

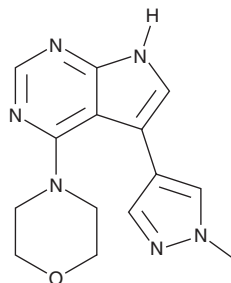
# PRODUCT INFORMATION



**PF-06454589**

Item No. 42957

**CAS Registry No.:** 1527473-30-8  
**Formal Name:** 5-(1-methyl-1H-pyrazol-4-yl)-4-(4-morpholinyl)-7H-pyrrolo[2,3-d]pyrimidine  
**MF:** C<sub>14</sub>H<sub>16</sub>N<sub>6</sub>O  
**FW:** 284.3  
**Purity:** ≥95%  
**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-06454589 is supplied as a solid. A stock solution may be made by dissolving the PF-06454589 in the solvent of choice, which should be purged with an inert gas. PF-06454589 is slightly soluble (0.1-1 mg/ml) in DMSO.

## Description

PF-06454589 is an inhibitor of leucine-rich repeat kinase 2 (LRRK2; IC<sub>50</sub> = 12 nM for the wild-type enzyme).<sup>1</sup> It also inhibits LRRK2 containing a glycine-to-serine substitution at position 2019 (LRRK2<sup>G2019S</sup>; IC<sub>50</sub> = 36 nM).

## Reference

1. Henderson, J.L., Kormos, B.L., Hayward, M.M., *et al.* Discovery and preclinical profiling of 3-[4-(morpholin-4-yl)-7H-pyrrolo[2,3-d]pyrimidin-5-yl]benzotrile (PF-06447475), a highly potent, selective, brain penetrant, and in vivo active LRRK2 kinase inhibitor. *J. Med. Chem* **58**(1), 419-432 (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, 04/01/2025

## 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