

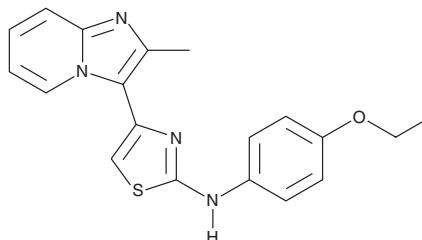
PRODUCT INFORMATION



JK 184

Item No. 15265

CAS Registry No.: 315703-52-7
Formal Name: N-(4-ethoxyphenyl)-4-(2-methylimidazo[1,2-a]pyridin-3-yl)-2-thiazolamine
MF: C₁₉H₁₈N₄O₂
FW: 350.4
Purity: ≥99%
Supplied as: A solid
Storage: 4°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

JK 184 is supplied as a solid. A stock solution may be made by dissolving the JK 184 in the solvent of choice, which should be purged with an inert gas. JK 184 is soluble in organic solvents such as ethanol and DMSO. The solubility of JK 184 in these solvents is approximately 25 mM and 100 mM, respectively.

Description

JK 184 is an imidazopyridine derivative that inhibits downstream hedgehog signaling, preventing Gli-dependent transcriptional activity (IC₅₀ = 30 nM).¹ It does so by inhibiting alcohol dehydrogenase 7 (IC₅₀ = 210 nM) and depolymerizing microtubules, preventing their assembly, which is crucial for Gli function.^{1,2} JK 184 exhibits antiproliferative activity in a range of cancer cell lines (L3.6pl, Panc 05.04, BxPC3, D283 med, Daoy; GI₅₀s = 3-21 nM) and in L3.6pl and BxPC3 pancreatic cancer xenograft models (0.2 mg/mouse/day).¹

References

1. Lee, J., Wu, X., Pasca de Magliano, M., *et al.* A small-molecule antagonist of the hedgehog signaling pathway. *ChemBioChem* **8(16)**, 1916-1919 (2007).
2. Cupido, T., Rack, P.G., Firestone, A.J., *et al.* The imidazopyridine derivative JK184 reveals dual roles for microtubules in hedgehog signaling. *Angew. Chem. Int. Ed. Engl.* **48(13)**, 2321-2324 (2009).

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, 11/04/2022

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