

PRODUCT INFORMATION

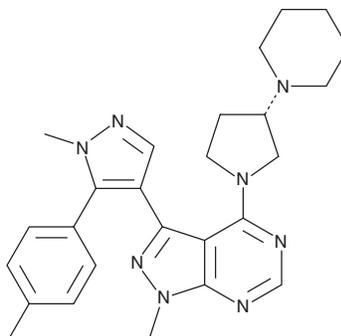


PF-04981517

Item No. 15019

CAS Registry No.: 1390637-82-7
Formal Name: 1-methyl-3-[1-methyl-5-(4-methylphenyl)-1H-pyrazol-4-yl]-4-[(3S)-3-(1-piperidinyl)-1-pyrrolidinyl]-1H-pyrazolo[3,4-d]pyrimidine

Synonym: CYP3cide
MF: C₂₆H₃₂N₈
FW: 456.6
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 295 nm
Supplied as: A crystalline 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-04981517 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-04981517 in the solvent of choice. PF-04981517 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of PF-04981517 in these solvents is approximately 2, 20, and 30 mg/ml, respectively.

PF-04981517 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-04981517 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PF-04981517 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PF-04981517 is a potent, time-dependent inhibitor of the cytochrome P450 (CYP450) isoform CYP3A4 (IC₅₀ = 30 nM).¹ It is less effective against CYP3A5 and CYP3A7 (IC₅₀s = 17 and 70 μM, respectively).¹ PF-04981517 will be useful for understanding the relative roles of CYP3A4 and CYP3A5 in drug metabolism.

Reference

1. Walsky, R.L., Obach, R.S., Hyland, R., *et al.* Selective mechanism-based inactivation of cytochrome P4503A4 by CYP3cide (PF-04981517) and its utility as an in vitro tool for delineating the relative roles of cytochrome P4503A4 versus P4503A5 in the metabolism of drugs. *Drug Metab. Dispos.* **40(9)**, 1686-1697 (2012).

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, 11/11/2022

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