

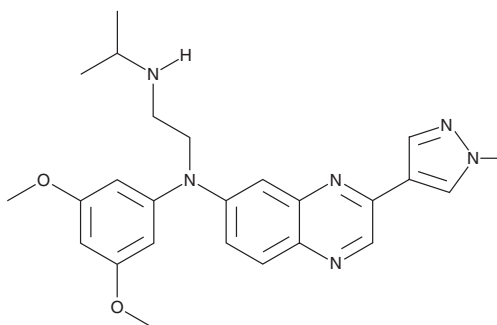
PRODUCT INFORMATION



JNJ-42756493

Item No. 21813

CAS Registry No.: 1346242-81-6
Formal Name: N¹-(3,5-dimethoxyphenyl)-N²-(1-methylethyl)-N¹-[3-(1-methyl-1H-pyrazol-4-yl)-6-quinoxaliny]-1,2-ethanediamine
Synonym: Erdafitinib
MF: C₂₅H₃₀N₆O₂
FW: 446.5
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 257, 297, 411 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-42756493 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-42756493 in the solvent of choice, which should be purged with an inert gas. JNJ-42756493 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JNJ-42756493 in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

JNJ-42756493 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNJ-42756493 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JNJ-42756493 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

JNJ-42756493 is a pan-FGFR inhibitor (IC₅₀s = 1.2, 2.5, 3, and 5.7 nM for FGFR1, -2, -3, and -4, respectively).¹ It is selective for FGFRs over VEGFR2 (IC₅₀ = 36.8 nM). JNJ-42756493 inhibits the proliferation of Ba/F3 cells expressing FGFR1, -3, or -4 (IC₅₀s = 22.1, 13.2, and 25 nM, respectively). It reduces tumor growth in a SNU-16 gastric cancer mouse xenograft model when administered at doses of 10 and 30 mg/kg.

Reference

1. Perera, T.P.S., Jovcheva, E., Mevellec, L., *et al.* Discovery and pharmacological characterization of JNJ-42756493 (erdafitinib), a functionally selective small molecule FGFR family inhibitor. *Mol. Cancer Ther.* **16**(6), 1010-1020 (2017).

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, 10/25/2022

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