

# PRODUCT INFORMATION

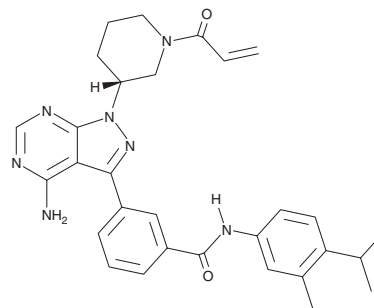


**PF-06465469**

Item No. 31118

**CAS Registry No.:** 1407966-77-1  
**Formal Name:** 3-[4-amino-1-[(3R)-1-(1-oxo-2-propen-1-yl)-3-piperidinyl]-1H-pyrazolo[3,4-d]pyrimidin-3-yl]-N-[3-methyl-4-(1-methylethyl) phenyl]-benzamide

**MF:** C<sub>30</sub>H<sub>33</sub>N<sub>7</sub>O<sub>2</sub>  
**FW:** 523.6  
**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-06465469 is supplied as a solid. A stock solution may be made by dissolving the PF-06465469 in the solvent of choice, which should be purged with an inert gas. PF-06465469 is soluble (≥10 mg/ml) in ethanol and DMSO.

## Description

PF-06465469 is a covalent inhibitor of IL-2-inducible T cell kinase (Itk; IC<sub>50</sub> = 2 nM).<sup>2</sup> It also inhibits Bruton's tyrosine kinase (BTK; IC<sub>50</sub> = 2 nM). PF-06465469 inhibits Itk phosphorylation of phospholipase Cγ (PLCγ) induced by an anti-CD3 antibody in Jurkat cells (IC<sub>50</sub> = 57 nM). It induces apoptosis in Jurkat cells, MOLT-4 and CCRF CEM T cell lymphoblasts, and Karpas-299 large cell lymphoma cells when used at a concentration of 16 μM.<sup>2</sup> PF-06465469 (2 μM) inhibits chemotaxis induced by chemokine (C-X-C motif) ligand 12 (CXCL12) in Jurkat cells.

## References

1. Zapf, C.W., Gerstenberger, B.S., Xing, L., *et al.* Covalent inhibitors of interleukin-2 inducible T cell kinase (Itk) with nanomolar potency in a whole-blood assay. *J. Med. Chem.* **55(22)**, 10047-10063 (2012).
2. Mamand, S., Allchin, R.L., Ahearne, M.J., *et al.* Comparison of interleukin-2-inducible kinase (ITK) inhibitors and potential for combination therapies for T-cell lymphoma. *Sci. Rep.* **8(1)**, 14216 (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, 07/24/2025

## 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