

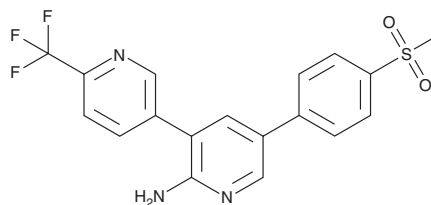
PRODUCT INFORMATION



MMV390048

Item No. 41835

CAS Registry No.: 1314883-11-8
Formal Name: 5-[4-(methylsulfonyl)phenyl]-6'-(trifluoromethyl)-[3,3'-bipyridin]-2-amine
MF: C₁₈H₁₄F₃N₃O₂S
FW: 393.4
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

MMV390048 is supplied as a solid. A stock solution may be made by dissolving the MMV390048 in the solvent of choice, which should be purged with an inert gas. MMV390048 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MMV390048 is an inhibitor of *P. vivax* PI4K (IC₅₀ = 3.4 nM).¹ It is active against *P. falciparum* (IC₅₀ = 28 nM). MMV390048 decreases the blood levels of *P. falciparum* in *P. falciparum*-infected mice (ED₉₀ = 0.57 mg/kg). It reduces the hepatic levels of hypnozoite- and schizont-stage *P. cynomolgi* parasites in *P. cynomolgi*-infected rhesus monkeys when administered at a dose of 20 mg/kg per day.

Reference

1. Paquet, T., Le Manach, C., Cabrera, D.G., *et al.* Antimalarial efficacy of MMV390048, an inhibitor of *Plasmodium* phosphatidylinositol 4-kinase. *Sci. Transl. Med.* **9**(387), eaad9735 (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, 10/30/2024

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