

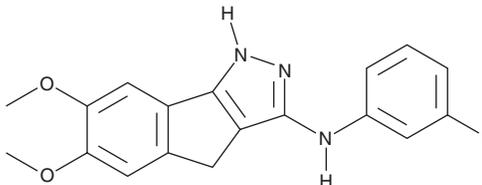
PRODUCT INFORMATION



JNJ-10198409

Item No. 10008131

CAS Registry No.: 627518-40-5
Formal Name: N-(3-fluorophenyl)-1,4-dihydro-6,7-dimethoxy-indeno[1,2-c]pyrazol-3-amine
MF: C₁₈H₁₆FN₃O₂
FW: 325.3
Purity: ≥98%
UV/Vis.: λ_{max}: 267, 302 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

JNJ-10198409 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-10198409 in the solvent of choice, which should be purged with an inert gas. JNJ-10198409 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JNJ-10198409 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Inhibition of the tyrosine kinase activity of growth factor receptors such as the platelet-derived growth factor (PDGF-BB) receptor can have potent antiangiogenic and antiproliferative activity.¹ JNJ-10198409 is a small molecule inhibitor of platelet-derived growth factor (PDGF-BB) tyrosine kinase with an IC₅₀ value of 4.2 nM when tested in human coronary artery smooth muscle cells.² JNJ-10198409 is a competitive antagonist of the ATP binding and hydrolysis at this receptor, resulting in a dose dependent inhibition of tumor growth and angiogenesis.

References

1. Bategay, E.J., Rupp, J., Iruela-Arispe, L., *et al.* PDGF-BB modulates endothelial proliferation and angiogenesis in vitro via PDGF β-receptors. *J. Cell Biol.* **125**(4), 917-928 (1994).
2. Ho, C.Y., Ludovici, D.W., Maharoo, U.S.M., *et al.* (6,7-Dimethoxy-2,4-dihydroindeno[1,2-c]pyrazol-3-yl)phenylamines: Platelet-derived growth factor receptor tyrosine kinase inhibitors with broad antiproliferative activity against tumor cells. *J. Med. Chem.* **48**(26), 8163-8173 (2005).

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, 11/17/2022

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