

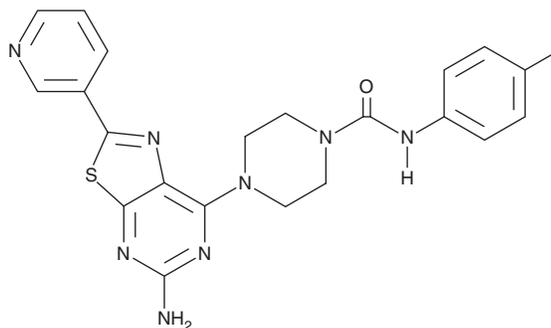
PRODUCT INFORMATION



PI4KIII β Inhibitor 3

Item No. 39530

CAS Registry No.: 1245319-54-3
Formal Name: 4-[5-amino-2-(3-pyridinyl)thiazolo[5,4-d]pyrimidin-7-yl]-N-(4-methylphenyl)-1-piperazinecarboxamide
MF: C₂₂H₂₂N₈OS
FW: 446.5
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

PI4KIII β inhibitor 3 is supplied as a solid. A stock solution may be made by dissolving the PI4KIII β inhibitor 3 in the solvent of choice, which should be purged with an inert gas. PI4KIII β inhibitor 3 is soluble in DMSO and methanol.

Description

PI4KIII β inhibitor 3 is an inhibitor of PI4KIII β (IC₅₀ = 19 nM).¹ It is approximately 100-fold selective for PI4KIII β over PI4KII α , PI4KII β , and PI4KIII α , PI3K α , PI3K β , PI3K γ , and PI3K δ , and PI3KC3, as well as 100-500-fold selective over PI3KC2 α , PI3KC2 β , and PI3KC2 γ .

Reference

1. Reuberson, J., Horsley, H., Franklin, R.J., *et al.* Discovery of a potent, orally bioavailable PI4KIII β inhibitor (UCB9608) able to significantly prolong allogeneic organ engraftment *in vivo*. *J. Med. Chem.* **61**(15), 6705-6723 (2018).

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, 10/26/2023

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