sub>6</sub>O<sub>4</sub> • XH<sub>2</sub>O  
**FW:** 432.4  
**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

PF-04937319 (hydrate) is supplied as a solid. A stock solution may be made by dissolving the PF-04937319 (hydrate) in the solvent of choice, which should be purged with an inert gas. PF-04937319 (hydrate) is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

### Description

PF-04937319 is an activator of glucokinase (EC<sub>50</sub>s = 174, 407, and 269 nM for the recombinant human, rat, and dog enzymes, respectively).<sup>1</sup> It increases glucose-induced insulin secretion in isolated rat islets when used at concentrations of 4 and 12 μM.<sup>2</sup> PF-04937319 (3-100 mg/kg) reduces postprandial plasma glucose levels in rats.

### References

1. Petterson, J.C., Litchfield, J., Neef, N., *et al.* The relationship of glucokinase activator-induced hypoglycemia with arteriopathy, neuronal necrosis, and peripheral neuropathy in nonclinical studies. *Toxicol. Pathol.* **42**(4), 696-708 (2014).
2. Pfefferkorn, J.A., Guzman-Perez, A., Oates, P.J., *et al.* Designing glucokinase activators with reduced hypoglycemia risk: Discovery of N,N-dimethyl-5-(2-methyl-6-((5-methylpyrazin-2-yl)-carbonyl) benzofuran-4-yloxy)pyrimidine-2-carboxamide as a clinical candidate for the treatment of type 2 diabetes mellitus. *Medchemcomm.* **2**, 828-839 (2011).

#### 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, 11/22/2022

#### 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](http://WWW.CAYMANCHEM.COM)