

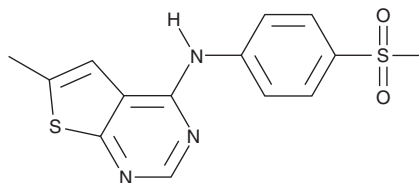
PRODUCT INFORMATION



ARUK2001607

Item No. 41527

CAS Registry No.: 2924824-56-4
Formal Name: 6-methyl-N-[4-(methylsulfonyl)phenyl]-thieno[2,3-d]pyrimidin-4-amine
MF: C₁₄H₁₃N₃O₂S₂
FW: 319.4
Purity: ≥98%
Supplied as: A 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

ARUK2001607 is supplied as a solid. A stock solution may be made by dissolving the ARUK2001607 in the solvent of choice, which should be purged with an inert gas. ARUK2001607 is soluble (≥10 mg/ml) in DMSO.

Description

ARUK2001607 is an inhibitor of phosphatidylinositol 5-phosphate 4-kinase γ (PI5P4K γ ; IC₅₀ = 0.079 μ M for PI5P4K γ +, a construct with mutations conferring increased functional activity).¹ It selectively binds to PI5P4K γ (K_d = 7.1 nM) over 22 other lipid kinases but does bind to phosphatidylinositol 4-phosphate 5-kinase type-1 γ (PIP5K1C; K_d = 230 nM). ARUK2001607 is selective for PI5P4K γ over PI5P4K α and PI5P4K β (IC₅₀s = >15.8 and >25.1 μ M, respectively). It is also selective for PI5P4K γ over 138 other kinases in a kinase panel and a variety of receptors, enzymes, and ion channels in a safety panel at 10 μ M but does inhibit Aurora B kinase and CDC-like kinase 2 (CLK2) by 31 and 37% and dopamine uptake by 59% at 10 μ M.

Reference

1. Rooney, T.P.C., Aldred, G.G., Boffey, H.K., *et al.* The identification of potent, selective, and brain penetrant PI5P4K γ inhibitors as in vivo-ready tool molecules. *J. Med. Chem.* **66**(1), 804-821 (2023).

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, 08/30/2024

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