

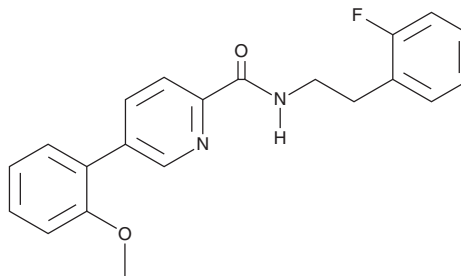
PRODUCT INFORMATION



N-(2-Fluorophenethyl)-5-(2-methoxyphenyl)picolinamide

Item No. 29740

CAS Registry No.: 2380261-53-8
Formal Name: N-[2-(2-fluorophenyl)ethyl]-5-(2-methoxyphenyl)-2-pyridinecarboxamide
MF: C₂₁H₁₉FN₂O₂
FW: 350.4
Purity: ≥98%
UV/Vis.: λ_{max}: 263 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

N-(2-Fluorophenethyl)-5-(2-methoxyphenyl)picolinamide is supplied as a crystalline solid. A stock solution may be made by dissolving the N-(2-fluorophenethyl)-5-(2-methoxyphenyl)picolinamide in the solvent of choice, which should be purged with an inert gas. N-(2-Fluorophenethyl)-5-(2-methoxyphenyl)picolinamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-(2-fluorophenethyl)-5-(2-methoxyphenyl)picolinamide in these solvents is approximately 30 mg/ml.

Description

N-(2-Fluorophenethyl)-5-(2-methoxyphenyl)picolinamide is an inhibitor of hypoxia-inducible factor (HIF-1) signaling (IC₅₀ = 0.32 μM in a cell-based hypoxia responsive element (HRE) reporter assay under hypoxic conditions).¹ It inhibits hypoxia-induced increases in HIF-1α levels in MDA-MB-231, SKOV3, and HeLa cells in a concentration-dependent manner. N-(2-Fluorophenethyl)-5-(2-methoxyphenyl)picolinamide also inhibits hypoxia-induced VEGF expression and capillary-like tube formation in human umbilical vein endothelial cells (HUVECs) and inhibits migration and invasion of MDA-MB-231 cells in wound healing and transwell assays under hypoxic conditions. It also reduces lung metastasis in an MDA-MB-231 mouse xenograft model when administered at doses of 15 and 30 mg/kg every other day.

Reference

1. Liu, M., Liang, Y., Zhu, Z., *et al.* Discovery of novel aryl carboxamide derivatives as hypoxia-inducible factor 1α signaling inhibitors with potent activities of anticancer metastasis. *J. Med. Chem.* **62**(20), 9299-9314 (2019).

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, 01/03/2020

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